

Technical Appendix

Restriction Enzyme Activity in Promega 10X Buffers, Reaction Temperature and Heat Inactivation.

The 10X Reaction Buffer supplied with each restriction enzyme is optimized to give 100% activity. In many cases good activity is also obtained using one of Promega's 4-CORE® 10X Buffers. Many commonly used cloning enzymes (i.e. restriction sites are found in vector multiple cloning sites) have buffers E and H as their optimal buffer and so we have determined the activity of many of our other restriction enzymes in those buffers as well. This table may be used to select the best buffer for digestion with multiple restriction enzymes. Enzyme activity is expressed as a percent of the activity obtained with the optimal buffer for each enzyme in a one-hour digest. Enzymes with 100% activity (green) perform as well as the optimal buffer. Enzymes with 50–75% or 75–100% (yellow) will give acceptable activity in that buffer. Enzymes with <10%, 10–25%, or 25–50% (pink) activity generally do not have acceptable activity in that buffer. Also, buffers leading to star activity of the enzyme (*; pink) should be avoided. If compatible buffers cannot be identified with acceptable activity for both enzymes, each digest should be performed separately in the optimal buffer for each enzyme.

Promega Enzyme	Cat. #	Buffer Supplied with Enzyme	Activity in							MULTI-CORE™	Heat Inactivation	Enzyme Assay Temperature
			A	B	C	D	E	H				
Aat II	R6541	J	50–75%	10–25%	<10%	<10%	10–25%	<10%	<10%	+	37°C	
Acc I	R6411	G	50–75%	25–50%	25–50%	10–25%	<10%	<10%	25–50%	–	37°C	
Acc III	R6581	F	<10%	10–25%	25–50%	25–50%	n.d.	n.d.	<10%	–	65°C	
Acc65 I	R6921	D	10–25%	50–75%	75–100%	100%	75–100%	100–125%**	100%	+	37°C	
AccB7 I	R7081	E	10–25%	50–75%	100%*	<10%	100%	n.d.	100%	+	37°C	
Age I	R7251	K	25–50%	25–50%	25–50%	50–75%	n.d.	n.d.	100%	+	37°C	
Alu I	R6281	B	75–100%	100%	75–100%	10–25%	n.d.	n.d.	10–25%	+	37°C	
Alw26 I	R6761	C	10–25%	25–50%	100%	10–25%	n.d.	n.d.	75–100%	+	37°C	
Alw44 I	R6771	C	<10%	25–50%	100%	25–50%	n.d.	n.d.	100%	+	37°C	
Apa I	R6361	A	100%	50–75%	50–75%	<10%	10–25%	<10%	75–100%	+	37°C	
Ava I	R6091	B	10–25%	100%	50–75%	25–50%	100%	10–25%	<10%	+/–	37°C	
Ava II	R6131	C	50–75%	50–75%	100%	25–50%	n.d.	n.d.	25–50%	+	37°C	
Bal I	R6691	G	10–25%	<10%	<10%	<10%	n.d.	n.d.	<10%	+	37°C	
BamH I	R6021	E	75–100%*	75–100%	75–100%	50–75%	100%	50–75%	75–100%	+	37°C	
Ban I	R6891	G	25–50%	25–50%	10–25%	<10%	n.d.	n.d.	100%	–	50°C	
Ban II	R6561	E	75–100%	75–100%	75–100%	25–50%	n.d.	n.d.	100%	+	37°C	
Bbu I	R6621	A	100%	75–100%	75–100%	<10%	10–25%	10–25%	100%	+	37°C	
Bcl I	R6651	C	10–25%	75–100%	100%	50–75%	50–75%	50–75%	10–25%	–	50°C	
Bgl I	R6071	D	10–25%	25–50%	75–100%	100%	25–50%	75–100%	100%	+	37°C	
Bgl II	R6081	D	25–50%	75–100%	75–100%	100%	n.d.	n.d.	<10%	–	37°C	
BsaM I	R6991	D	10–25%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	–	65°C	
Bsp1286 I	R6741	A	100%	50–75%	25–50%	10–25%	n.d.	n.d.	75–100%	+	37°C	
BsrS I	R7241	D	10–25%	25–50%	10–25%	100%	n.d.	n.d.	100%	–	65°C	
BssH II	R6831	H	75–100%	50–75%	75–100%	50–75%	n.d.	100%	75–100%	–	50°C	
Bst98 I	R7141	D	<10%	10–25%	10–25%	100%	n.d.	n.d.	25–50%	–	37°C	
BstE II	R6641	D	25–50%	50–75%	50–75%	100%	n.d.	n.d.	100%	–	60°C	
BstO I	R6931	C	10–25%	25–50%	100%	25–50%	n.d.	n.d.	<10%	–	60°C	
BstX I	R6471	D	<10%	10–25%	25–50%	100%	100%	75–100%	10–25%	+/–	50°C	
BstZ I	R6881	D	<10%	<10%	10–25%	100%	10–25%	75–100%	10–25%	–	50°C	
Bsu36 I	R6821	E	<10%	25–50%	50–75%	25–50%	100%	n.d.	50–75%	–	37°C	
Cfo I	R6241	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	100%	+/–	37°C	
Cla I	R6551	C	75–100%	75–100%	100%	75–100%	100%	50–75%	100%	+	37°C	
Csp I	R6671	K	<10%	10–25%	25–50%	50–75%	100%	100–125%**	10–25%	+	30°C	
Csp45 I	R6571	B	25–50%	100%	50–75%	25–50%	100%	25–50%	50–75%	+	37°C	
Dde I	R6291	D	25–50%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	+/–	37°C	
Dpn I	R6231	B	50–75%	100%	75–100%	50–75%	n.d.	n.d.	100%	+	37°C	
Dra I	R6271	B	75–100%	100%	75–100%	50–75%	n.d.	n.d.	25–50%	+	37°C	
EclHK I	R7111	E	<10%	<10%	75–100%	10–25%	100%	n.d.	50–75%	+	37°C	
Eco47 III	R6731	D	<10%	25–50%	50–75%	100%	n.d.	n.d.	25–50%	+	37°C	
Eco52 I	R6751	L	<10%	<10%	10–25%	25–50%	25–50%	50–75%	<10%	+	37°C	
EcoICR I	R6951	B	10–25%	100%	75–100%	<10%	25–50%	n.d.	100%	+	37°C	
EcoR I	R6011	H	25–50%	50–75%	50–75%	50–75%	75–100%	100%	100%*	+	37°C	
EcoR V	R6351	D	10–25%	25–50%	50–75%	100%	25–50%	50–75%	100%	+	37°C	
Fok I	R6781	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	50–75%	+	37°C	
Hae II	R6661	B	50–75%	100%	50–75%	10–25%	n.d.	n.d.	100%	–	37°C	
Hae III	R6171	C	75–100%	75–100%	100%	50–75%	n.d.	n.d.	100%	–	37°C	
Hha I	R6441	C	50–75%	75–100%	100%	50–75%	n.d.	n.d.	75–100%	+	37°C	
Hinc II	R6031	B	25–50%	100%	25–50%	50–75%	75–100%	50–75%	100%	+	37°C	
Hind III	R6041	E	25–50%	100%	75–100%	10–25%	100%	25–50%	50–75%	+	37°C	
Hint I	R6201	B	50–75%	100%	75–100%	75–100%	n.d.	n.d.	50–75%	–	37°C	

Technical Appendix

Restriction Enzyme Activity in Promega 10X Buffers, Reaction Temperature and Heat Inactivation (continued).

