siCHECKTM Vectors

Instructions for Use of Products **C8011 and C8021**



Revised 4/16 TB329



siCHECKTM Vectors

All technical literature is available at: www.promega.com/protocols/ Visit the web site to verify that you are using the most current version of this Technical Bulletin. E-mail Promega Technical Services if you have questions on use of this system: techserv@promega.com

1.	Description	2
	Product Components and Storage Conditions	
	General Considerations	
	3.A. siCHECK™ Vector Features	3
	3.B. How the siCHECK™ Vectors Work	
	3.C. Sample Experiments Using the siCHECK™ Vectors	
4.	siCHECK™ Vector Maps	9
5.	siCHECK™ Vector Restriction Enzyme Tables	11
	5.A. Restriction Enzyme Sites for the psiCHECK™-1 Vector	11
	5.B. Restriction Enzyme Sites for the psiCHECK™-2 Vector	
6.	siCHECK™ Vector Backbones and Components	15
7.	References	16
8.	Related Products	18
9.	Summary of Changes	18



1. Description

The psiCHECKTM-1 Vector^(a,b,c) (Cat.# C8011) and psiCHECKTM-2 Vector^(a,b,c) (Cat.# C8021) are designed to provide a quantitative and rapid approach for optimization of RNA interference (RNAi). The vectors enable the monitoring of changes in expression of a target gene fused to the reporter gene. In both vectors, *Renilla* luciferase is used as a primary reporter gene, and the gene of interest can be cloned into the multiple cloning region located downstream of the *Renilla* luciferase translational stop codon. Initiation of the RNAi process toward a gene of interest results in cleavage and subsequent degradation of fusion mRNA. Measurement of decreased *Renilla* luciferase activity is a convenient indicator of RNAi effect (1).

RNAi is a phenomenon by which double-stranded RNA complementary to a target mRNA can specifically inactivate gene function by stimulating the degradation of the target mRNA (2–4). Because of the ability to inactivate genes, RNAi has emerged as a powerful tool for analyzing gene function.

In mammalian systems, including cultured mammalian cells, chemically synthesized double-stranded short interfering RNA molecules (<30 nucleotides; siRNA) result in dsRNA duplexes <30 base pairs in length that induce RNAi (5–10). RNAi duplexes >30bp induce the interferon response and nonspecific degradation of mRNA and cannot be used as tools for specific gene silencing (11, 12).

Interestingly, a significant percentage of the siRNA designed for a specific gene are not effective (5, 13–16). On average only 1 in 5 of the siRNAs selected for targeting a specific region show efficient gene silencing (16, 17). Possible causes for the failure of a particular siRNA may be instability of an siRNA probe in vivo, inability to interact with components of the RNAi machinery or the inaccessibility of the target mRNA due to local secondary structural constraints. Analysis of nucleotide sequences, melting temperatures and secondary structures have not revealed any obvious difference between effective and ineffective siRNA (18).

At present, one of the most serious limitations for the RNAi technology is the lack of a rapid, reliable, quantitative target-site screening method. Various algorithm programs exist that aid in the design of potential siRNA targets. However, an experimental method is needed to screen these siRNAs. Current screening technologies include such semi-quantitative, time-consuming methods as fluorescence change for GFP-target fusions, Western blot analysis, monitoring phenotypic changes or RT-PCR. In addition, the current screening technologies are not easily modified for the rapid, simultaneous screening of multiple siRNA.

2. Product Components and Storage Conditions

2

PRODUCT	SIZE	CAT.#
psiCHECK™-1 Vector	20μg	C8011
psiCHECK™-2 Vector	20μg	C8021

Storage Conditions: Store the psiCHECKTM-1 and psiCHECKTM-2 Vectors at -20° C.



3. General Considerations

3.A. siCHECKTM Vector Features

Current methods to monitor changes in gene expression as the result of RNAi are either semi-quantitative, time-consuming or not applicable to high-throughput screening. The siCHECKTM Vectors are easier to use than currently available methods, allow optimal quantitative target site selection and can be adapted for use in high-throughput methodologies.

There are two siCHECKTM Vectors, the psiCHECKTM-1 Vector and the psiCHECKTM-2 Vector. Both vectors contain as the primary reporter gene the synthetic version of *Renilla* luciferase, hRluc, which is used to monitor changes in expression as the result of RNAi induction. This synthetic gene is engineered for more efficient expression in mammalian cells and for reduced anomalous transcription.

To aid in fusion of the target gene to the synthetic *Renilla* luciferase reporter gene, a region of restriction sites (i.e., the multiple cloning region) has been added 3′ to the *Renilla* translational stop. The restriction sites present in the multiple cloning region can be used to create genetic fusions between the gene of interest and the *Renilla* reporter gene. Because no fusion protein is expressed, there is no need to be concerned about whether you have cloned into a proper translational reading frame.

