

# Sequencing Primers

## Sequencing Primers.

Primer Sequence	Cat.#	T <sub>m</sub> (°C)*	Applicable Vectors
<b>RNA Polymerase Promoter Primers</b>			
<b>SP6</b>			
5'-d(TATTTAGGTGACACTATAG)-3'	Q5011	42	pGEM <sup>®</sup> -4Z, -3Zf(+/-), -5Zf(+/-), -7Zf(+/-), -9Zf(-), -11Zf(+/-), -13Zf(+); pGEM <sup>®</sup> - <i>luc</i> ; pSP- <i>luc</i> +; pSP- <i>luc</i> +NF; pALTER <sup>®</sup> -1; pGEM <sup>®</sup> -T, -T Easy; PinPoint <sup>™</sup> Xa-1, -2, -3; pTnT <sup>™</sup> Vector; pCMVTnT <sup>™</sup> Vector; pGeneClip <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin, hMGFP
<b>T7</b>			
5'-d(TAATACGACTCACTATAGGG)-3'	Q5021	47	pGEM <sup>®</sup> -4Z, -3Zf(+/-), -5Zf(+/-), -7Zf(+/-), -9Zf(-), -11Zf(+/-), -13Zf(+); pGEM <sup>®</sup> - <i>luc</i> ; pSP- <i>luc</i> +NF; pALTER <sup>®</sup> -1; pGEM <sup>®</sup> -T, -T Easy; pTARGET <sup>™</sup> ; pGeneClip <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin, hMGFP; pF1A, pF1K T7 Flexi <sup>®</sup> Vector; pFN2A, pFN2K (GST) Flexi <sup>®</sup> Vector; pF4A, pF4K CMV Flexi <sup>®</sup> Vector; pF5A, pF5K CMV- <i>neo</i> Flexi <sup>®</sup> Vector; pFC8A, pFC8K (HaloTag <sup>®</sup> ) CMV Flexi <sup>®</sup> Vector; pF9 CMV <i>Rluc</i> - <i>neo</i> Flexi <sup>®</sup> Vector; pFN10A (ACT) Flexi <sup>®</sup> Vector; pFN11A (BIND) Flexi <sup>®</sup> Vector
<b>T3</b>			
5'-d(ATAACCCCTCACTAAAGGGA)-3'	Q5741	50	pACT, pALTER <sup>®</sup> -MAX, pBIND; pCI- <i>neo</i> ; HaloTag <sup>®</sup> pHT2
<b>T7 EEV</b>			
5'-d(AAGGCTAGAGTACTTAATACGA)-3'	Q6700	50	pALTER <sup>®</sup> -MAX, pSI, pCI, pCI- <i>neo</i> , pCMVTnT <sup>™</sup> , pTnT <sup>™</sup> , pHMGFP Vector; HaloTag <sup>®</sup> pHT2; psiCHECK <sup>™</sup> -1, -2
<b>pUC/M13 Primers</b>			
<b>Forward (17mer) (-40)</b>			
5'-d(GTTTCCAGTCACGAC)-3'	Q5391	50	pGEM <sup>®</sup> -3Z, -4Z, -3Zf(+/-), -5Zf(+/-), -7Zf(+/-), -9Zf(-), -11Zf(+/-), -13Zf(+); pGEM <sup>®</sup> - <i>luc</i> ; pGEM <sup>®</sup> -T, -T Easy; pALTER <sup>®</sup> -1; pCAT <sup>®</sup> -Basic, -Promoter, -Enhancer, -Control; pGeneClip <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin, hMGFP
<b>Forward (24mer) (-47)</b>			
5'-d(CGCCAGGGTTTCCAGTCACGAC)-3'	Q5601	64	[same as Forward (17mer)]
<b>Reverse (17mer)</b>			
5'-d(CAGGAAACAGCTATGAC)-3'	Q5401	47	pGEM <sup>®</sup> -3Z, -4Z, -3Zf(+/-), -5Zf(+/-), -7Zf(+/-), -9Zf(-), -11Zf(+/-), -13Zf(+); pGEM <sup>®</sup> - <i>luc</i> ; pSP64, -64 Poly(A), -65; pGEM <sup>®</sup> -T, -T Easy; pCAT <sup>®</sup> -Basic, -Promoter, -Enhancer, -Control; pALTER <sup>®</sup> -1, <i>Ex</i> -1, <i>Ex</i> -2; pGeneClip <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin, hMGFP; psiSTRIKE <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin, hMGFP; psiLentGene <sup>™</sup> Basic, Puromycin, Hygromycin, Neomycin
<b>Reverse (22mer)</b>			
5'-d(TCACACAGGAAACAGCTATGAC)-3'	Q5421	55	[same as Reverse (17mer)]
<b>Luciferase Primers</b>			
<b>GLprimer1</b>			
5'-d(TGTATCTTATGGTACTGTAAGT)-3'	E1651	50	pGL2-Basic, -Promoter, -Enhancer, -Control