Promega Enzyme	Cat.#	Buffer Supplied with Enzyme	Activity in							MULTI-CORE™	Heat Inactivation	Enzyme Assay Temperature
			A	B	C	D	E	H				
<i>Hpa</i> I	R6301	J	25–50%	50–75%	25–50%	10–25%	n.d.	n.d.	100%	—	37°C	
<i>Hpa</i> II	R6311	A	100%	50–75%	50–75%	10–25%	n.d.	n.d.	100%	—	37°C	
<i>Hsp92</i> I	R7151	F	10–25%	75–100%	50–75%	25–50%	n.d.	n.d.	10–25%	+	37°C	
<i>Hsp92</i> II	R7161	K	10–25%	25–50%	25–50%	<10%	n.d.	n.d.	<10%	+	37°C	
<i>I-Ppo</i> I	R7031	NA	10–25%	25–50%	25–50%	25–50%	n.d.	n.d.	—	+	37°C	
<i>Kpn</i> I	R6341	J	100%*	25–50%	25–50%	<10%	25–50%	<10%	75–100%	+/–	37°C	
<i>Mbo</i> I	R6711	C	10–25%	75–100%	100%	50–75%	n.d.	n.d.	<10%	+	37°C	
<i>Mbo</i> II	R6723	B	10–25%	100%	50–75%	75–100%	n.d.	n.d.	100%	+	37°C	
<i>Mlu</i> I	R6381	D	10–25%	25–50%	50–75%	100%	25–50%	100–125%**	10–25%	+/–	37°C	
<i>Msp</i> I	R6401	B	75–100%	100%	75–100%	25–50%	n.d.	n.d.	25–50%	+	37°C	
<i>MspA1</i> I	R7021	C	25–50%	100%*	100%	10–25%	n.d.	n.d.	100%	+	37°C	
<i>Nae</i> I	R7131	A	100%	50–75%	25–50%	<10%	n.d.	n.d.	50–75%	+	37°C	
<i>Nar</i> I	R6861	G	75–100%	50–75%	75–100%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Nci</i> I	R7061	B	100%*	100%	25–50%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Nco</i> I	R6513	D	50–75%	75–100%	75–100%	100%	100%	100–125%**	75–100%	+	37°C	
<i>Nde</i> I	R6801	D	<10%	<10%	25–50%	100%	n.d.	n.d.	25–50%	+	37°C	
<i>Nde</i> II	R7291	D	<10%	<10%	10–25%	100%	n.d.	n.d.	25–50%	+	37°C	
<i>Ngo</i> M IV	R7171	MULTI-CORE™	100%*	100%*	100%*	<10%	n.d.	n.d.	100%	+	37°C	
<i>Nhe</i> I	R6501	B	75–100%	100%	75–100%	10–25%	75–100%	10–25%	100%	+	37°C	
<i>Not</i> I	R6431	D	<10%	10–25%	25–50%	100%	25–50%	100–125%**	25–50%	+	37°C	
<i>Nru</i> I	R7091	K	<10%	<10%	<10%	50–75%	n.d.	n.d.	10–25%	+	37°C	
<i>Nsi</i> I	R6531	D	10–25%	50–75%	50–75%	100%	25–50%	>125%**	10–25%	+/–	37°C	
<i>Pst</i> I	R6111	H	10–25%	50–75%	50–75%	50–75%	25–50%	100%	25–50%	+	37°C	
<i>Pvu</i> I	R6321	D	10–25%	25–50%	50–75%	100%	n.d.	n.d.	<10%	–	37°C	
<i>Pvu</i> II	R6331	B	25–50%	100%	50–75%	25–50%	n.d.	n.d.	50–75%	+	37°C	
<i>Rsa</i> I	R6371	C	75–100%	75–100%	100%	<10%	n.d.	n.d.	<10%	+	37°C	
<i>Sac</i> I	R6061	J	75–100%	25–50%	25–50%	<10%	100%	25–50%	100%	+	37°C	
<i>Sac</i> II	R6221	C	100%	50–75%	100%	50–75%	25–50%	>125%**	<10%	+	37°C	
<i>Sal</i> I	R6051	D	<10%	10–25%	25–50%	100%	25–50%	25–50%	<10%	+	37°C	
<i>Sau</i> 3A I	R6191	B	25–50%	100%	75–100%	<10%	n.d.	n.d.	100%	+	37°C	
<i>Sca</i> I	R6211	K	<10%	100%*	50–75%	75–100%	n.d.	n.d.	10–25%	+	37°C	
<i>Sfi</i> I	R6391	B	75–100%	100%	75–100%	25–50%	75–100%	50–75%	75–100%	–	50°C	
<i>Sgf</i> I	R7103	C	25–50%	25–50%	100%	<10%	n.d.	n.d.	<10%	+/–	37°C	
<i>Sin</i> I	R6141	A	100%	75–100%	50–75%	10–25%	n.d.	n.d.	100%	+	37°C	
<i>Sma</i> I	R6121	J	<10%	<10%	<10%	<10%	<10%	<10%	100%	+	25°C	
<i>Sna</i> B I	R6791	B	50–75%	100%	50–75%	<10%	n.d.	n.d.	100%	–	37°C	
<i>Spe</i> I	R6591	B	75–100%	100%	75–100%	75–100%	100%	25–50%	100%	+	37°C	
<i>Sph</i> I	R6261	K	75–100%	75–100%	100%*	75–100%	100%	>125%**	10–25%	+	37°C	
<i>Ssp</i> I	R6601	E	10–25%	50–75%	50–75%	75–100%	100%	100–125%**	50–75%	+	37°C	
<i>Stu</i> I	R6421	B	75–100%	100%	75–100%	50–75%	n.d.	n.d.	50–75%	+	37°C	
<i>Sty</i> I	R6481	F	25–50%	75–100%	75–100%	75–100%	10–25%	50–75%	<10%	+	37°C	
<i>Taq</i> I	R6151	E	10–25%	25–50%	50–75%	50–75%	100%	n.d.	100%	–	65°C	
<i>Tru</i> 9 I	R7011	F	75–100%	50–75%	75–100%	25–50%	n.d.	n.d.	25–50%	–	65°C	
<i>Tth</i> 111 I	R6841	B	50–75%	100%	75–100%	25–50%	n.d.	n.d.	100%	–	65°C	
<i>Vsp</i> I	R6851	D	<10%	25–50%	75–100%	100%	n.d.	n.d.	<10%	+	37°C	
<i>Xba</i> I	R6181	D	50–75%	75–100%	75–100%	100%	100%	100–125%**	100%	–	37°C	
<i>Xho</i> I	R6161	D	25–50%	75–100%	75–100%	100%	25–50%	100–125%**	10–25%	+	37°C	
<i>Xho</i> II	R6811	C	25–50%	25–50%	100%	10–25%	n.d.	n.d.	<10%	+	37°C	
<i>Xma</i> I	R6491	B	50–75%	100%	25–50%	<10%	25–50%	<10%	50–75%	+	37°C	
<i>Xmn</i> I	R7271	B	75–100%	100%	75–100%	10–25%	n.d.	n.d.	75–100%	+	37°C	

* Not recommended due to potential star activity.

** Unit activity is based on recommended buffer. In Buffer H, some enzymes have enhanced activity.

n.d. = Not determined.

Heat Inactivation Key:

+	=greater than 95% inactivation (DNA is undigested)
—	=less than 95% inactivation (DNA digest is complete, i.e., ≥5% of the initial 20 activity units [≥1 unit] remains)
+/-	=partial inactivation (DNA is partially digested)

Heat Inactivation Conditions:

Twenty units of enzyme in 50µl of its optimal buffer were heated at 65°C for 15 minutes.

One microgram of DNA was added and incubated for 1 hour in accordance with the unit definition, then analyzed by agarose gel electrophoresis.

Technical Appendix

Isoschizomers. The enzymes in boldface type are available from Promega.

Enzyme	Isoschizomer(s)	Recognition Sequence
<i>mAat</i> I	Stu I, <i>Eco</i> 147 I, <i>Pme</i> 55 I, <i>Sse</i> B I	AGG▼CCT
Aaf II	—	GACGT▼C
Acc I	<i>Fbl</i> I, <i>Xmi</i> I	GT▼(A/C)(G/T)AC
Acc III	<i>Bsp</i> E II, <i>Mro</i> I	T▼CCGGA
Acc65 I	<i>Asp</i> 718 I Kpn I*	G▼GTACC GGTAC▼C
<i>Acc</i> B1 I	Ban I, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(C/T)(G/A)CC
AccB7 I	<i>Pfi</i> M I, <i>Van</i> 91 I	CCAN ₄ ▼NTGG
<i>Acc</i> N I	Spe I	A▼CTAGT
<i>Acc</i> W I	<i>Alw</i> I	GGATCNNNN▼
<i>Acy</i> I	<i>Bbi</i> II, <i>Hin</i> 1 I, Hsp92 I, <i>Bsa</i> H I, <i>Msp</i> 171 I	G(A/G)▼CG(T/C)C
<i>Acs</i> I	<i>Apo</i> I	(G/A)▼AATT(C/T)
<i>Afa</i> I	<i>Csp</i> 6 I*, Rsa I	GT▼AC
<i>Afe</i> I	Eco47 III	AGC▼GCT
<i>Afl</i> II	Bst 98 I	C▼TTAAGG
Age I	<i>Pin</i> A I	A▼CCGGT
<i>Aha</i> III	Dra I	TTT▼AAA
<i>Ahd</i> I	Ecl HK I	GACNNN▼NNGTC
Alu I	—	AG▼CT
<i>Alw</i> I	<i>Acc</i> W I	GGATCNNNN▼
Alw26 I ¹	<i>Bsm</i> A I	GTCTC(1/5)
Alw44 I	<i>Apa</i> L I	G▼TGCAC
<i>Aoc</i> I	Bsu 36 I, <i>Cvn</i> I	CC▼TNAGG
Apa I	<i>Bsp</i> 120 I	GGGCC▼C
<i>Apa</i> L I	Alw44 I, <i>Vne</i> I	G▼TGCAC
<i>Apo</i> I	<i>Acs</i> I	(G/A)▼AATT(C/T)
<i>Ase</i> I	Vsp I, <i>Asn</i> I	AT▼TAAT
<i>Asn</i> I	Vsp I, <i>Ase</i> I	AT▼TAAT
<i>Asp</i> I	Tth 111 I	GACN▼NNGTC
<i>Asp</i> E I	<i>Ahd</i> I, <i>Eam</i> 1105 I, Ecl HK I	GACNNN▼NNGTC
<i>Asp</i> 700 I	Xmn I	GAANN▼NNTTC
<i>Asp</i> 718 I	Acc65 I Kpn I*	G▼GTACC GGTAC▼C
<i>Asu</i> I	<i>Sau</i> 96 I, <i>Cfr</i> 13 I	G▼GNCC
<i>Asu</i> II	<i>Csp</i> 45 I, <i>Bst</i> B I	TT▼CGAA
<i>Asu</i> HP I	<i>Hph</i> I	GGTGAN ₈ ▼
Ava I	<i>Ama</i> 87 I, <i>Bco</i> I, <i>Bso</i> B I, <i>Eco</i> 88 I	C▼(C/T)CG(G/A)G
Ava II	Sin I, <i>Eco</i> 47 I, <i>Hgi</i> E I	G▼G(A/T)CC
<i>Axy</i> I	Bsu 36 I	CC▼TNAGG
Bal I	<i>Msc</i> I, <i>Mlu</i> N I	TGG▼CCA
Bam H I	—	G▼GATCC
Ban I	<i>Acc</i> B1, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(T/C)(A/G)CC
Ban II	<i>Eco</i> 24 I	G(A/G)GC(T/C)▼C
<i>Bbe</i> I	— Nar I*	GGCGC▼C GG▼CGCC
<i>Bbr</i> P I	<i>Eco</i> 72 I, <i>Pml</i> I	CAC▼GTG
<i>Bbs</i> I ¹	<i>Bsc</i> 91 I, <i>Bpi</i> I	GAAGAC(2/6)
Bbu I	<i>Pae</i> I, Sph I	GCATG▼C
Bcl I	<i>Bsi</i> Q I, <i>Fba</i> I	T▼GATCA
<i>Bcn</i> I	Nci I	CC▼(C/G)GG
<i>Bfr</i> I	Bst 98 I	C▼TTAAG
Bgl I	—	GCCNNNN▼NGGC
Bgl II	—	A▼GATCT
<i>Bmy</i> I	Bsp1286 I	G(G/A/T)GC(C/A/T)▼C
<i>Bpm</i> I	<i>Gsu</i> I	CTGGAG(16/14)
<i>Bsa</i> H I	Hsp92 I	G(A/G)▼CG(T/C)C
Bsa M I	<i>Bsm</i> I	GAATGC(1/–1)
<i>Bsa</i> O I	<i>Bsh</i> 1285 I, <i>Bsi</i> E I	CG(A/G)(T/C)▼C
<i>Bse</i> A I	Acc III	T▼CCGGA
<i>Bse</i> N I	Bsr S I, <i>Bsr</i> I	ACTGGN (1/–1)
<i>Bse</i> P I	Bss H II, <i>Pau</i> I	G▼CGCGC
<i>Bsh</i> 1285 I	<i>Bsa</i> O I	CG(A/G)(T/C)▼CG
<i>Bsh</i> N I	Ban I, <i>Acc</i> B1 I, <i>Eco</i> 64 I	G▼G(T/C)(A/G)CC
<i>Bsh</i> 1365 I	<i>Bsr</i> BR I	GATNN▼NATC
<i>Bsi</i> E I	<i>Bsa</i> O I	CG(A/G)(T/C)▼CG

