

Datasheet of GLuc-ON™ TCF/LEF Transcriptional Response Element (TRE) Lentiviral Clone (TR202)

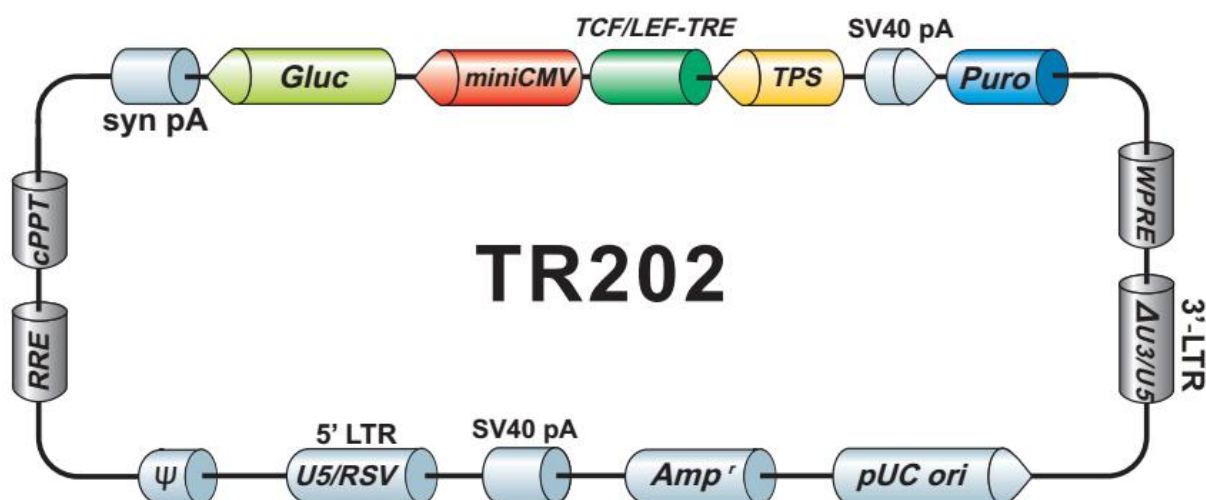


Figure 1. GLuc-ON™ TCF/LEF Transcriptional Response Element (TRE) Lentiviral Clone (TR202)

Cat. No TR202

Description

The TCF/LEF transcriptional response element (TRE) lentiviral clone assays the status of the Wnt signaling pathway through the activity of the transcription factor TCF/LEF. The TCF/LEF transcriptional TRE clone contains a minimal promoter and tandem repeats of the TCF/LEF transcriptional response element upstream of a secreted Gaussia luciferase reporter gene. The number of response elements, as well as the intervening sequence between the elements has been optimized to maximize the sensitivity and signal to noise ratio.

Delivery 500ng/μL × 40μL

Storage Conditions -20°C

QC

- Functional validation in common cell line
- Full sequence verification
- Restriction Digestion

Validation Data

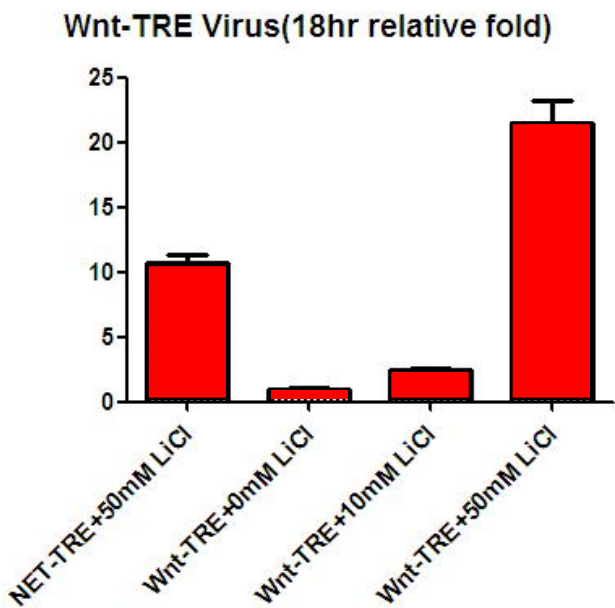


Figure 2. Activation of the GLuc-ON™ Wnt TRE by lithium chloride. HEK293a cells were transduced with negative control lentiviral particles or the GLuc-ON™ TCF/LEF TRE lentiviral particles. Transduced cells were incubated for 48 hours, followed by 6 days incubation in the presence of puromycin. Puromycin-resistant cells were then treated with lithium chloride for 18 hours. A Gaussia Luciferase assay was performed, and transcriptional response activity values are expressed as luminescence fold activation.