The multiple cloning region of the psiCHECK $^{\text{TM}}$ -1 Vector contains unique restriction sites SgfI, XhoI, SmaI, EcoRI, PmeI and NotI. Due to the presence of the firefly expression cassette, the psiCHECK $^{\text{TM}}$ -2 Vector contains fewer unique restriction sites. The restriction sites in the psiCHECK $^{\text{TM}}$ -2 Vector multiple cloning region are SgfI, XhoI, PmeI and NotI.

The promoter used for *Renilla* luciferase expression in the siCHECK™ Vectors is the SV40 promoter. Experimental results (data not shown) demonstrate that the SV40 promoter results in the best balance between *Renilla* luciferase expression and the detection of RNAi activity when used with siRNA.

The difference between the two siCHECKTM Vectors is that the psiCHECKTM-2 Vector possesses a secondary firefly reporter expression cassette. The firefly expression cassette consists of an HSV-TK promoter, a synthetic firefly luciferase gene and an SV40 late poly(A) signal. To reduce the potential for recombination events, the *Renilla* luciferase reporter gene in the psiCHECKTM-2 Vector uses a synthetic poly(A). This firefly reporter cassette has been specifically designed to be an intraplasmid transfection normalization reporter; thus when using the psiCHECKTM-2 Vector, the *Renilla* luciferase signal can be normalized to the firefly luciferase signal.

If no transfection normalization is required or one would prefer to have the transfection normalization reporter on a second plasmid, the psi $CHECK^{TM}$ -1 Vector is the vector of choice.



3.A. siCHECK™ Vector Features (continued)

The psiCHECKTM-1 Vector is recommended for use in monitoring RNAi effects in live cells. The changes in *Renilla* luciferase activity are measured with EnduRenTM Live Cell Substrate (Cat.# E6481), which allows continuous monitoring of intracellular *Renilla* luminescence (19; Figure 2). EnduRenTM Live Cell Substrate is for use only with *Renilla* luciferase.

Promega offers several reagents that can be used in conjunction with the siCHECK™ Vectors to monitor *Renilla* and firefly luciferase signals. For the psiCHECK™-1 Vector, which only contains the *Renilla* luciferase reporter gene, the *Renilla* Luciferase Assay System (Cat.# E2810, E2820) can be used. The psiCHECK™-2 Vector, which contains *Renilla* and firefly luciferase reporter genes, requires the use of either the Dual-Luciferase® Reporter Assay System (Cat.# E1910) or the Dual-Glo™ Luciferase Assay System (Cat.# E2920) to generate the firefly and *Renilla* luciferase signals.

3.B. How the siCHECK™ Vectors Work

Figure 1 provides a basic description of how the siCHECKTM Vectors work. Using the unique restriction sites, the gene of interest is cloned into the multiple cloning region located 3 to the synthetic Renilla luciferase gene and its translational stop codon. After cloning, the vector is transfected into the mammalian cell line of choice, and a fusion of the Renilla luciferase gene and the gene of interest is transcribed. Vectors expressing potential siRNA can be cotransfected simultaneously or sequentially, depending on your experimental design. If a specific siRNA binds to the target mRNA and initiates the RNAi process, the fused Renilla luciferase:gene of interest mRNA will be cleaved and subsequently degraded, decreasing the Renilla luciferase signal.



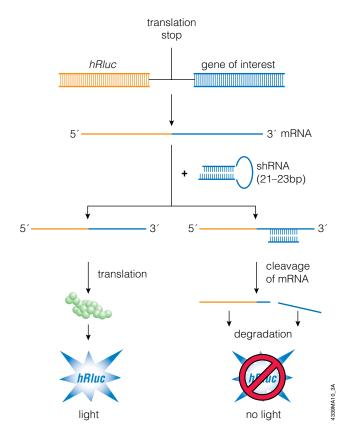


Figure 1. Mechanism of action of the siCHECK™ Vectors.



6

3.C. Sample Experiments Using the siCHECK™ Vectors

To demonstrate the utility of the siCHECK[™] Vectors, two experiments are detailed in this Technical Bulletin. In the first experiment, human p53 cDNA was subcloned into the psiCHECK[™]-1 and the psiCHECK[™]-2 Vectors using the SgfI and NotI restriction sites located in the multiple cloning region of both vectors. Note the SgfI and NotI restriction sites are located 3′ to the *Renilla* luciferase translational stop codon. As shown in Figure 2, the psiCHECK[™]-1 Vector containing the human p53 cDNA was cotransfected into HEK-293T cells with the psiLentGene[™] Basic Vector expressing either a *Renilla* luciferase (*hRluc*). The negative control was the psiLentGene[™] Basic Vector with a nonspecific 19bp insert. (A BLAST search using this 19bp sequence and a threshold >90% revealed no homology to any known mammalian gene or to the synthetic *Renilla* luciferase gene.) This nonspecific sequence was used for all RNAi experiments in this Technical Bulletin.