Enzyme	Isoschizomer(s)	Recognition Sequence
<i>Bsm</i> I	Bsa M I	GAATGCN▼
<i>Bsm</i> A I ¹	Alw26 I	GTCTC(1/5)
<i>Bso</i> B I	Ava I, <i>Ama</i> 87 I, <i>Bco</i> I, <i>Eco</i> 88 I	C(C/T)CG(G/A)G
<i>Bsp</i> 19 I	Nco I	C▼CATGG
<i>Bsp</i> 68 I	Nru I	TCG▼CGA
<i>Bsp</i> 106 I	Cla I, <i>Bsa</i> D I	AT▼CGAT
<i>Bsp</i> 119 I	Csp45 I, <i>Nsp</i> V, <i>Bst</i> B I	TT▼CGAA
<i>Bsp</i> 120 I	Apa I	G▼GGCCC
<i>Bsp</i> 143 I	Mbo I, Sau3A I, Nde II	▼GATC
<i>Bsp</i> 143 II	Hae II	(A/G)GCGC▼(T/C)
Bsp1286 I	<i>Bmy</i> I, <i>Sdu</i> I	G(G/A/T)GC(C/A/T)▼C
<i>Bsp</i> C I	Pvu I	CGAT▼CG
<i>Bsp</i> D I	Cla I	AT▼CGAT
<i>Bsp</i> E I	Acc III	T▼CCGGA
<i>Bsr</i> I ¹	Bsr S I, <i>Bse</i> N I	ACTGGN(1/–1)
Bsr S I ¹	<i>Bse</i> N I, <i>Bsr</i> I	ACTGGN(1/–1)
Bss H II	<i>Bse</i> P I, <i>Pau</i> I	G▼CGCGC
Bst 98 I	<i>Afl</i> II, <i>Bfr</i> I	C▼TTAAG
<i>Bst</i> B I	Csp45 I, <i>Nsp</i> V, <i>Bsp</i> 119 I	TT▼CGAA
Bst E II	<i>Bst</i> P I, <i>Eco</i> 91 I, <i>Psp</i> E I	G▼GTNACC
<i>Bst</i> N I	Bst O I, <i>Mva</i> I, <i>Eco</i> R II	CC▼(A/T)GG
Bst O I	<i>Bst</i> N I, <i>Eco</i> R II, <i>Mva</i> I	CC▼(A/T)GG
Bst X I	—	CCANNNNN▼NTGG
<i>Bst</i> Y I	Xho II, <i>Mil</i> I	(A/G)▼GATC(T/C)
Bst Z I	Eco52 I, <i>Eag</i> I, <i>Xma</i> III, <i>Ecl</i> X I	C▼GGCCG
<i>Bsu</i> 15 I	Cla I	AT▼CGAT
Bsu 36 I	<i>Cvn</i> I, <i>Aoc</i> I, <i>Eco</i> 81 I	CC▼TNAGG
<i>Bsu</i> R I	Hae III, <i>Pal</i> I	GG▼CC
Cfo I	Hha I <i>Hin</i> 6 I <i>Hin</i> P1 I*	GCG▼C GCG▼C G▼CGC
<i>Cfr</i> 9 I	Xma I Sma I*	C▼CCGGG CCC▼GGG
<i>Cfr</i> 13 I	<i>Sau</i> 96 I	G▼GNCC
<i>Cfr</i> 42 I	Sac II	CCGC▼GG
Cla I	<i>Ban</i> III, <i>Bsp</i> 106 I, <i>Bsp</i> D I, <i>Bsu</i> 15 I	AT▼CGAT
<i>Cpo</i> I	Csp I, <i>Rsr</i> II	CG▼G(A/T)CCG
Csp I	<i>Cpo</i> I, <i>Rsr</i> II	CG▼G(A/T)CCG
<i>Csp</i> 6 I	Rsa I*, <i>Afa</i> I*	GT▼AC
Csp45 I	<i>Bst</i> B I, <i>Nsp</i> V, <i>Bsp</i> 119 I	TT▼CGAA
<i>Cvn</i> I	Bsu 36 I	CC▼TNAGG
Dde I	<i>Bst</i> DE I	C▼TNAG
Dpn I ²	<i>Dpn</i> II*	GmeA▼TC
<i>Dpn</i> II	Mbo I, Sau3A I, Nde II, Dpn I*	▼GATC
Dra I	—	TTT▼AAA
<i>Eag</i> I	Eco52 I, Bst Z I, <i>Ecl</i> X I, <i>Xma</i> III	C▼GGCCG
<i>Eam</i> 1105 I	Ecl HK I, <i>Ahd</i> I, <i>Asp</i> E I	GACNNN▼NNGTC
<i>Ecl</i> 136 II	Ecol CR I Sac I*	GAG▼CTC GAGCT▼C
Ecl HK I	<i>Ahd</i> I, <i>Eam</i> 1105 I, <i>Asp</i> E I	GACNNN▼NNGTC
<i>Ecl</i> X I	Bst Z I, <i>Eag</i> I, Eco52 I, <i>Xma</i> III	C▼GGCCG
<i>Eco</i> 24 I	Ban II, <i>Fri</i> O I	G(AG)GC(TC)▼C
<i>Eco</i> 32 I	Eco R V	GAT▼ATC
<i>Eco</i> 47 I	Ava II, Sin I	G▼G(A/T)CC
Eco47 III	<i>Afe</i> I	AGC▼GCT
Eco52 I	Bst Z I, <i>Xma</i> III, <i>Eag</i> I, <i>Ecl</i> X I	C▼GGCCG
<i>Eco</i> 64 I	Ban I, <i>Bsh</i> N I, <i>Eco</i> 64 I	G▼G(TC)(AG)CC
<i>Eco</i> 81 I	Bsu 36 I	CC▼TNAGG
<i>Eco</i> 88 I	Ava I	C▼(TC)CG(AG)G
<i>Eco</i> 91 I	Bst E II	G▼GTNACC
<i>Eco</i> 105 I	Sna B I	TAC▼CTA
<i>Eco</i> 130 I	Sty I	C▼C(A/T)(T/A)GG
<i>Eco</i> 147 I	Stu I	AGG▼CCT
Ecol CR I	<i>Ecl</i> 136 II	GAG▼CTC
	Sac I*	GAGCT▼C
	<i>Sst</i> I*	GAGCT▼C

Technical Appendix

Isoschizomers (continued). The enzymes in boldface type are available from Promega.

