

Datasheet of GLuc-ONTM ERSE Transcriptional Response Element (TRE)
Clone (TR110)

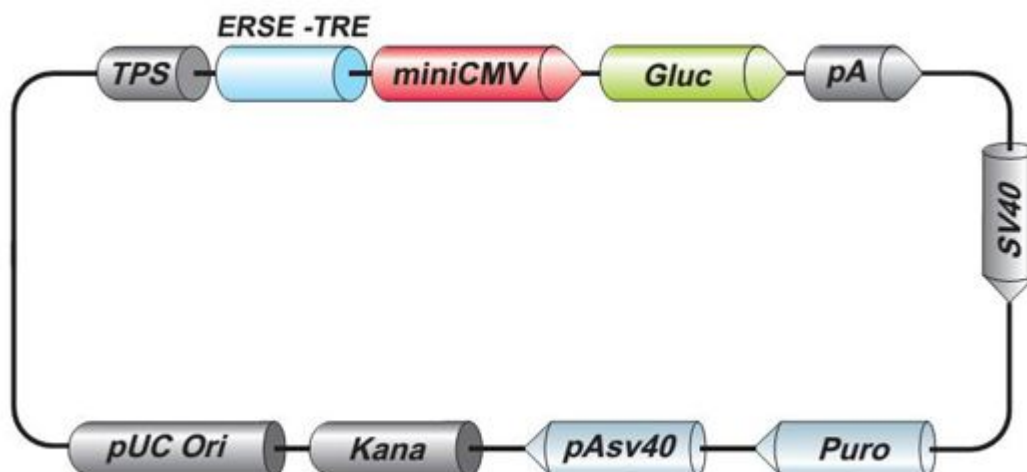


Figure 1. GLuc-ONTM ERSE Transcriptional Response Element (TRE) Clone(TR110)

Catalog:

TR110

Description:

The ERSE transcriptional response element (TRE) clone assays the status of the ER stress pathway through the activity of the transcription factor ERSE. The ERSE transcriptional TRE clone contains a minimal promoter and tandem repeats of the ERSE 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 × 40 μL

Storage Conditions:

-20°C

QC:

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

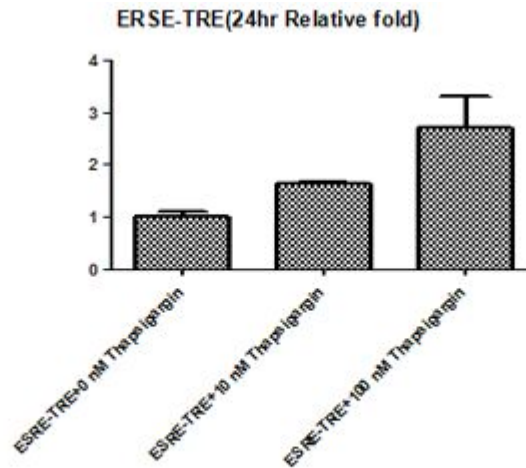


Figure 2. Validation data of GLuc-ON™ ERSE Transcriptional Response Element (TRE) Clone(TR110).

Restriction Enzyme Information for TR110

Table 1. Restriction Enzymes That Do Not Cleave TR110

AfeI	AgeI	AscI	BbvCI	BlpI	BsaAI	BstXI	BstZ17I
Bsu36I	BtrI	EcoRV	FspI	MluI	NdeI	PacI	PmeI
PmlI	PpuMI	PshAI	PvuII	SanDI	SapI	SbfI	SgrAI
SnaBI	SpeI	SrfI	Swal				