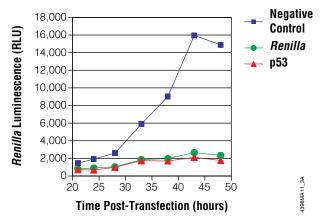


Figure 2. Inhibition of *Renilla* luciferase expression by targeting either the *Renilla* luciferase or p53 gene. The human p53 cDNA was subcloned into the psiCHECK[™]-1 Vector using the SgfI and NotI restriction sites located 3′ to the *Renilla* luciferase translational stop codon. To begin the transfection assay, HEK-293T cells were plated in a 96-well plate at 3,000 cells/well. After an overnight incubation, the cells were treated with a transfection mixture consisting of 35μl of serum-free medium, 0.3μl of TransFast[™] Transfection Reagent (Cat.# E2431), 0.02μg of psiCHECK[™]-1:p53 vector and 0.08μg of psiLentGene[™] Basic Vector per well. For this experiment, the psiLentGene[™] Vector expressed shRNAs directed against human p53, *Renilla* luciferase or the nonspecific 19bp sequence, which serves as a negative control, (Section 3.C). After a one-hour incubation, 100μl of serum-containing medium was added to the wells. At 21 hours post-transfection, EnduRen[™] Live Cell Substrate (Cat.# E6481) was added to a final concentration of 60μM, and *Renilla* luciferase activity was monitored. *Renilla* luciferase activities were normalized to the number of viable cells using the CellTiter-Glo[®] Luminescent Cell Viability Assay (Cat.# G7573; 20).



At 21 hours post-transfection, nonlytic EnduRenTM Live Cell Substrate was added to the wells; luminescence was monitored for the next 27 hours until 48 hours post-transfection. The data in Figure 2 show that the psiLentGeneTM Basic Vector expressing either *Renilla* luciferase inhibits the expression of the *Renilla* luciferase reporter gene from the psiCHECKTM-1:p53 vector. Interestingly, using either *Renilla* luciferase results in virtually identical inhibition of *Renilla* luciferase expression.

In a second experiment, the human p53 cDNA used in Figure 2 was subcloned into the psiCHECK™-2 Vector using the SgfI and NotI restriction sites. Five potential p53 shRNAs designed to bind to five different target sites were cloned into the psiLentGene™ Basic Vector; the resulting vectors were named Site 1 through Site 5. The control is a psiLentGene™ Vector containing the nonspecific 19bp sequence. The psiCHECK™-2 Vector containing the p53 cDNA was cotransfected with the psiLentGene™ Vector expressing either a p53 shRNA (Figure 3, Sites 1−5) or the nonspecific shRNA into HEK-293T cells as described in Figure 3. Forty-eight hours after transfection, the medium was removed and cells were lysed in Passive Lysis Buffer (Cat.# E1941). The firefly and *Renilla* luciferase signals were generated using the Dual-Luciferase® Reporter 1000 Assay System (21).

Figure 3, Panel A, displays the *Renilla* luciferase signal, while Figure 3, Panel B, shows the *Renilla* luciferase signal normalized (corrected for transfection efficiency to the firefly luciferase signal). The data in Figure 3, Panel A, is difficult to interpret due to transfection variations. The *Renilla* luciferase positive control, which should demonstrate inhibition of reporter expression, is not statistically different (i.e., overlapping error bars) from the negative control (no effect on reporter expression was detected). The inability to distinguish between the positive and negative controls renders any conclusion regarding the effectiveness of potential shRNAs suspect.

However, when the *Renilla* luciferase signals are normalized (see Figure 3, Panel B) to the internal firefly luciferase transfection control, the data interpretation is different, as the *Renilla* luciferase positive control is statistically different from the negative control. In addition, the normalized data allow the ability to distinguish the effectiveness of the various target site shRNAs.



3.C. Sample Experiments Using the siCHECK™ Vectors (continued)