Enzyme	Isoschizomer(s)	Recognition Sequence	Enzyme	Isoschizomer(s)	Recognition Sequence
EcoR I	—	G▼AATTC	Rsa I	<i>Afa I</i>	GT▼AC
<i>EcoR II</i>	BstO I , <i>BstN I</i> , <i>Mva I</i>	CC▼(A/T)GG	<i>Rsr II</i>	Csp I , <i>Cpo I</i>	CG▼G(A/T)CCG
EcoR V	<i>Eco32 I</i>	GAT▼ATC	Sac I	<i>Sst I</i>	GAGCT▼C
<i>EcoT14 I</i>	Sty I	C▼C(A/T)(A/T)GG		<i>EcI136 II*</i> , EcoICR I*	GAG▼CTC
<i>EcoT22 I</i>	Nsi I	ATGCA▼T	Sac II	<i>Sst II</i> , <i>Ksp I</i> , <i>Cfr42 I</i>	CCGC▼GG
<i>Ehe I</i>	Nar I*	GG▼CGCC	Sal I	—	G▼TCGAC
Fok I ²	—	GGATG(9/13)	Sau3A I	Mbo I , Nde II , <i>Dpn II</i>	▼GATC
Hae II	<i>Bsp143 II</i>	(A/G)GCGC▼(T/C)	<i>Sau96 I</i>	<i>Cfr13 I</i>	G▼GNCC
Hae III	<i>BsuR I</i> , <i>Pal I</i>	GG▼CC	Sca I	—	AGT▼ACT
<i>Hap II</i>	Hpa II , Msp I	C▼CGG	<i>Sdu I</i>	Bsp1286 I	G(G/A/T)GC(C/A/T)▼C
<i>HgiE I</i>	<i>Eco47 I</i> , Sin I , Ava II	G▼G(A/T)CC	Sfi I	—	GGCCNNNN▼NGGCC
Hha I	Cfo I	GCG▼C	<i>Sfu I</i>	Csp45 I	TT▼CGAA
	<i>HinP1 I*</i> , <i>Hin6 I*</i>	G▼CGC	Sgf I	—	GCGAT▼CGC
<i>Hin1 I</i>	<i>Acy I</i> , Hsp92 I	G(A/G)▼CG(T/C)C	Sin I	Ava II , <i>Eco47 I</i>	G▼G(A/T)CC
Hinc II	<i>Hind II</i>	GT(T/C)▼(A/G)AC	Sma I	—	CCC▼GGG
<i>Hind II</i>	Hinc II	GT(T/C)▼(A/G)AC		Xma I* , <i>Cfr9 I*</i>	C▼CCGGG
Hind III	—	A▼AGCTT	SnaB I	<i>Eco105 I</i>	TAC▼GTA
Hinf I	—	G▼ANTC	Spe I	<i>AccI</i>	A▼CTAGT
<i>HinP1 I</i>	—	G▼CGC	Sph I	Bbu I , <i>Pae I</i>	GCATG▼C
	Hha I* , Cfo I*	GCG▼C	Ssp I	—	AAT▼ATT
Hpa I	<i>KspA I</i>	GTT▼AAC	<i>Sst I</i>	Sac I	GAGCT▼C
Hpa II ³	Msp I , <i>Hap II</i>	C▼CGG		EcoICR I*	GAG▼CTC
Hsp92 I	<i>Acy I</i> , <i>BsaH I</i> , <i>Hin1 I</i>	G(A/G)▼CG(C/T)	<i>Sst II</i>	Sac II	CCGC▼GG
Hsp92 II	<i>Nla III</i>	CATG▼	Stu I	<i>Aat I</i> , <i>Eco147 I</i>	AGG▼CCT
I-Ppo I	—	CTCTCTTAA▼GGTAGC	Sty I	<i>EcoT14 I</i>	C▼C(A/T)(A/T)GG
<i>Kas I</i>	Nar I*	GG▼CGCC	Taq I	<i>TthHB8 I</i>	T▼CGA
Kpn I	—	GGTAC▼C	Tru9 I	<i>Mse I</i>	T▼TAA
	Acc65 I* , <i>Asp718 I*</i>	G▼GTACC	Tth111 I	<i>Asp I</i>	GACN▼NNGTC
<i>Ksp I</i>	Sac II	CCGC▼GG	<i>TthHB8 I</i>	Taq I	T▼CGA
Mbo I	Sau3A I , Nde II , <i>Dpn II</i>	▼GATC	<i>Van91 I</i>	AccB7 I , <i>PfiM I</i>	CCAN ₄ ▼NTGG
Mbo II ⁴	—	GAAGA(8/7)	<i>Vne I</i>	<i>ApaL I</i> , Alw44 I	G▼TGCAC
<i>Mil I</i>	Xho II	(A/G)▼GATC(T/C)	Vsp I	<i>Ase I</i> , <i>Asn I</i>	AT▼TAAT
Mlu I	—	A▼CGCGT	Xba I	—	T▼CTAGA
<i>MluN I</i>	Bal I , <i>Msc I</i>	TGG▼CCA	Xho I	<i>PaeR7 I</i>	C▼TCGAG
<i>Mro I</i>	Acc III	T▼CCGGA	Xho II	<i>BstY I</i> , <i>Mil I</i>	(A/G)▼GATC(T/C)
<i>Msc I</i>	Bal I , <i>MluN I</i>	TGG▼CCA	Xma I	<i>Cfr9 I</i> , <i>XmaC I</i> , Sma I*	C▼CCGGG CCC▼GGG
<i>Mse I</i>	Tru9 I	T▼TAA	<i>Xma III</i>	Eco52 I , BstZ I , <i>Eag I</i> , <i>EcIX I</i>	C▼GGCCG
Msp I ³	Hpa II , <i>Hap II</i>	C▼CGG	<i>XmaC I</i>	Xma I	C▼CCGGG
MspA1 I	<i>NspB II</i>	C(A/C)G▼C(G/T)G		Sma I*	CCC▼GGG
<i>Mst II</i>	Bsu36 I	CC▼TNAGG	Xmn I	<i>Asp700 I</i>	GAANN▼NNTTC
<i>Mva I</i>	BstO I , <i>EcoR II</i> , <i>BstN I</i>	CC▼(A/T)GG	Key: N = A, C, G or T * = neoschizomer Notes: 1. The locations of cleavage sites falling outside the recognition site are indicated in parentheses. For example, GTCTC(1/5) indicates cleavage at: 5'...GTCTCN▼...3' 3'...CAGAGNNNN▼...5' 2. <i>Dpn I</i> is unique among commercially available restriction enzymes in requiring methylation of a nucleotide (adenine) in its recognition sequence in order to cut. Therefore, <i>Dpn I</i> cannot be substituted for other enzymes recognizing the GATC sequence (e.g., <i>Mbo I</i> and <i>Sau3A I</i>). 3. Although <i>Hpa II</i> and <i>Msp I</i> recognize the same nucleotide sequence, <i>Hpa II</i> is sensitive to methylation of either cytosine in its recognition sequence, while <i>Msp I</i> is sensitive only to methylation of the external cytosine. These enzymes may not be interchanged for all applications.		
Nae I	NgoM IV	G▼CCGGC			
Nar I	—	GG▼CGCC	Reference Roberts, R.J. (1991) <i>Nucl. Acids Res.</i> 19 (supp), 2077–109.		
	<i>Ehe I*</i>	GGC▼GCC			
	<i>Kas I*</i>	G▼GCGCC			
	<i>Bbe I*</i>	GGCGC▼C			
Nci I	<i>Bcn I</i>	CC▼(C/G)GG			
Nco I	<i>Bsp19 I</i>	C▼CATGG			
Nde I	—	CA▼TATG			
Nde II	Mbo I , Sau3A I , <i>Dpn II</i>	▼GATC			
NgoM IV	Nae I	G▼CCGGC			
Nhe I	—	G▼CTAGC			
<i>Nla III</i>	Hsp92 II	CATG▼			
Not I	—	GC▼GGCCGC			
Nru I	<i>Bsp68 I</i>	TCG▼CGA			
Nsi I	<i>EcoT22 I</i> , <i>Mph1103 I</i>	ATGCA▼T			
<i>Nsp V</i>	Csp45 I , <i>BstB I</i> , <i>Bsp119 I</i>	TT▼CGAA			
<i>NspB II</i>	MspA1 I	C(A/C)G▼C(G/T)G			
<i>Pae I</i>	Bbu I , Sph I	GCATG▼C			
<i>PaeR7 I</i>	Xho I	C▼TCGAG			
<i>Pal I</i>	Hae III , <i>BsuR I</i>	GG▼CC			
<i>PfiM I</i>	AccB7 I , <i>Vau91 I</i>	CCAN ₄ ▼NTGG			
<i>PinA I</i>	Age I	A▼CCGGT			
Pst I	—	CTGCA▼G			
Pvu I	<i>BspC I</i>	CGAT▼CG			
Pvu II	—	CAG▼CTG			

Key:

N = A, C, G or T
* = neoschizomer

Notes:

- The locations of cleavage sites falling outside the recognition site are indicated in parentheses. For example, GTCTC(1/5) indicates cleavage at:
5'...GTCTCN▼...3'
3'...CAGAGNNNN▼...5'
- Dpn I* is unique among commercially available restriction enzymes in requiring methylation of a nucleotide (adenine) in its recognition sequence in order to cut. Therefore, *Dpn I* cannot be substituted for other enzymes recognizing the GATC sequence (e.g., *Mbo I* and *Sau3A I*).
- Although *Hpa II* and *Msp I* recognize the same nucleotide sequence, *Hpa II* is sensitive to methylation of either cytosine in its recognition sequence, while *Msp I* is sensitive only to methylation of the external cytosine. These enzymes may not be interchanged for all applications.

Reference

Roberts, R.J. (1991) *Nucl. Acids Res.* **19** (supp), 2077–109.

Technical Appendix

Compatible Ends.

