

Datasheet of GLuc-ON™ EGR1 Transcriptional Response Element (TRE) Lentiviral Clone (TR207)

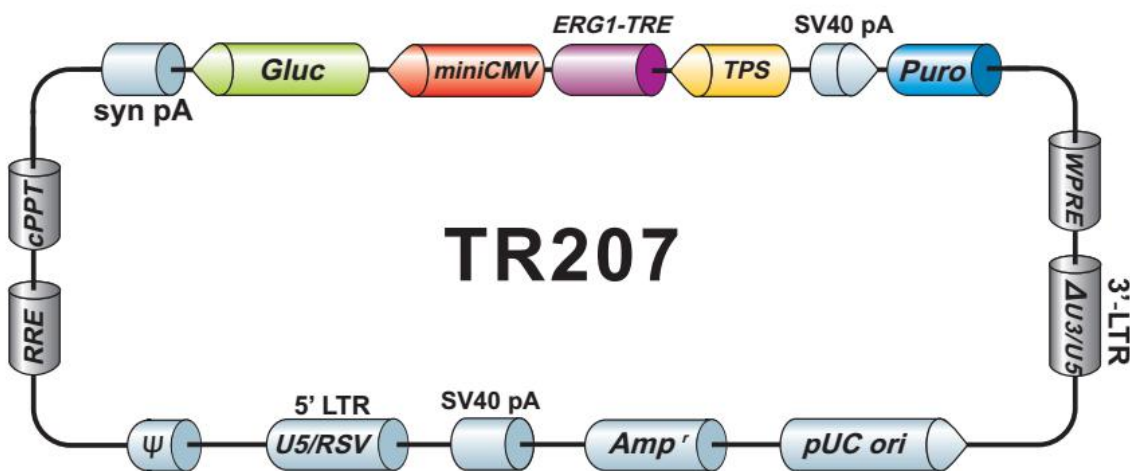


Figure 1. GLuc-ON™ EGR1 Transcriptional Response Element (TRE) Lentiviral Clone (TR207)

Cat. No	TR207
Description	The EGR1 transcriptional response element (TRE) clone assays the status of the EGR1 signaling pathway through the activity of the transcription factor EGR1. The EGR1 transcriptional TRE clone contains a minimal promoter and tandem repeats of the EGR1 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

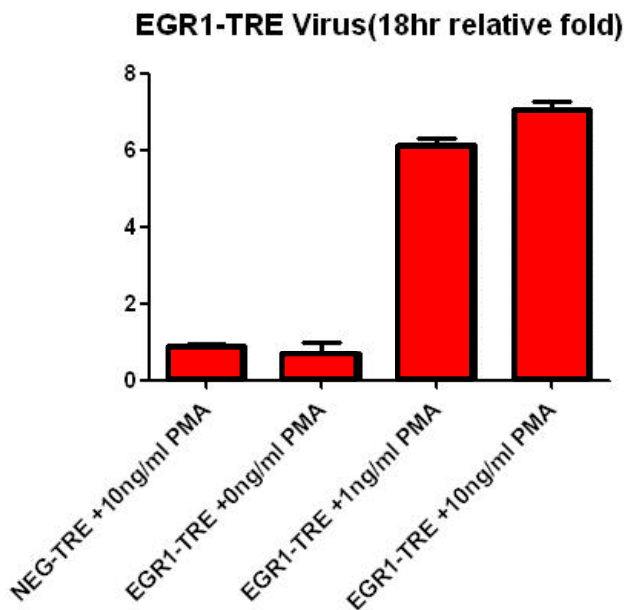


Figure 2. Activation of the GLuc-ON™ EGR1 TRE by phorbol 12-myristate 13-acetate (PMA). HEK293a cells were transduced with negative control lentiviral particles or the GLuc-ON™ EGR1 TRE lentiviral particles. Transduced cells were incubated for 48 hours, followed by 6 days incubation in the presence of puromycin. Puromycin-resistant cells were then treated with PMA for 18 hours. A Gaussia Luciferase assay was performed, and transcriptional response activity values are expressed as luminescence fold activation.

Restriction Enzyme Information for TR207

Table 1. Restriction Enzymes That Do Not Cleave TR207

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 TR207 Once

AasI	3216	Acc16I	4220	AcvI	715	AjiI	942	AleI	6556	Alol	7848	Apal	516	AspI	1047
AsuII	8260	AviII	4220	AxyI	2339	BaeI	5338	BamHI	7879	BbrPI	715	BbvCI	6403	BmgBI	942
Bpu14I	8260	Bsa29I	8171	BsaAI	715	BsaBI	7840	Bse21I	2339	Bse8I	7840	BseCI	8171	BseJI	7840
BshVI	8171	BsiWI	1061	Bsp119I	8260	Bsp120I	512	BspDI	8171	BspT104I	8260	BstBAI	715	BstBI	8260
BstEII	1139	BstENI	6145	BstPI	1139	Bsu15I	8171	Bsu36I	2339	BtrI	942	ClaI	8171	CpoI	1121
CspI	1121	DrdI	3216	DseDI	3216	Eco32I	8492	Eco72I	715	Eco81I	2339	Eco91I	1139	EcoNI	6145
EcoO65I	1139	EcoRI	8233	EcoRV	8492	FseI	6125	FspI	4220	KfiI	6913	MluI	1605	NsbI	4220
NspV	8260	OliI	6556	PaeR7I	7308	Pfi23II	1061	PfiFI	1047	PmaCI	715	PmlI	715	Ppu21I	715
PpuMI	6913	Psp5II	6913	PspCI	715	PspEI	1139	PspLI	1061	PspOMI	512	PspPPI	6913	PspXI	7308
PstI	2223	PsyI	1047	RigI	6125	Rsr2I	1121	RsrII	1121	Sfr274I	7308	SfuI	8260	SgrDI	4919
Slal	7308	Sspl	4802	StrI	7308	Tth111I	1047	XagI	6145	XcmI	5144	XhoI	7308		