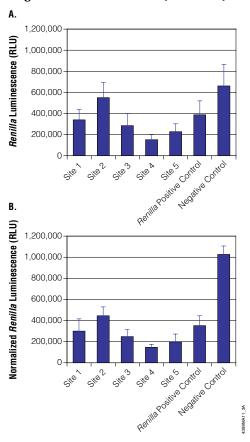


Figure 3. Target site selection using the psiCHECK™-2 Vector. HEK-293T cells were seeded into a 96-well plate at a density of 3,000 cells/well. Human p53 cDNA was subcloned into the psiCHECK™-2 Vector using the SgfI and NotI restriction sites. After an overnight incubation, the cells were treated with a transfection mixture consisting of 35μl of serum-free medium, 0.3μl of TransFast™ Transfection Reagent (Cat.# E2431), 0.02μg of psiCHECK™-2 Vector: p53 and 0.08μg of psiLentGene™ Basic Vector per well. The psiLentGene™ Basic Vector expressed one of five different shRNAs directed against human p53, *Renilla* luciferase or a nonspecific 19bp sequence (Section 3.C) as a negative control. After a one-hour incubation, 100μl of serum-containing medium was added to the wells. Forty-eight hours post-transfection *Renilla* and firefly luciferase activities were measured using the Dual-Luciferase® Reporter 1000 Assay System (Cat.# E1980; 21). Panel A displays the raw *Renilla* luciferase data, while in Panel B, the *Renilla* luciferase data has been normalized to firefly luciferase data. The data represent the mean of 12 wells plus or minus the standard deviation. Note that in other experiments the ability of different shRNAs to inhibit gene expression might vary more dramatically.



4. siCHECK™ Vector Maps

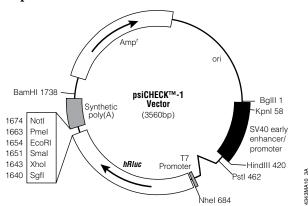


Figure 4. psiCHECK™-1 Vector map. -^- denotes the intron.

psiCHECK™-1 Vector sequence reference points:

SV40 early enhancer/promoter	7-425
Chimeric intron	489-621
T7 RNA polymerase promoter	666-684
Synthetic Renilla luciferase gene (hRluc)	694-1629
Multiple cloning region	1636-1680
Synthetic poly(A)	1688-1736
β-lactamase (Amp ^r) coding region	1874-2734

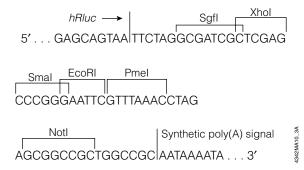


Figure 5. psiCHECK™-1 Vector multiple cloning region.



4. siCHECKTM Vector Maps (continued)

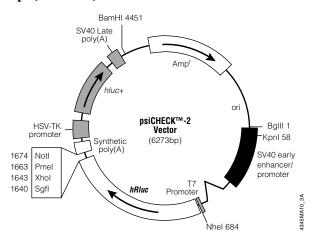


Figure 6. psiCHECK™-2 Vector map. -^- denotes the intron.

psiCHECK™-2 Vector sequence reference points:

SV40 early enhancer/promoter	7-425
Chimeric intron	489-621
T7 RNA polymerase promoter	666-684
Synthetic Renilla luciferase gene (hRluc)	694-1629
Multiple cloning region	1636-1680
Synthetic poly(A)	1688-1736
HSV-TK promoter	1744-2496
Synthetic firefly luciferase gene (hluc+)	2532-4184
SV40 late poly(A)	4219-4440
β-lactamase (Amp ^r) coding region	4587-5447

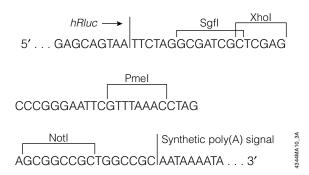


Figure 7. psiCHECK™-2 Vector multiple cloning region.

10



5. siCHECK[™] Vector Restriction Enzyme Tables

5.A. Restriction Enzyme Sites for the psiCHECK™-1 Vector

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3′-end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or to report a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. Vector sequences are available from the GenBank® database (GenBank®/EMBL accession number AY535006) and online at: www.promega.com/vectors/

Table 1. Restriction Enzymes That Cut the psiCHECK™-1 Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	1391	Cfr10I	1	2576
Acc65I	1	54	DraI	4	1663, 2083, 2775, 2794
AcyI	2	1388, 2121	DraII	1	1539
AflII	2	452, 649	DraIII	1	882
Alw44I	2	1989, 3235	DrdI	2	441, 3447
AlwNI	1	3140	DsaI	4	15, 311, 692, 899
AspHI	4	1091, 1993, 2078, 3239	EaeI	3	1674, 1681, 2268
AvaI	3	715, 1643, 1649	EagI	1	1674
AvaII	2	2297, 2519	EarI	2	1193, 1862
AvrII	1	404	EclHKI	1	2661
BamHI	1	1738	Eco52I	1	1674
BanI	3	54, 575, 2708	Eco81I	1	1280
BanII	3	759, 899, 1650	EcoRI	1	1654
BbsI	1	560	EcoRV	1	1179
BbuI	2	152, 224	FspI	2	8, 2438
BclI	2	734, 1187	HaeII	1	3309
BglI	3	357, 694, 2543	HgaI	4	1570, 2129, 2859, 3437
BglII	1	1	HindIII	1	420
BsaI	3	514, 1234, 2595	Hsp92I	2	1388, 2121
BsaOI	5	1640, 1677, 2143, 2292, 3215	KpnI	1	58
BsaBI	1	1453	MspA1I	5	80, 1679, 2025, 2966, 3211
BsaHI	2	1388, 2121	NciI	5	1650, 1651, 2125, 2476, 3172
BspHI	2	1821, 2829	NcoI	3	15, 311, 692
BspMI	1	476	NheI	1	684
BssSI	2	1992, 3376	NotI	1	1674
Bst98I	2	452, 649	NruI	1	1355
BstZI	1	1280	NsiI	3	154, 226, 913