Promega Restriction Enzymes That Generate 5' Overhangs

Overhang	Definite Compatible Ends	Possible Compatible Ends
5'-N		<i>Tth</i> 111 I
5'-S		<i>Nci</i> I
5'-W		<i>Bst</i> O I
5'-AT	<i>Acc</i> I	
5'-CG	<i>Nar</i> I, <i>Msp</i> I, <i>Hsp</i> 92 I, <i>Taq</i> I, <i>Cla</i> I, <i>Csp</i> 45 I, <i>Hpa</i> II	
5'-GN		<i>Bsr</i> S I
5'-MK		<i>Acc</i> I
5'-TA	<i>Vsp</i> I, <i>Nde</i> I, <i>Tru</i> 9 I	
5'-ANT		<i>Hint</i> I
5'-GNC		<i>Sau</i> 96 I
5'-GWC		<i>Ava</i> II, <i>Csp</i> I, <i>Sin</i> I
5'-TNA		<i>Dde</i> I, <i>Bsu</i> 36 I
5'-AATT	<i>Eco</i> R I	
5'-AGCT	<i>Hind</i> III	
5'-CATG	<i>Nco</i> I	<i>Sty</i> I
5'-CCGG	<i>Age</i> I, <i>Xma</i> I, <i>Acc</i> III, <i>Ngo</i> M IV	<i>Ava</i> I
5'-CGCG	<i>Mlu</i> I, <i>Bss</i> H I	
5'-CTAG	<i>Spe</i> I, <i>Nhe</i> I, <i>Xba</i> I	<i>Sty</i> I
5'-CWWG		<i>Sty</i> I
5'-GATC	<i>Mbo</i> I, <i>Sau</i> 3A I, <i>Bam</i> H I, <i>Bgl</i> II, <i>Xho</i> II, <i>Bcl</i> I, <i>Nde</i> II	
5'-GCGC	<i>Ban</i> I	
5'-GGCC	<i>Not</i> I, <i>Bst</i> Z I, <i>Eco</i> 52 I	
5'-GTAC	<i>Acc</i> 65 I	<i>Ban</i> I
5'-GTNAC		<i>Bst</i> E II
5'-GYRC		<i>Ban</i> I
5'-TCGA	<i>Sal</i> I, <i>Xho</i> I	<i>Ava</i> I
5'-TGCA	<i>Alw</i> 44 I	
5'-TTAA	<i>Bst</i> 98 I	
5'-YCGR		<i>Ava</i> I

Promega Restriction Enzymes That Generate 3' Overhangs

Overhang	Definite Compatible Ends	Possible Compatible Ends
N-3'		<i>Ecl</i> HK I
AT-3'	<i>Sgf</i> I, <i>Pvu</i> I	
CG-3'	<i>Cfo</i> I, <i>Hha</i> I	
CN-3'		<i>Bsa</i> M I
GC-3'	<i>Sac</i> II	<i>Bsa</i> O I
NNN-3'		<i>Acc</i> B7 I, <i>Bgl</i> I, <i>Sfi</i> I
ACGT-3'	<i>Aat</i> II	
AGCT-3'	<i>Sac</i> I	<i>Ban</i> II, <i>Bsp</i> 1286 I
CATG-3'	<i>Hsp</i> 92 II, <i>Sph</i> I, <i>Bbu</i> I	
DGCH-3'		<i>Bsp</i> 1286 I
GCGC-3'	<i>Hae</i> II	
GGCC-3'	<i>Apa</i> I	<i>Ban</i> II, <i>Bsp</i> 1286 I
GTAC-3'	<i>Kpn</i> I	
NNNN-3'		<i>Bst</i> X I
RGCY-3'		<i>Ban</i> II, <i>Bsp</i> 1286 I
TGCA-3'	<i>Nsi</i> I, <i>Pst</i> I	<i>Bsp</i> 1286 I
TTAA-3'	<i>I-Ppo</i> I	

Key:

D = A or G or T
 H = A or C or T
 K = G or T
 M = A or C
 N = A or C or G or T
 R = A or G
 S = C or G
 W = A or T
 Y = C or T

Technical Appendix

Site-Specific Methylation Sensitivity of Promega Restriction Enzymes.

This table lists the sensitivities of several Promega restriction enzymes to site-specific methylation at *dam*, *dcm*, CpG and CpNpG sites (p = phosphoryl group). These four modifications are frequently found in DNA of bacteria, eukaryotes or their viruses. Many strains of *E. coli* contain the site-specific *dam* and *dcm* DNA methylases. Higher eukaryotes contain the site-specific CpG and CpNpG DNA methylases. In mammalian genomes, methylation occurs mainly at the CG dinucleotide. In plant genomes, methylation may occur at both the CG and CNG sequences.

Prokaryotic Methylation

dcm Cytosine methylase mutation—methylates the C5 position of the internal cytosine residue in the sequence 5'...CCTGG...3'.

dam Adenine methylase mutation—methylates the N6 position of the adenine residue in the sequence 5'...GATC...3'.

For further information regarding site-specific methylation, refer to McClelland, M., Nelson, M. and Raschke, E. (1994) *Nucl. Acids Res.* **22**, 3640–59.

Key:

- s = sensitive to this methylation
- i = insensitive to this methylation
- s(ol) = overlapping (sensitive when restriction site overlaps methylation sequence)
- n/a = information not available

Eukaryotic Methylation

CpG Methylates the C5 position of the cytosine residue in the dinucleotide recognition sequence 5'...CG...3'.

CpNpGp Methylates the C5 position of the cytosine residue in the trinucleotide recognition sequence 5'...CNG...3' (N = any base).

Enzyme	Recognition Sequence	<i>dam</i>	<i>dcm</i>	CpG	CpNpG	Enzyme	Recognition Sequence	<i>dam</i>	<i>dcm</i>	CpG	CpNpG
<i>Aat</i> II	GACGTC	i	i	s	i	<i>Kpn</i> I	GGTACC	i	i	i	i
<i>Acc</i> B7 I	CCANNNNTGG	i	s(ol)	i	i	<i>Mbo</i> II	GAAGA(8/7)	s(ol)	i	i	i
<i>Acc</i> III	TCCGGA	s(ol)	i	i	i	<i>Mlu</i> I	ACGCGT	i	i	s	i
<i>Acc</i> 65 I	GGTACC	i	s(ol)	i	i	<i>Msp</i> I	CCGG	i	i	i	s
<i>Apa</i> I	GGGCCC	i	s(ol)	s(ol)	i	<i>Nae</i> I	GCCGGC	i	i	s	s
<i>Ava</i> I	CYCGRG	i	i	s	i	<i>Nar</i> I	GGCGCC	i	i	s	i
<i>Ava</i> II	GGWCC	i	s(ol)	s(ol)	s(ol)	<i>Nde</i> II	GATC	s	i	i	i
<i>Bal</i> I	TGGCCA	i	s(ol)	i	s(ol)	<i>Ngo</i> M IV	GCCGGC	i	i	s	s
<i>Bam</i> HI	GGATCC	i	i	i	s(ol)	<i>Nhe</i> I	GCTAGC	i	i	s(ol)	s(ol)
<i>Ban</i> II	GRGCTC	i	i	i	i	<i>Not</i> I	GCGGCCGC	i	i	s	s
<i>Bbu</i> I	GCATGC	i	i	i	i	<i>Nru</i> I	TCGCGA	s(ol)	i	s	i
<i>Bcl</i> I	TGATCA	s	i	i	i	<i>Pst</i> I	CTGCAG	i	i	i	s
<i>Bgl</i> I	GCCNNNNNGGC	i	i	s(ol)	s(ol)	<i>Pvu</i> I	CGATCG	i	i	s	s(ol)
<i>Bgl</i> II	AGATCT	i	i	i	s(ol)	<i>Pvu</i> II	CAGCTG	i	i	i	s
<i>Bsp</i> 1286 I	GDGCHC	i	i	i	i	<i>Sac</i> I	GAGCTC	i	i	i	i
<i>Bss</i> HI	GCGCGC	i	i	s	i	<i>Sac</i> II	CCGCGG	i	i	s	s
<i>Bst</i> E II	GGTNACC	i	i	i	i	<i>Sal</i> I	GTCGAC	i	i	s	n/a
<i>Bst</i> O I	CCWGG	i	i	i	n/a	<i>Sau</i> 3A I	GATC	i	i	s(ol)	s(ol)
<i>Bst</i> X I	CCANNNNNNTGG	i	i	i	i	<i>Sau</i> 96 I	GGNCC	i	s(ol)	s(ol)	s(ol)
<i>Bst</i> Z I	CGGCCG	i	i	s(ol)	s(ol)	<i>Sca</i> I	AGTACT	i	i	i	i
<i>Cfo</i> I	GCGC	i	i	s	n/a	<i>Sfi</i> I	GGCCNNNNNGGCC	i	s(ol)	s(ol)	s(ol)
<i>Cla</i> I	ATCGAT	s(ol)	i	s	i	<i>Sgf</i> I	GCGATCGC	i	i	s	n/a
<i>Csp</i> I	CGGWCCG	i	i	i	s	<i>Sin</i> I	GGWCC	i	i	i	s(ol)
<i>Csp</i> 45 I	TTCGAA	i	i	s	i	<i>Sma</i> I	CCCGGG	i	i	s	s
<i>Dde</i> I	CTNAG	i	i	i	s(ol)	<i>Sna</i> B I	TACGTA	i	i	s	i
<i>Eco</i> 47 III	AGCGCT	i	i	s	i	<i>Sph</i> I	GCATGC	i	i	i	i
<i>Eco</i> 52 I	CGGCCG	i	i	s	i	<i>Stu</i> I	AGGCCT	i	s(ol)	i	s(ol)
<i>Eco</i> R I	GAATTC	i	i	s(ol)	i	<i>Taq</i> I	TCGA	s(ol)	i	i	i
<i>Fok</i> I	GGATC	i	i	i	i	<i>Xba</i> I	TCTAGA	s(ol)	i	i	i
<i>Hae</i> III	GGCC	i	i	i	s(ol)	<i>Xho</i> I	CTCGAG	i	i	s	i
<i>Hha</i> I	GCGC	i	i	s	s(ol)	<i>Xho</i> II	RGATCY	i	i	i	s(ol)
<i>Hinc</i> II	GTYRAC	i	i	i	i	<i>Xma</i> I	CCCGGG	i	i	i	n/a
<i>Hind</i> III	AAGCTT	i	i	i	i	<i>Xmn</i> I	GAANNNN	i	i	n/a	n/a
<i>Hpa</i> II	CCGG	i	i	s	s						

Technical Appendix

Restriction Enzyme Buffer Composition.