Restriction Enzyme Information for TR202

Table 1. Restriction Enzymes That Do Not Cleave TR202

AfeI	AgeI	AscI	BclI	BlpI	BspEI	BsrGI	BstXI
BstZ17I	HpaI	MfeI	NdeI	NotI	PacI	PmeI	PshAI
SbfI	SgfI	SgrAI	SnaBI	SpeI	SrfI	Swal	XbaI

Table 2. Restriction Enzymes That Cleave TR202 Once

AasI	3216	Acc16I	4220	AcvI	715	AjiI	942	AleI	6556	Alol	7848	AspI	1047	AsuII	8290
AviII	4220	AxyI	2339	BaeI	5338	BamHI	7879	BbrPI	715	BbvCI	6403	BmgBI	942	Bpu14I	8290
Bsa29I	8201	BsaAI	715	BsaBI	7840	Bse21I	2339	Bse8I	7840	BseCI	8201	BseJI	7840	BshVI	8201
BsiWI	1061	Bsp119I	8290	BspDI	8201	BspT104I	8290	BstBAI	715	BstBI	8290	BstEII	1139	BstENI	6145
BstPI	1139	Bsu15I	8201	Bsu36I	2339	BtrI	942	Clal	8201	CpoI	1121	CspI	1121	DrdI	3216
DseDI	3216	Eco32I	8522	Eco72I	715	Eco81I	2339	Eco91I	1139	EcoNI	6145	EcoO65I	1139	EcoRI	8263
EcoRV	8522	FseI	6125	FspI	4220	KfiI	6913	MluI	1605	NsbI	4220	NspV	8290	OliI	6556
PaeR7I	7308	Pfi23II	1061	PfiFI	1047	PmaCI	715	PmlI	715	Ppu21I	715	PpuMI	6913	Psp5II	6913
PspCI	715	PspEI	1139	PspLI	1061	PspPPI	6913	PspXI	7308	PstI	2223	PsyI	1047	RigI	6125
Rsr2I	1121	RsrII	1121	Sfr274I	7308	SfuI	8290	SgrDI	4919	Slal	7308	Sspl	4802	StrI	7308
Tth111I	1047	XagI	6145	XcmI	5144	XhoI	7308								

Table 3. Restriction Enzymes That Cleave TR202 Twice

AanI	2684	AanI	4957	AarI	738	AarI	8127	Adel	762	Adel	1576	AhdI	3998	AhdI	7569
AjuI	7228	AjuI	8293	ApaI	516	ApaI	8027	AspA2I	324	AspA2I	550	AspEI	3998	AspEI	7569
AssI	4478	AssI	8196	AsuNHI	2720	AsuNHI	8070	AvrII	324	AvrII	550	Ball	1380	Ball	7777
BauI	3281	BauI	4662	Bcgl	1446	Bcgl	4503	BlnI	324	BlnI	550	BmcAI	4478	BmcAI	8196
BmeRI	3998	BmeRI	7569	BmtI	2724	BmtI	8074	Bpu10I	6403	Bpu10I	8079	BsePI	1460	BsePI	5686
BsmBI	986	BsmBI	1397	Bsp120I	512	Bsp120I	8023	Bsp68I	5808	Bsp68I	7765	BspHI	3825	BspHI	4833
BspOI	2724	BspOI	8074	BspQI	2992	BspQI	6077	BssHII	1460	BssHII	5686	BssSI	3281	BssSI	4662
Bst2BI	3281	Bst2BI	4662	BtuMI	5808	BtuMI	7765	CciI	3825	CciI	4833	Cfr9I	345	Cfr9I	1054
CsiI	91	CsiI	1572	DrallI	762	DrallI	1576	Dril	3998	Dril	7569	Eam1105I	3998	Eam1105I	7569
Eco147I	323	Eco147I	1336	EcoT22I	174	EcoT22I	146	Esp3I	986	Esp3I	1397	KroI	2130	KroI	6121
LguI	2992	LguI	6077	MabI	91	MabI	1572	MlsI	1380	MlsI	7777	MluNI	1380	MluNI	7777
Mph1103I	74	Mph1103I	146	MroNI	2130	MroNI	6121	MscI	1380	MscI	7777	Msp20I	1380	Msp20I	7777
NaeI	2132	NaeI	6123	NgoMIV	2130	NgoMIV	6121	NheI	2720	NheI	8070	NruI	5808	NruI	7765
NsiI	74	NsiI	146	PagI	3825	PagI	4833	Paul	1460	Paul	5686	PceI	323	PceI	1336
PciI	889	PciI	3108	PciSI	2992	PciSI	6077	PdiI	2132	PdiI	6123	PfoI	1480	PfoI	1891
PscI	889	PscI	3108	Psil	2684	Psil	4957	PspOMI	512	PspOMI	8023	Ptel	1460	Ptel	5686
PvuII	2376	PvuII	2932	RcaI	3825	RcaI	4833	RruI	5808	RruI	7765	SapI	2992	SapI	6077
Scal	4478	Scal	8196	SexAI	91	SexAI	1572	SfiI	7968	SfiI	8096	SmaI	347	SmaI	1056
SseBI	323	SseBI	1336	StuI	323	StuI	1336	TspMI	345	TspMI	1054	XmaI	345	XmaI	1054
XmaJI	324	XmaJI	550	Zrml	4478	Zrml	8196	Zsp2I	74	Zsp2I	146				

Table 4. Restriction Enzymes That Cleave TR202 Three times

[illegible]