5.A. Restriction Enzyme Sites for the psiCHECK™-1 Vector (continued)

Table 1. Restriction Enzymes That Cut the psiCHECK™-1 Vector Between 1 and 5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
NspI	2	152, 224	SmaI	1	1651
PaeR7I	1	1643	SphI	2	152, 224
PmeI	1	1663	SspI	1	1856
Ppu10I	3	150, 222, 909	StuI	1	403
PspAI	1	1649	StyI	5	15, 311, 404, 692, 701
PstI	1	462	TfiI	2	426, 805
PvuI	2	1640, 2292	Tth111I	1	1390
PvuII	1	80	VspI	1	2486
ScaI	2	662, 2180	XhoI	1	1643
SfiI	1	357	XmaI	1	1649
SgfI	1	1640	XmnI	2	1228, 2061
SinI	2	2297, 2519			

Table 2. Restriction Enzymes That Do Not Cut the psiCHECK™-1 Vector.

AccB7I	BsaAI	Eco47III	NaeI	SacI
AccI	BsaMI	Eco72I	NarI	SacII
AccIII	BsmI	EcoICRI	NdeI	SalI
AflIII	Bsp120I	EcoNI	NgoMIV	SgrAI
AgeI	BsrGI	EheI	PacI	SnaBI
ApaI	BssHII	FseI	PflMI	SpeI
AscI	Bst1107I	HincII	PinAI	SplI
BalI	BstEII	HindII	PmlI	SrfI
BbeI	BstXI	HpaI	PpuMI	Sse8387I
BbrPI	ClaI	I-PpoI	PshAI	SwaI
BlpI	CspI	KasI	Psp5II	XbaI
Bpu1102I	Csp45I	MluI	RsrII	XcmI

Table 3. Restriction Enzymes That Cut the psiCHECK™-1 Vector 6 or More Times.

AciI	BstOI	HhaI	MboII	Sau3AI
AluI	BstUI	HinfI	MnlI	Sau96I
Alw26I	CfoI	HpaII	MseI	ScrFI
BbvI	DdeI	HphI	MspI	SfaNI
BsaJI	DpnI	Hsp92II	NdeII	TaqI
Bsp1286I	DpnII	MaeI	NlaIII	Tru9I
BsrI	Fnu4HI	MaeII	NlaIV	XhoII
BsrSI	FokI	MaeIII	PleI	
Bst71I	HaeIII	MboI	RsaI	

Note: The enzymes listed in boldface type are available from Promega.



5.B. Restriction Enzyme Sites for the psiCHECK™-2 Vector

The following restriction enzyme tables were constructed using DNASTAR® sequence analysis software. Please note that we have not verified this information by restriction digestion with each enzyme listed. The location given specifies the 3´-end of the cut DNA (the base to the left of the cut site). For more information on the cut sites of these enzymes, or to report a discrepancy, please contact your local Promega Branch or Distributor. In the U.S., contact Promega Technical Services at 800-356-9526. Vector sequences are available from the GenBank® database (GenBank®/EMBL accession number AY535007) and online at: www.promega.com/vectors/

Table 4. Restriction Enzymes That Cut the psiCHECK™-2 Vector Between 1 and 5 Times.

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
AatII	1	1391	BssHII	1	1978
AccI	2	2079, 3132	BssSI	3	3459, 4705, 6089
Acc65I	1	54	Bst1107I	1	2080
AflII	4	452, 649, 1773, 1897	Bst98I	4	452, 649, 1773, 1897
AflIII	1	2450	BstXI	1	3650
Alw44I	2	4702, 5948	BstZI	3	1674, 4202, 4206
AlwNI	2	2094, 5853	Bsu36I	3	1280, 3145, 3745
ApaI	1	2562	ClaI	1	4444
AvrII	2	404, 2059	Csp45I	1	2390
BalI	3	1865, 3513, 4038	DraI	5	1663, 4410, 4796, 5488, 5507
BamHI	1	4451	DraIII	1	882
BanII	5	759, 899, 1650, 2050, 2562	DrdI	2	441,6160
BbeI	4	2030, 2815, 3481, 3613	EagI	3	1674, 4202, 4206
BbsI	2	560, 1743	EarI	5	1193, 1874, 2616, 2727, 4575
BbuI	2	152, 224	EclHKI	1	5374
BclI	5	734, 1187, 3112, 3853, 4147	Eco47III	1	3519
BglII	1	1	Eco52I	3	1674, 4202, 4206
BsaI	4	514, 1234, 2123, 5308	Eco81I	3	1280, 3145, 3745
BsaAI	2	2083, 3734	EcoNI	3	2721, 3144, 4149
BsaBI	4	1453, 2979, 4146, 4450	EcoRI	2	1654, 2386
BsaMI	3	2504, 4270, 4363	EcoRV	1	1179
BsmI	3	2504, 4270, 4363	EheI	4	2028, 2813, 3479, 3611
Bsp120I	1	2558	FseI	2	3943, 4208
BspHI	3	3115, 4534, 5542	FspI	3	8, 3354, 5151
BspMI	2	476, 3463	HincII	1	4349
BsrGI	1	3022	HindII	1	4349