Buffer	pH (at 37°C)	Tris-HCl (mM)	MgCl ₂ (mM)	NaCl (mM)	KCl (mM)	DTT (mM)
A	7.5	6	6	6	—	1
B	7.5	6	6	50	—	1
C	7.9	10	10	50	—	1
D	7.9	6	6	150	—	1
E	7.5	6	6	100	—	1
F	8.5	10	10	100	—	1
G	8.2	50	5	—	—	—
H	7.5	90	10	50	—	—
J	7.5	10	7	—	50	1
K	7.4	10	10	—	150	—
L	9.0	10	3	100	—	—

MULTI-CORE™ Buffer (1X) = 25mM Tris-acetate (pH 7.5 at 37°C), 100mM potassium acetate, 10mM magnesium acetate, 1mM DTT.

Notes:

- For each 10°C rise in temperature between 0°C and 25°C, the pH of Tris buffers decreases 0.31 pH units.
- For each 10°C rise in temperature between 25°C and 37°C, the pH of Tris buffers decreases 0.25 pH units.
- All of Promega's Restriction enzymes are supplied with 10mg/ml Acetylated BSA. Although BSA is not absolutely required for activity, it has been shown to enhance activity of many restriction enzymes. We recommend adding BSA to all restriction digests at a final concentration of 0.1mg/ml.

Copy Number of Commonly Used Plasmids.

Plasmid	Plasmid Size (approx.)	Origin of Replication*	Copy Number	**Yield per ml of Culture	Reference
pGEM®	2,700bp	mutated pMB1	300–700	1.8–4.1µg	1
pUC	2,700bp	mutated pMB1	500–700	2.9–4.1µg	1
pBR322	4,400bp	pMB1	>25	>0.23µg	2
ColE1	4,500bp	ColE1	>15	>0.15µg	3
pACYC	4,000bp	p15A	~10	~0.09µg	4
pSC101	9,000bp	pSC101	~6	~0.12µg	5
pGL Series	5,000bp	mutated pMB1	300–700	3.3–7.6µg	1
pRL Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
phRL Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
phRG Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
pGEM®-T/ T easy	3,000bp	mutated pMB1	300–700	2.0–4.6µg	1
psiLentGene™ Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
psiCHECK™ 1/2	3,500bp	mutated pMB1	300–700	2.7–5.3µg	1
psiSTRIKE™ Series	4,000bp	mutated pMB1	300–700	2.7–6.0µg	1
pALTER®-1/Ex1	5,800bp	pMB1	>25	>0.3µg	3
pALTER®-Ex2	5,800bp	p15A	~10	~0.13µg	4
pSP	2,500bp	mutated pMB1	300–700	1.6–3.8µg	1
pCI, pSI	3,600bp	mutated pMB1	300–700	2.4–5.5µg	1

* Plasmids carrying the pMB1, mutated pMB1 and ColE1 belong to the same incompatibility group, so they are not compatible with one another, but they are fully compatible with those carrying p15A and pSC101 replicons.

** Theoretical plasmid yields were calculated from the reported copy number and size of each plasmid assuming 2.0×10^9 cells per milliliter of culture grown for 16 hours at 37°C.

References

- Summerton, J., Atkins, T. and Bestwick, R. (1983) *Anal. Biochem.* **133**, 79–84.
- Holmes, D.S. and Quigley, M. (1981) *Anal. Biochem.* **114**, 193–7.
- Jansz, H.S., Pouwels, P.H. and Schiphorst, J. (1966) *Biochem. Biophys. Acta* **123**, 626.
- Birnboim, H.C. and Doly, J. (1979) *Nucl. Acids Res.* **7**, 1513–23.
- Birnboim, H.C. (1983) *Meth. Enzymol.* **100**, 243–55.

Star Activity.

Restriction enzymes, under nonstandard conditions, can demonstrate the ability to cleave DNA at sequences different from their defined recognition sites. The term "star activity" has been given to this nonsequence-specific cleavage of DNA under nonoptimal reaction conditions. The most common types of altered activity are single-base substitutions, truncation of the outer bases in the recognition sequence and single-strand nicking (1). In general, star activity is not a concern if restriction endonucleases are used in the recommended buffers at the appropriate temperatures. Star activity is evident with a number of restriction enzymes when the following parameters are altered in the reaction environment (2):

- High enzyme concentration (generally >100 units/µg).
- High glycerol content (>5% v/v).
- Substitution of Mn²⁺ for Mg²⁺ (or substitution of other divalent cations).
- Low salt concentration (generally <25mM).
- Extremes of pH, especially pH>8.0.
- Presence of DMSO, ethanol or other organic solvents.

References

- Barany, F. (1988) *Gene* **65**, 149–65.
- Brown, T.A., Hames, B.D. and Rickwood, D. (1991) In: *Molecular Biology Lab Fax*, BIOS Scientific Publishers Limited, Oxford, United Kingdom.

Technical Appendix

Genotypes of Frequently Used Bacterial Strains.

All genes in the bacterium are presumed to be in the wildtype state, except for those listed, which are mutant alleles carried by that bacterium. Genes listed on the F' episome, however, represent wildtype alleles unless specified otherwise. Strains are λ^- unless specified otherwise. *Strains available from Promega as competent cells are indicated by an asterisk. Strains shown in **bold** are available from Promega as glycerol freezer stocks.

Strain	Genotype
BL21(DE3)	F ⁻ , <i>ompT</i> , <i>hsdS_B</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>dcm</i> , <i>gal</i> , λ (DE3)
*BL21(DE3)pLysS	F ⁻ , <i>ompT</i> , <i>hsdS_B</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>dcm</i> , <i>gal</i> , λ (DE3), pLysS (Cm ^r)
*BMH 71-18 mutS	<i>thi</i> , <i>supE</i> , Δ (<i>lac-proAB</i>), [<i>mutS</i> ::Tn10(<i>tet^r</i>)] [F', <i>tra</i> D36, <i>proAB</i> , <i>laqI^a</i> Δ M15]
C600 (1)	<i>thi-1</i> , <i>thr-1</i> , <i>leuB6</i> , <i>lacY1</i> , <i>tonA21</i> , <i>supE44</i>
C600 <i>hfl</i> (1)	<i>thi-1</i> , <i>thr-1</i> , <i>leuB6</i> , <i>lacY1</i> , <i>tonA21</i> , <i>supE44</i> , <i>hflA150</i> ::Tn10(<i>tet^r</i>)
DH1 (2)	<i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i>
DH10B	F ⁻ , <i>mcrA</i> Δ (<i>mrr-hsdRMS-mcrBC</i>) ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>endA1</i> , <i>araD139</i> , Δ (<i>ara, leu</i>)7697, <i>galU</i> , <i>galK</i> , λ^- , <i>rpsL</i> (<i>str^r</i>), <i>nupG</i>
DH5 α TM	ϕ 80d <i>lacZ</i> Δ M15, <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i> , <i>deoR</i> , Δ (<i>lacZYA-argF</i>) U169, <i>phoA</i>
DM1 (3)	F', <i>dam-13</i> ::Tn9(Cm ^r) <i>dcm</i> , <i>mcrB</i> , <i>hsdR-M^r</i> , <i>gal1</i> , <i>gal2</i> , <i>ara-</i> , <i>lac-</i> , <i>thr-</i> , <i>leu-</i> , <i>ton^R</i> , <i>tsx^R</i> , <i>Su^o</i>
ES1301 mutS	<i>lacZ53</i> , <i>thyA36</i> , <i>rha-5</i> , <i>metB1</i> , <i>deoC</i> , IN(<i>rrnD-rrnE</i>), [<i>mutS201</i> ::Tn5]
*HB101 (4)	<i>thi-1</i> , <i>hsdS20</i> (<i>r_B⁻</i> , <i>m_B⁻</i>), <i>supE44</i> , <i>recA13</i> , <i>ara-14</i> , <i>leuB6</i> , <i>proA2</i> , <i>lacY1</i> , <i>galK2</i> , <i>rpsL20</i> (<i>str^r</i>), <i>xyl-5</i> , <i>mtl-1</i>
JM101 (5)	<i>supE</i> , <i>thi</i> , Δ (<i>lac-proAB</i>), F' (<i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15)
*JM109 (5)	<i>endA1</i> , <i>recA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>relA1</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15]
JM109(DE3) (5)	<i>endA1</i> , <i>recA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>relA1</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15], λ (DE3)
JM110 (5)	<i>rpsL</i> (<i>str^r</i>), <i>thr</i> , <i>leu</i> , <i>thi</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>lacY</i> , <i>galK</i> , <i>galT</i> , <i>ara</i> , <i>tonA</i> , <i>tsx</i> , <i>dam</i> , <i>dcm</i> , <i>supE44</i> , Δ (<i>lac-proAB</i>), [F', <i>traD36</i> , <i>proAB</i> , <i>lacI^a</i> Δ M15]
KW251	<i>supE44</i> , <i>galK2</i> , <i>galT22</i> , <i>metB1</i> , <i>hsdR2</i> , <i>mcrB1</i> , <i>mcrA</i> , [<i>argA81</i> ::Tn10(<i>tet^r</i>)], <i>recD1014</i>
LE392 (6)	<i>hsdR514</i> , (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>supF58</i> , <i>lacY1</i> or Δ (<i>lacIZY</i>)6, <i>galK2</i> , <i>galT22</i> , <i>metB1</i> , <i>trpR55</i>
NM522 (7)	<i>supE</i> , <i>thi</i> , Δ (<i>lac-proAB</i>), Δ <i>hsd5</i> (<i>r_K⁻</i> , <i>m_K⁻</i>), [F', <i>proAB</i> , <i>lacI^a</i> Δ M15]
NM538 (8)	<i>supF</i> , <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>trpR</i> , <i>lacY</i>
NM539 (8)	<i>supF</i> , <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>lacY</i> , (P2)
*Select96 TM	<i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>recA1</i> , <i>araD139</i> (<i>ara-leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> , <i>endA1</i> , <i>nupG</i>
Stb12 TM	F ⁻ , <i>mcrA</i> , Δ (<i>mcrBC-hsdRMS-mrr</i>), <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>supE44</i> , <i>relA1</i> , λ^- , Δ (<i>lac-proAB</i>)
Stb14 TM	<i>mcrA</i> , Δ (<i>mcrBC-hsdRMS-mrr</i>), <i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>supE44</i> , <i>relA1</i> , λ^- , Δ (<i>lac-proAB</i>), <i>gal</i> , F' { <i>proAB⁺</i> , <i>lacI^a</i> , Δ M15, <i>Tn10</i> (<i>tet^R</i>)}
SURE [®]	<i>e14-</i> , (<i>mcrA-</i>) Δ (<i>mcrCB-hsdSMR-mrr</i>)171, <i>endA1</i> , <i>supE44</i> , <i>thi-1</i> , <i>gyrA96</i> , <i>relA1</i> , <i>lac</i> , <i>recB</i> , <i>recJ</i> , <i>sbcC</i> , <i>umuC</i> ::Tn5 (<i>kan^r</i>), <i>uvrC</i> , [F' <i>proAB</i> , <i>lacI^a</i> Δ M15::Tn10 (<i>tet^r</i>)]
TOP10	F ⁻ , <i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>araD139</i> , Δ (<i>ara, leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> (<i>str^R</i>), <i>endA1</i> , <i>nupG</i>
TOP10F'	F' { <i>lacI^a</i> Tn10 (<i>tet^R</i>)}, <i>mcrA</i> , Δ (<i>mrr-hsdRMS-mcrBC</i>), ϕ 80 <i>lacZ</i> Δ M15, Δ <i>lacX74</i> , <i>deoR</i> , <i>recA1</i> , <i>araD139</i> , Δ (<i>ara-leu</i>)7697, <i>galU</i> , <i>galK</i> , <i>rpsL</i> (<i>str^r</i>), <i>endA1</i> , <i>nupG</i>
XL1-Blue	<i>recA1</i> , <i>endA1</i> , <i>gyrA96</i> , <i>thi-1</i> , <i>hsdR17</i> (<i>r_K⁻</i> , <i>m_K⁺</i>), <i>supE44</i> , <i>relA1</i> , <i>lac</i> , [F', <i>proAB</i> , <i>lacI^a</i> Δ M15::Tn10(<i>tet^r</i>)]
Y1089 (9)	Δ (<i>lacU169</i>), <i>proA⁺</i> , Δ (<i>lon</i>), <i>araD139</i> , <i>strA</i> , <i>hflA150</i> , [<i>chr</i> ::Tn10(<i>tet^r</i>)], (pMC9)
Y1090 (9)	Δ (<i>lacU169</i>), <i>proA⁺</i> , Δ (<i>lon</i>), <i>araD139</i> , <i>strA</i> , <i>supF</i> , <i>rpsL</i> (<i>str^r</i>), [<i>trpC22</i> ::Tn10 (<i>tet^r</i>)], (pMC9), <i>hsdR</i> (<i>r_K⁻</i> , <i>m_K⁺</i>)