Table 3. Restriction Enzymes That Cleave TR207 Twice

AanI	2684	AanI	4957	AarI	738	AarI	8097	Adel	762	Adel	1576	AfIII	2532	AfIII	5492
AhdI	3998	AhdI	7569	AjuI	7228	AjuI	8263	AspA2I	324	AspA2I	550	AspEI	3998	AspEI	7569
AssI	4478	AssI	8166	AsuNHI	2720	AsuNHI	8040	AvrII	324	AvrII	550	Ball	1380	Ball	7777
BauI	3281	BauI	4662	Bcgl	1446	Bcgl	4503	BfrI	2532	BfrI	5492	BinI	324	BinI	550
BmcAI	4478	BmcAI	8166	BmeRI	3998	BmeRI	7569	BmtI	2724	BmtI	8044	Bpu10I	6403	Bpu10I	8049
BsmBI	986	BsmBI	1397	Bsp68I	5808	Bsp68I	7765	BspHI	3825	BspHI	4833	BspOI	2724	BspOI	8044
BspQI	2992	BspQI	6077	BspTI	2532	BspTI	5492	BssSI	3281	BssSI	4662	Bst2BI	3281	Bst2BI	4662
BstAFI	2532	BstAFI	5492	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	MisI	1380	MisI	7777	MluNI	1380	MluNI	7777
Mph1103I	74	Mph1103I	146	MroNI	2130	MroNI	6121	MscI	1380	MscI	7777	Msp20I	1380	Msp20I	7777
MspCI	2532	MspCI	5492	NaeI	2132	NaeI	6123	NgoMIV	2130	NgoMIV	6121	NheI	2720	NheI	8040
NruI	5808	NruI	7765	NsiI	74	NsiI	146	PagI	3825	PagI	4833	PceI	323	PceI	1336
PciI	889	PciI	3108	PciSI	2992	PciSI	6077	PdII	2132	PdII	6123	PfoI	1480	PfoI	1891
PscI	889	PscI	3108	PsiI	2684	PsiI	4957	PvuII	2376	PvuII	2932	RcaI	3825	RcaI	4833
RruI	5808	RruI	7765	SapI	2992	SapI	6077	Scal	4478	Scal	8166	SexAI	91	SexAI	1572
SfiI	7968	SfiI	8066	SmaI	347	SmaI	1056	SseBI	323	SseBI	1336	StuI	323	StuI	1336
TspMI	345	TspMI	1054	Vha464I	2532	Vha464I	5492	XmaI	345	XmaI	1054	XmaJI	324	XmaJI	550
Zrml	4478	Zrml	8166	Zsp2I	74	Zsp2I	146								

Table 4. Restriction Enzymes That Cleave TR207 Three times

Acc36I	738	Acc36I	7394	Acc36I	8097	Acc65I	840	Acc65I	2345	Acc65I	8053	AccI	422	AccI	4224
AccI	4597	Asp718I	840	Asp718I	2345	Asp718I	8053	BciVI	1687	BciVI	3311	BciVI	4835	BfuAI	738
BfuAI	7394	BfuAI	8097	BfuI	1687	BfuI	3311	BfuI	4835	BpvUI	4368	BpvUI	5320	BpvUI	7538
BseX3I	1453	BseX3I	6119	BseX3I	6123	Bsp19I	231	Bsp19I	2026	Bsp19I	7868	BspMI	738	BspMI	7394
BspMI	8097	BstZI	1453	BstZI	6119	BstZI	6123	BsuI	1687	BsuI	3311	BsuI	4835	BveI	738
BveI	7394	BveI	8097	Cfr42I	1219	Cfr42I	2121	Cfr42I	7741	DinI	1214	DinI	1345	DinI	5614
EagI	1453	EagI	6119	EagI	6123	EclXI	1453	EclXI	6119	EclXI	6123	Eco52I	1453	Eco52I	6119
Eco52I	6123	EgeI	1214	EgeI	1345	EgeI	5614	EheI	1214	EheI	1345	EheI	5614	KasI	1212
KasI	1343	KasI	5612	KpnI	844	KpnI	2349	KpnI	8057	KspI	1219	KspI	2121	KspI	7741
Mly113I	1213	Mly113I	1344	Mly113I	5613	MvrI	4368	MvrI	5320	MvrI	7538	NarI	1213	NarI	1344
NarI	5613	NcoI	231	NcoI	2026	NcoI	7868	Ple19I	4368	Ple19I	5320	Ple19I	7538	PluTI	1216
PluTI	1347	PluTI	5616	Psp1406I	422	Psp1406I	4224	Psp1406I	4597	PvuI	4368	PvuI	5320	PvuI	7538
SacII	1219	SacII	2121	SacII	7741	SfoI	1214	SfoI	1345	SfoI	5614	Sfr303I	1219	Sfr303I	2121
Sfr303I	7741	SgrBI	1219	SgrBI	2121	SgrBI	7741	SspDI	1212	SspDI	1343	SspDI	5612		