5.B. Restriction Enzyme Sites for the psiCHECK™-2 Vector (continued)

Table 4. Restriction Enzymes That Cut the psiCHECK™-2 Vector Between 1 and 5 Times (continued).

Enzyme	# of Sites	Location	Enzyme	# of Sites	Location
HindIII	2	420, 2497	PspAI	2	1649, 2019
HpaI	1	4349	PvuI	2	1640, 5005
KasI	4	2026, 2811, 3477, 3609	PvuII	3	80, 2268, 2606
KpnI	1	58	SacII	1	2036
MluI	1	2450	ScaI	3	662, 2697, 4893
NaeI	3	3941, 3962, 4206	SfiI	1	357
NarI	4	2027, 2812, 3478, 3610	SgfI	1	1640
NcoI	5	15, 311, 692, 2067, 2530	SmaI	2	1651, 2021
NgoMIV	3	3939, 3960, 4204	SphI	2	152, 224
NheI	1	684	SspI	1	4569
NotI	1	1674	StuI	1	403
NruI	1	1355	TfiI	2	426, 805
NsiI	3	154, 226, 913	Tth111I	1	1390
NspI	5	152, 224, 2336, 3023, 3278	VspI	1	5199
PaeR7I	1	1643	XbaI	1	4189
PmeI	1	1663	XhoI	1	1643
Ppu10I	3	150, 222, 909	XmaI	2	1649, 2019
PpuMI	1	2056	XmnI	2	1228, 4774
Psp5II	1	2056			

Table 5. Restriction Enzymes That Do Not Cut the psiCHECK™-2 Vector.

AccB7I	Bpu1102I	NdeI	RsrII	SplI
AccIII	BstEII	PacI	SacI	SrfI
AgeI	CspI	PflMI	SalI	Sse8387I
AscI	Eco72I	PinAI	SgrAI	SwaI
BbrPI	EcoICRI	PmlI	SnaBI	XcmI
BlpI	I-PpoI	PshAI	SpeI	

Note: The enzymes listed in boldface type are available from Promega.



Table 6. Restriction Enzymes That Cut the psiCHECK™-2 Vector 6 or More Times.

AciI	Bsp1286I	EaeI	MaeII	PstI	
AcyI	BsrI	Fnu4HI	MaeIII	RsaI	
AluI	BsrSI	FokI	MboI	Sau3AI	
Alw26I	Bst71I	HaeII	MboII	Sau96I	
AspHI	BstOI	HaeIII	MnlI	ScrFI	
AvaI	BstUI	HgaI	MseI	SfaNI	
AvaII	CfoI	HhaI	MspI	SinI	
BanI	Cfr10I	HinfI	MspA1I	StyI	
BbvI	DdeI	HpaII	NciI	TaqI	
BglI	DpnI	HphI	NdeII	Tru9I	
BsaOI	DpnII	Hsp92I	NlaIII	XhoII	
BsaHI	DraII	Hsp92II	NlaIV		
BsaJI	DsaI	MaeI	PleI		

Note: The enzymes listed in boldface type are available from Promega.

6. siCHECK™ Vector Backbones and Components

The vector backbones of the psiCHECK™-1 and psiCHECK™-2 Vectors are based on the phRL-SV40 Vector (Cat.# E6261). Both the psiCHECK™-1 Vector and psiCHECK™-2 Vector contain the synthetic *Renilla* luciferase reporter gene. The psiCHECK™-2 Vector also contains a synthetic firefly luciferase gene. These synthetic luciferase genes have been codon optimized for more efficient mammalian expression and have been designed with a greatly reduced number of consensus transcription factor binding sites for reduced risk of anomalous transcriptional behavior.

SV40 Early Enhancer/Promoter

The psiCHECK™-1 Vector and psiCHECK™-2 Vector contain the SV40 early enhancer/promoter region, which provides strong, constitutive expression of *Renilla* luciferase in a variety of cell types.

Chimeric Intron

Downstream of the SV40 enhancer/promoter region is a chimeric intron composed of the 5´-donor site from the first intron of the human β -globin and the branch and 3´-acceptor site from the intron that is between the leader and the body of an immunoglobin gene heavy chain variable region (22). The sequences of the donor and acceptor sites, along with the branch point site, have been changed to match the consensus sequence for splicing (23). Transfection studies have demonstrated that the presence of an intron flanking the cDNA insert frequently increases the level of gene expression (24–27).