Miscellaneous

F' Host contains an F' episome with the stated features

λ (DE3) Bacteriophage λ carrying the gene for T7 RNA polymerase is integrated into the host genome.

pMC9 is pBR322 with *lacI^a* inserted and confers amp and tet resistance.

References

- Jendrisak, J., Young, R.A. and Engel, J. (1987) In: *Guide to Molecular Cloning Techniques*, Berger, S. and Kimmel, A., eds., Academic Press, San Diego, CA.
- Hanahan, D. (1983) *J. Mol. Biol.* **166**, 557–80.
- Lorow-Murray, D. and Bloom, F. (1991) *Focus* **13**, 20.
- Lacks, S. and Greenberg, J.B. (1977) *J. Mol. Biol.* **114**, 153–60.
- Yanisch-Perron, C., Vieira, J. and Messing, J. (1985) *Gene* **33**, 103–19.
- Murray, N. *et al.* (1977) *Mol. Gen. Genet.* **150**, 53–61.
- Gough, J. and Murray, N. (1983) *J. Mol. Bio.* **166**, 1–19.
- Frischauf, A. *et al.* (1983) *J. Mol. Biol.* **170**, 827–42.
- Huynh, T., Young, R.A. and Davis, R. (1985) In: *DNA Cloning*, Vol. 1, Glover, D., ed., IRL Press Ltd., Oxford, UK.

Technical Appendix

Genetic Markers in *E. coli*.

Symbol	Description	Effect of Mutation
<i>ara-14</i>	Mutation in arabinose metabolism	Blocks arabinose catabolism.
<i>araD</i>	L-ribulose phosphate 4-epimerase mutation; part of an inducible operon <i>araBAD</i> repressed by L-arabinose	Blocks arabinose catabolism.
<i>argA</i>	N-Acetylglutamate synthase mutation; inhibited by the presence of arginine	Arginine required from growth in minimal media.
<i>cycA</i>	Involved in D-alanine, glycine, D-serine and D-cycloserine transport, and an L-alanine carrier	Mutants cannot use D-alanine as a carbon source.
<i>dam</i>	DNA adenine methylase mutation	Blocks methylation of adenine residues in the sequence 5'...G ^m ATC...3'.
<i>dapD</i>	Succinyl-diaminopimelate aminotransferase mutation	Mutant reflects impaired synthesis of succinyl CoA and needs to be supplemented with succinate or lysine + methionine.
<i>dcm</i>	DNA cytosine methylase mutation	Blocks methylation of cytosine in the sequence 5'...C ^m CAGG...3' or 5'...C ^m CTGG...3'.
<i>deoC</i>	Deoxyribose-phosphate aldolase mutation	
<i>deoR</i>	Regulatory gene mutation allowing constitutive expression of genes for deoxyribose synthesis	Allows efficient propagation of large plasmids.
<i>dut1</i>	Mutation of deoxyuridine triphosphatase, which catalyzes dUTP the conversion to dUMP and PPi	Mutants are impaired in conversion of dUTP to dUMP, leading to higher dUTP pools that can lead to misincorporation of uracil instead of thymidine. Stable incorporation of dUTP needs mutation in <i>ung</i> gene.
<i>endA1</i>	DNA-specific endonuclease I mutation	Improves quality of plasmid DNA isolations.
<i>galE</i>	Part of the <i>galETK</i> operon that encodes UDP galactose-4-epimerase	Mutant is more resistant to bacteriophage P1 infection.
<i>galK</i>	Galactokinase mutation	Blocks catabolism of galactose.
<i>galT</i>	Galactose-1-phosphate uridylyltransferase mutation	Blocks catabolism of galactose.
<i>gyrA96</i>	DNA gyrase mutation	Confers resistance to nalidixic acid.
<i>hflA150</i>	Protease mutation that leads to stabilization of <i>cll</i> gene products	Leads to high frequency of lysogeny by λ phages (1).
<i>hflB</i>	Gene encodes a possible protease component	Mutations lead to high frequency of bacteriophage lambda lysogenization.
<i>hsdR</i> (r_K^- , m_K^+)	Host DNA restriction and methylation system mutation: Restriction minus, modification positive for the <i>E. coli</i> K strain methylation system	Allows cloning without cleavage of transformed DNA by endogenous restriction endonucleases. DNA prepared from this strain can be used to transform r_K^+ <i>E. coli</i> strains.
<i>hsdS20</i> (r_B^- , m_B^-)	Mutation of specificity determinant for host DNA restriction and methylation system. Restriction minus, modification minus for the <i>E. coli</i> B strain methylation system	Allows cloning without cleavage of transformed DNA by endogenous restriction endonucleases. DNA prepared from this strain is unmethylated by the <i>hsdS20</i> methylases.
<i>lacI^q</i>	Overproduction of the <i>lac</i> repressor protein	Leads to high levels of the <i>lac</i> repressor protein, inhibiting transcription from the <i>lac</i> promoter.
<i>lacY</i>	Galactoside permease mutation	Blocks lactose utilization.
<i>lacZΔM15</i>	Partial deletion of β -D-galactosidase gene	Allows complementation of β -galactosidase activity by α -complementation sequence in pGEM [®] -Z Vectors. Allows blue/white selection for recombinant colonies when plated on X-Gal.
<i>leuB</i>	β -isopropylmalate dehydrogenase mutation	Requires leucine for growth on minimal media.
Δ (<i>lon</i>)	Deletion of <i>lon</i> protease	Reduces proteolysis of expressed proteins.
<i>LysS</i>	pLysS plasmid is integrated into the host genome	Strains carrying this plasmid will be tet resistant and produce T7 lysozyme, a natural inhibitor of T7 RNA polymerase, thus lowering background transcription of sequences under the control of the T7 RNA polymerase promoter (2).
<i>mcrA</i>	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'...G ^m CGC...3'.
<i>mcrB</i>	Mutation in methylcytosine restriction system	Blocks restriction of DNA methylated at the sequence 5'...AG ^m CT...3'.
<i>metB</i>	Cystathionine γ -synthase mutation	Requires methionine for growth on minimal media.
<i>metC</i>	Cystathionine beta-lyase mutation; involved in methionine biosynthesis	Methionine required from growth in minimal media.
<i>mtl</i>	Mutation in mannitol metabolism	Blocks catabolism of mannitol.
<i>mutS</i>	Methyl-directed mismatch repair mutation	Prevents repair of the newly synthesized, unmethylated strand.
<i>ompT</i>	Mutation of protease VII, an outer membrane protein	Reduces proteolysis of expressed proteins.
P2	P2 bacteriophage lysogen present in host	λ phages containing the <i>red</i> and <i>gam</i> genes of λ are growth inhibited by P2 lysogens (3).
<i>proA</i>	γ -glutamyl phosphate reductase mutation	<i>proA/argD</i> mutant will not block proline synthesis, but will be repressed by arginine. Mutants excrete proline on minimal media and are resistant to proline analogs. <i>proA/argD/argR</i> triple mutant grows slowly on minimal media + arginine.
<i>proAB</i>	Mutations in proline metabolism	Requires proline for growth in minimal media.