<b>GLprimer2</b>			
5'-d(CTTTATGTTTTGGCGTCTCCA)-3'	E1661	55	pSP- <i>luc</i> +; pGL3-Basic, -Promoter, -Enhancer, -Control; pGL2-Basic, -Promoter, -Enhancer, -Control
<b>RVprimer3</b>			
5'-d(CTAGCAAATAGGCTGTCCC)-3'	E4481	53	pGL3-Basic, -Promoter, -Enhancer, -Control; pCAT <sup>®</sup> 3-Basic, -Promoter, -Enhancer, -Control; pGL3(R2.1)-Basic; pGL3(R2.2)-Basic; pGL4.10[ <i>luc</i> 2]; pGL4.11[ <i>luc</i> 2P]; pGL4.12[ <i>luc</i> 2CP]; pGL4.13[ <i>luc</i> 2/SV40]; pGL4.14[ <i>luc</i> 2/Hygro]; pGL4.15[ <i>luc</i> 2P/Hygro]; pGL4.16[ <i>luc</i> 2CP/Hygro]; pGL4.17[ <i>luc</i> 2/Neo]; pGL4.18[ <i>luc</i> 2P/Neo]; pGL4.19[ <i>luc</i> 2CP/Neo]; pGL4.20[ <i>luc</i> 2Puro]; pGL4.21[ <i>luc</i> 2P/Puro]; pGL4.22[ <i>luc</i> 2CP/Puro]; pGL4.70[ <i>hRluc</i> ]; pGL4.71[ <i>hRluc</i> P]; pGL4.72[ <i>hRluc</i> CP]; pGL4.73[ <i>hRluc</i> /SV40]; pGL4.74[ <i>hRluc</i> /TK]; pGL4.75[ <i>hRluc</i> /CMV]; pGL4.76[ <i>hRluc</i> /Hygro]; pGL4.77[ <i>hRluc</i> P/Hygro]; pGL4.78[ <i>hRluc</i> CP/Hygro]; pGL4.79[ <i>hRluc</i> /Neo]; pGL4.80[ <i>hRluc</i> P/Neo]; pGL4.81[ <i>hRluc</i> CP/Neo]; pGL4.82[ <i>hRluc</i> /Puro]; pGL4.83[ <i>hRluc</i> P/Puro]; pGL4.84[ <i>hRluc</i> CP/Puro]; pCBR-Basic, -Control; pCBG68-Basic, -Control; pCBG99-Basic, -Control
<b>RVprimer4</b>			
5'-d(GACGATAGTCATGCCCGCG)-3'	E4491	60	pGL3-Basic, -Promoter, -Enhancer, -Control; pCAT <sup>®</sup> 3-Basic, -Promoter, -Enhancer, -Control; pGL3(R2.1)-Basic; pGL3(R2.2)-Basic; pGL4.10[ <i>luc</i> 2]; pGL4.11[ <i>luc</i> 2P]; pGL4.12[ <i>luc</i> 2CP]; pGL4.13[ <i>luc</i> 2/SV40]; pGL4.14[ <i>luc</i> 2/Hygro]; pGL4.15[ <i>luc</i> 2P/Hygro]; pGL4.16[ <i>luc</i> 2CP/Hygro]; pGL4.17[ <i>luc</i> 2/Neo]; pGL4.18[ <i>luc</i> 2P/Neo]; pGL4.19[ <i>luc</i> 2CP/Neo]; pGL4.20[ <i>luc</i> 2Puro]; pGL4.21[ <i>luc</i> 2P/Puro]; pGL4.22[ <i>luc</i> 2CP/Puro]; pGL4.70[ <i>hRluc</i> ]; pGL4.71[ <i>hRluc</i> P]; pGL4.72[ <i>hRluc</i> CP]; pGL4.73[ <i>hRluc</i> /SV40]; pGL4.74[ <i>hRluc</i> /TK]; pGL4.75[ <i>hRluc</i> /CMV]; pGL4.76[ <i>hRluc</i> /Hygro]; pGL4.77[ <i>hRluc</i> P/Hygro]; pGL4.78[ <i>hRluc</i> CP/Hygro]; pGL4.79[ <i>hRluc</i> /Neo]; pGL4.80[ <i>hRluc</i> P/Neo]; pGL4.81[ <i>hRluc</i> CP/Neo]; pGL4.82[ <i>hRluc</i> /Puro]; pGL4.83[ <i>hRluc</i> P/Puro]; pGL4.84[ <i>hRluc</i> CP/Puro]; pCBR-Basic, -Control; pCBG68-Basic, -Control; pCBG99-Basic, -Control
<b>Miscellaneous Sequencing Primers</b>			
<b>PinPoint<sup>™</sup> Sequencing Primer</b>			
5'-d(CGTGACGCGGTGCAGGGCG)-3'	V4211	66	PinPoint <sup>™</sup> Xa-1, -2, -3; PinPoint <sup>™</sup> Xa-1 T-Vector
<b>pTARGET<sup>™</sup> Sequencing Primer</b>			
5'-d(TTAGCCAAAGTTATTTAGGTGACA)-3'	Q4461	55	pTARGET <sup>™</sup> Vector

\*Melting temperatures (T<sub>m</sub>) for each oligonucleotide determined by the base-stacking free-energy method are listed on Promega's technical resources Web site at: [www.promega.com/biomath/](http://www.promega.com/biomath/)  
 N/A = Not available from Promega.