6. siCHECK[™] Vector Backbones and Components (continued)

T7 Promoter

A T7 RNA polymerase promoter is located downstream of the chimeric intron and immediately precedes the synthetic *Renilla* luciferase reporter gene. This promoter can be used to synthesize RNA transcripts in vitro using T7 RNA Polymerase (Cat.# P2075). Note that the T7 promoter has been verified by sequence only; there has been no functional testing of the T7 promoter.

Polyadenylation Signals (SV40 Late and Synthetic)

Polyadenylation signals are coupled to the termination of transcription by RNA polymerase II and signal the addition of approximately 200–250 adenosine residues to the 3´-end of the RNA transcript (28). Polyadenylation has been shown to enhance RNA stability and translation (29, 30). The late SV40 polyadenylation signal is extremely efficient and has been shown to increase the steady-state level of RNA to approximately fivefold more than that of the early SV40 polyadenylation signal (31). The synthetic poly(A) was cloned from our pCI-neo Vector (Cat.# E1841). The synthetic poly(A) signal is based on the highly efficient polyadenylation signal of the rabbit β -globin gene (32).

7. References

- 1. Kumar, R., Conklin, D.S. and Mittal, V. (2003) High-throughput selection of effective RNAi probes for gene silencing. *Genome Res.* **13**, 2333–40.
- 2. Bass, B.L. (2000) Double-stranded RNA as a template for gene silencing. *Cell* **101**, 235–8.
- 3. Zamore, P.D. (2001) RNA interference: Listening to the sound of silence. Nature Struct. Biol. 8, 746-50.
- 4. Sharp, P.A. (2001) RNA interference—2001. *Genes Dev.* **15**, 485–90.
- 5. Gil, J. and Esteban, M. (2000) Induction of apoptosis by the dsRNA-dependent protein kinase (PKR): Mechanism of action. *Apoptosis* **5**, 107–14.
- 6. Marcus, P.I. and Sekellick, M.J. (1985) Interferon induction by viruses. XIII. Detection and assay of interferon induction-suppressing particles. *Virology* **142**, 411–5.
- 7. Elbashir, S.M. *et al.* (2001) Duplexes of 21-nucleotide RNAs mediate RNA interference in cultured mammalian cells. *Nature* **411**, 494–8.
- 8. Brummelkamp, T.R., Bernards, R. and Agami, R. (2002) A system for stable expression of short interfering RNAs in mammalian cell. *Science* **296**, 550–3.
- 9. Elbashir, S.M. *et al.* (2002) Analysis of gene function in somatic mammalian cells using small interfering RNAs. *Methods* **26**, 199–213.
- 10. Paddison, P.J. *et al.* (2002) Short hairpin RNAs (shRNAs) induce sequence-specific silencing in mammalian cells. *Genes Dev.* **16**, 948–58.
- 11. Paul, C.P. et al. (2002) Effective expression of small interfering RNA in human cells. Nature Biotechnol. 20, 505–8.
- 12. Sui, G. *et al.* (2002) A DNA vector-based RNAi technology to suppress gene expression in mammalian cells. *Proc. Natl. Acad. Sci. USA* **99**, 5515–20.