Technical Appendix

Genetic Markers in *E. coli* (continued).

Symbol	Description	Effect of Mutation
<i>recA1</i> , <i>recA13</i>	Mutation in recombination	Minimizes recombination of introduced DNA with host DNA, increasing stability of inserts. Inserts are more stable in <i>recA1</i> than <i>recA13</i> hosts.
<i>recB</i> , <i>recC</i> <i>recD</i>	Exonuclease V mutations The Rec BCD trimer (exonuclease V) progressively degrades ssDNA and dsDNA in an ATP-dependent manner to form oligonucleotides; implicated in homologous recombination	Reduces general recombination and affects repair of radiation damage. Allows easier propagation of sequences with inverted repeats.
<i>recF</i>	Recombination and repair mutation	Mutant cannot repair daughter strand gaps (post-replicative repair).
<i>relA</i>	ppGpp synthetase I mutation, a novel nucleotide guanosine 5'-diphosphate-3'-diphosphate produced in response to starvation by <i>relA</i> ribosomal protein sensing uncharged tRNA	Allows RNA synthesis in the absence of protein synthesis.
<i>rha</i>	Utilization of L-rhamnose, a methylpentose	Blocks rhamnose catabolism.
<i>rpsL</i>	Mutation in subunit S12 of 30S ribosome	Confers resistance to streptomycin.
<i>sbcB</i>	Exonuclease I mutation	Allows general recombination in <i>recBC</i> mutant strains.
<i>strA</i>	Mutant alters ribosome protein S12	Confers resistance to streptomycin
<i>supB</i> , <i>supC</i> , <i>supG</i> , <i>supL</i> , <i>supM</i> , <i>supN</i> , <i>supO</i>	Suppressor mutations	Suppresses ochre (UAA) and amber (UAG) mutations.
<i>supD</i> , <i>supE</i> , <i>supF</i>	Suppressor mutations	Suppresses amber (UAG) mutations.
<i>thi-1</i>	Mutation in thiamine metabolism	Thiamine required for growth in minimal media.
<i>thr</i>	Threonine biosynthesis mutation	Mutants are obligate threonine auxotrophs.
<i>thyA</i>	Thymidylate synthase; dTTP biosynthesis	Mutants are obligate thymidine auxotrophs.
Tn5	Transposon	Encodes resistance to kanamycin.
Tn10	Transposon	Encodes resistance to tetracycline.
<i>tonA</i>	Mutation in outer membrane protein	Confers resistance to bacteriophage T1.
<i>traD36</i>	Transfer factor mutation	Prevents transfer of F' episome.
<i>trpC</i>	Phosphoribosyl anthranilate isomerase mutation; part of tryptophan biosynthesis pathway	
<i>trpR</i>	<i>trpR</i> aporepressor; regulates the biosynthesis of tryptophan and its transport	
<i>tsx</i>	T6 and colicin K phage receptor; outer membrane protein involved in specific diffusion of nucleosides; transports the antibiotic albicidin	Resistant to bacteriophage T6 and colicin K.
<i>ung1</i>	Uracil-DNA N-glycosylase	Allows uracil to exist in plasmid DNA.
<i>xyl-5</i>	Mutation in xylose metabolism	Blocks catabolism of xylose.

References

1. Hoyt, M.A. *et al.* (1982) *Cell* **31**, 565–73.
2. Studier, F.W. (1991) *J. Mol. Biol.* **219**, 37–44.
3. Kaiser, K. and Murray, N. (1985) In: *DNA Cloning*, Vol. 1, Glover, D., ed., IRL Press Ltd., Oxford, UK.
4. Neidhardt, F. ed. (1996) *Escherichia coli and Salmonella Cellular and Molecular Biology* 2nd ed, ASM Press, Washington, D.C.

Technical Appendix

Nucleic Acids and Proteins: Calculations.

An online calculator for these values is available in the "tools" section of Promega's Web site at: www.promega.com/techserv/tools/

Metric Prefixes

Prefix	Symbol	Factor
kilo	k	10 ³
centi	c	10 ⁻²
milli	m	10 ⁻³
micro	μ	10 ⁻⁶
nano	n	10 ⁻⁹
pico	p	10 ⁻¹²
femto	f	10 ⁻¹⁵
atto	a	10 ⁻¹⁸
zepto	z	10 ⁻²¹

Spectrophotometric Conversions

- 1 A₂₆₀ unit of double-stranded DNA = 50 μg/ml
- 1 A₂₆₀ unit of single-stranded DNA = 33 μg/ml
- 1 A₂₆₀ unit of single-stranded RNA = 40 μg/ml

DNA Molar Conversions

- 1 μg of 1,000bp DNA = 1.52 pmol (3.03 pmol of ends)
- 1 μg of pBR322 DNA = 0.36 pmol DNA
- 1 pmol of 1,000bp DNA = 0.66 μg
- 1 pmol of pBR322 DNA = 2.8 μg

Formulas for DNA Molar Conversions

For dsDNA:

To convert pmol to μg:

$$\text{pmol} \times N \times \frac{660 \text{ pg}}{\text{pmol}} \times \frac{1 \mu\text{g}}{10^6 \text{ pg}} = \mu\text{g}$$

To convert μg to pmol:

$$\mu\text{g} \times \frac{10^6 \text{ pg}}{1 \mu\text{g}} \times \frac{\text{pmol}}{660 \text{ pg}} \times \frac{1}{N} = \text{pmol}$$

where N is the number of nucleotide pairs and 660 pg/pmol is the average MW of a nucleotide pair.

For ssDNA:

To convert pmol to μg:

$$\text{pmol} \times N \times \frac{330 \text{ pg}}{\text{pmol}} \times \frac{1 \mu\text{g}}{10^6 \text{ pg}} = \mu\text{g}$$

To convert μg to pmol:

$$\mu\text{g} \times \frac{10^6 \text{ pg}}{1 \mu\text{g}} \times \frac{\text{pmol}}{330 \text{ pg}} \times \frac{1}{N} = \text{pmol}$$

where N is the number of nucleotides and 330 pg/pmol is the average MW of a nucleotide.

Dalton (Da) is an alternate name for the atomic mass unit, and kiloDalton (kDa) is 1,000 Daltons. Thus a protein with a mass of 64 kDa has a molecular weight of 64,000 grams per mole.

- (a) The PCR process is covered by patents issued and applicable in certain countries*. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process. *In the U.S., effective March 29, 2005, U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 will expire. In Europe, effective March 28, 2006, European Pat. Nos. 201,184 and 200,362 will expire.
- (b) Purchase of this product is accompanied by a limited license under U.S. Pat. Nos. 5,082,784 and 5,192,675 for the internal research use of the computer.
- (c) Turbo™ Nae I and Turbo™ Nar I are the subjects of U.S. Pat. Nos. 5,248,600 and 5,418,150 and are licensed exclusively to Promega Corporation, as well as a license under DD 264 231.
- (d) Licensed using U.S. Pat. No. 4,935,361.
- (e) Licensed under U.S. Pat. No. 5,7075.
- (f) Licensed using one or more of U.S. Pat. Nos. 5,487,993 and 5,827,657 and European Pat. No. 0 550 693.
- (g) U.S. Pat. No. 4,766,072.
- (h) The PCR process, which is the subject of European Pat. Nos. 201,184 and 200,362 and U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 owned by Hoffmann-LaRoche*, is covered by patents issued and applicable in certain countries. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process. Use of this product is recommended for persons that either have a license to perform PCR or are not required to obtain a license. *In the U.S., effective March 29, 2005, the above primary U.S. Pat. Nos. 4,683,195, 4,965,188 and 4,683,202 will expire. In Europe, effective March 28, 2006, the above primary European Pat. Nos. 201,184 and 200,362 will expire.
- (i) Certain applications of this product are covered by patents issued and applicable in certain countries. Because purchase of this product does not include a license to perform any patented application, users of this product may be required to obtain a patent license depending upon the particular application and country in which the product is used.
- (j) Australian Pat. No. 730718 and other patents and patents pending.
- (k) U.S. Pat. No. 5,981,235, Australian Pat. No. 729932 and other patents pending.

©2004 Promega Corporation. All Rights Reserved.

All prices and specifications are subject to change without prior notice. Product claims are subject to change. Please contact Promega Technical Services or access Promega online for the most up-to-date information on Promega products.

GeneEditor, LigaFast, MULTICORE, Select96, TagBead, and Turbo are trademarks of Promega Corporation. 4-CORE, Altered Sites, GoTaq, pALTER, pGEM, VacMan and Wizard are trademarks of Promega Corporation and are registered with the U.S. Patent and Trademark Office.

ABLE and SURE are registered trademarks of Stratagene. Bacto is a registered trademark of Difco Laboratories, Detroit, Michigan. DH5α is a trademark of Life Technologies, Inc. Ficoll is a registered trademark of Amersham Biosciences Ltd. Stb12 adn Stb14 are trademarks of Invitrogen Corporation.