- 13. Holen, T. et al. (2002) Positional effects of short interfering RNAs targeting the human coagulation trigger tissue factor. Nucl. Acids Res. **30**, 1757–66.
- 14. Lee, N.S. *et al.* (2002) Expression of small interfering RNAs targeted against HIV-1 rev transcripts in human cells. *Nature Biotechnol.* **20**, 500–5.
- 15. Yu, J-Y., DeRuiter, S.L. and Turner, D.L. (2002) RNA interference by expression of short-interfering RNAs and hairpin RNAs in mammalian cells. *Proc. Natl. Acad. Sci. USA* **99**, 6047–52.
- 16. Kapadia, S.B., Brideau-Andersen, A. and Chisari, F.V. (2003) Interference of hepatitis C virus RNA replication by short interfering RNAs. *Proc. Natl. Acad. Sci. USA* **100**, 2014–8.
- 17. McManus, M.T. et al. (2002) Gene silencing using micro-RNA designed hairpins. RNA 8, 842-50.
- 18. Hohjoh, H. (2002) RNA interference (RNA(i)) induction with various types of synthetic oligonucleotide duplexes in cultured human cells. *FEBS Lett.* **521**, 195–9.
- 19. EnduRen™ Live Cell Substrate Technical Manual, #TM244, Promega Corporation.
- 20. CellTiter-Glo® Luminescent Cell Viability Assay Technical Bulletin, #TB288, Promega Corporation.
- 21. Dual-Luciferase® Reporter 1000 Assay System Technical Manual, #TM046, Promega Corporation.
- 22. Bothwell, A.L. *et al.* (1981) Heavy chain variable region contribution to the NPb family of antibodies: Somatic mutation evident in a gamma 2a variable region. *Cell* **24**, 625–37.
- 23. Senapathy, P., Shapiro, M.B. and Harris, N.L. (1990) Splice junctions, branch point sites, and exons: Sequence statistics, identification, and applications to genome project. *Methods Enzymol.* **183**, 252–78.
- 24. Gross, M.K., Kainz, M.S. and Merrill, G.F. (1987) Introns are inconsequential to efficient formation of cellular thymidine kinase mRNA in mouse L cells. *Mol. Cell. Biol.* 7, 4576–81.
- 25. Buchman, A.R. and Berg, P. (1988) Comparison of intron-dependent and intron-independent gene expression. *Mol. Cell. Biol.* **8**, 4395–405.
- 26. Evans, M.J. and Scarpulla, R.C. (1989) Introns in the 3′-untranslated region can inhibit chimeric CAT and beta-galactosidase gene expression. *Gene* **84**, 135–42.
- 27. Huang, M.T. and Gorman, C.M. (1990) Intervening sequences increase efficiency of RNA 3′ processing and accumulation of cytoplasmic RNA. *Nucl Acids Res.* **18**, 937–47.
- 28. Proudfoot, N. (1991) Poly(A) signals. Cell 64, 671-4.
- 29. Bernstein, P. and Ross, J. (1989) Poly(A), poly(A) binding protein and the regulation of mRNA stability. *Trends Biochem. Sci.* **14**, 373–7.
- 30. Jackson, R.J. and Standart, N. (1990) Do the poly(A) tail and 3´ untranslated region control mRNA translation? *Cell* **62**, 15–24.
- 31. Carswell, S. and Alwine, J.C. (1989) Efficiency of utilization of the simian virus 40 late polyadenylation site: Effects of upstream sequences. *Mol. Cell. Biol.* **9**, 4248–58.
- 32. Levitt, N. et al. (1989) Definition of an efficient poly(A) site. Genes Dev. 3, 1019–25.



8. Related Products

Firefly and Renilla Luciferase Reagents

Product	Size	Cat.#
EnduRen™ Live Cell Substrate	0.34mg	E6481
	3.4mg	E6482
	34mg	E6485
Passive Lysis 5X Buffer	30ml	E1941
Glo Lysis Buffer, 1X	100ml	E2661

9. Summary of Changes

The following changes were made to the 4/16 revision of this document:

1. Removed discussion of using this product with shRNA.



(a) BY USE OF THIS PRODUCT, RESEARCHER AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE LABEL LICENSE. If the researcher is not willing to accept the terms of this label license, and the product is unused, Promega will accept return of the unused product and provide the researcher with a full refund.

Researchers may use this product for research use only, no commercial use is allowed. "Commercial use" means any and all uses of this product and derivatives by a party for money or other consideration and may include but is not limited to use in: (1) product manufacture; and (2) to provide a service, information or data; and/or resale of the product or its derivatives, whether or not such product or derivatives are resold for use in research. Researchers shall have no right to modify or otherwise create variations of the nucleotide sequence of the luciferase gene except that researchers may: (1) create fused gene sequences provided that the coding sequence of the resulting luciferase gene has no more than four deoxynucleotides missing at the affected terminus compared to the intact luciferase gene sequence, and (2) insert and remove nucleic acid sequences in splicing research predicated on the inactivation or reconstitution of the luminescence of the encoded luciferase. No other use or transfer of this product or derivatives is authorized without the prior express written consent of Promega. In addition, researchers must either: (1) use luminescent assay reagents purchased from Promega for all determinations of luminescence activity of this product and its derivatives; or (2) contact Promega to obtain a license for use of the product and its derivatives. Researchers may transfer derivatives to others for research use provided that at the time of transfer a copy of this label license is given to the recipients and recipients agree to be bound by the terms of this label license. With respect to any uses outside this label license, including any diagnostic, therapeutic or prophylactic uses, please contact Promega for supply and licensing information. PROMEGA MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING FOR MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THE PRODUCT. The terms of this label license shall be governed under the laws of the State of Wi

 $^{\mbox{\tiny (b)}}\mbox{U.S.}$ Pat. No. 8,008,006 and European Pat. No. 1341808

(c)Certain applications of this product may require licenses from others.

© 2002, 2004, 2005, 2007, 2009, 2015, 2016 Promega Corporation. All Rights Reserved.

CellTiter-Glo, Dual-Glo and Dual-Luciferase are registered trademarks of Promega Corporation. EnduRen, psiCHECK, psiLentGene, siCHECK and TransFast are trademarks of Promega Corporation.

DNASTAR is a registered trademark of DNASTAR, Inc. GenBank is a registered trademark of the U.S. Department of Health and Human Services.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.