Table 2. Restriction Enzymes That Cleave TR110 Once

AarI	619	AasI	4636	Acc65I	663	AccI	1301	AccIII	2277	AcII	209	AcuI	4196	Adel	2735
AflII	343	AflIII	4738	AhdI	1266	Alol	991	Aor13HI	2277	AseI	3819	AsiSI	3620	Asp718I	663
AspA2I	2087	AspEI	1266	AspI	2206	AssI	554	AsuII	458	AvrII	2087	BaeI	2845	BamHI	951
BauI	4565	BciVI	4540	BclI	2138	BfrI	343	BfuI	4540	BlnI	2087	BmcAI	554	BmeRI	1266
Bpu14I	458	Bsa29I	547	BsaBI	994	BsaI	2638	Bse8I	994	BseAI	2277	BseCI	547	BseJI	994
BsePI	2619	BseRI	2083	BshVI	547	BsiWI	2220	Bso31I	2638	Bsp119I	458	Bsp13I	2277	BspDI	547
BspEI	2277	BspHI	3144	BspT104I	458	BspTI	343	BssHII	2619	BssSI	4565	Bst2BI	4565	BstAFI	343
BstBI	458	BstEII	2298	BstENI	3532	BstPI	2298	Bsu15I	547	BsuI	4540	CciI	3144	CciNI	1523
Cfr9I	2108	Clal	547	CpoI	2280	Cspl	2280	DraI	3044	DraIII	2735	DrdI	4636	Dril	1266
DseDI	4636	Eam1105I	1266	Ecl136II	672	Eco31I	2638	Eco53kl	672	Eco57I	4196	Eco91I	2298	EcoCRI	672
EcoNI	3532	EcoO65I	2298	EcoRI	483	FbaI	2138	FblI	1301	FseI	2842	HindIII	875	HpaI	2983
Kpn2I	2277	KpnI	667	KroI	2838	Ksp22I	2138	KspAI	2983	MfeI	2992	MroI	2277	MroNI	2838
MspCI	343	MunI	2992	NaeI	2840	NgoMIV	2838	NotI	1523	NspV	458	PaeR7I	1529	PagI	3144
PasI	2213	Paul	2619	PciI	4738	PdII	2840	Pfi23II	2220	PfiFI	2206	PfoI	2639	PscI	4738
PshBI	3819	Psp124BI	674	Psp1406I	209	PspEI	2298	PspLI	2220	PspXI	1529	PstI	3085	Psyl	2206
Ptel	2619	RcaI	3144	RgaI	3620	RigI	2842	Rsr2I	2280	RsrII	2280	SacI	674	Sall	1300
Scal	554	SfaAI	3620	Sfr274I	1529	SfuI	458	SgfI	3620	Slal	1529	Smal	2110	Sspl	3545
SstI	674	StrI	1529	TspMI	2108	Tth111I	2206	Vha464I	343	VspI	3819	XagI	3532	XbaI	1541
XcmI	1762	XhoI	1529	XmaI	2108	XmaJI	2087	XmiI	1301	Zrml	554				

Table 3. Restriction Enzymes That Cleave TR110 Twice

Acc36I	619	Acc36I	1436	AccB7I	479	AccB7I	3883	Ajul	462	Ajul	3648	Alw44I	1206	Alw44I	4424
AlwNI	1178	AlwNI	4329	Apal	356	Apal	4046	ApaLI	1206	ApaLI	4424	Asp700I	1225	Asp700I	1549
AsuNHI	44	AsuNHI	676	Ball	1057	Ball	2539	BasI	479	BasI	3883	BbsI	134	BbsI	1040
Bcgl	71	Bcgl	2605	BfuAI	619	BfuAI	1436	BgIII	475	BgIII	856	BmtI	48	BmtI	680
Bpil	134	Bpil	1040	BpII	2022	BpII	2054	Bpu10I	668	Bpu10I	3637	BpuAI	134	BpuAI	1040
BpvUI	1298	BpvUI	3620	BsaXI	989	BsaXI	2542	BsmBI	2556	BsmBI	3636	Bsp120I	352	Bsp120I	4042
Bsp1407I	440	Bsp1407I	1663	Bsp19I	962	Bsp19I	1994	Bsp68I	1069	Bsp68I	3277	BspMI	619	BspMI	1436
BspOI	48	BspOI	680	BsrGI	440	BsrGI	1663	BstAUI	440	BstAUI	1663	BstV2I	134	BstV2I	1040
BtgZI	981	BtgZI	1338	BtuMI	1069	BtuMI	3277	BveI	619	BveI	1436	Cail	1178	Cail	4329
Cfr42I	1095	Cfr42I	2378	Csil	1854	Csil	2731	DinI	2373	DinI	2504	Drall	353	Drall	4043
Eco147I	2086	Eco147I	2495	EcoO109I	353	EcoO109I	4043	EgeI	2373	EgeI	2504	EheI	2373	EheI	2504
Esp3I	2556	Esp3I	3636	HincII	1302	HincII	2983	HindII	1302	HindII	2983	KasI	2371	KasI	2502
KspI	1095	KspI	2378	MabI	1854	MabI	2731	MisI	1057	MisI	2539	MluNI	1057	MluNI	2539
Mly113I	2372	Mly113I	2503	MroXI	1225	MroXI	1549	MscI	1057	MscI	2539	MslI	524	MslI	1000
Msp20I	1057	Msp20I	2539	Mvrl	1298	Mvrl	3620	NarI	2372	NarI	2503	NcoI	962	NcoI	1994
NheI	44	NheI	676	Nrul	1069	Nrul	3277	PceI	2086	PceI	2495	Pdml	1225	Pdml	1549
PfiMI	479	PfiMI	3883	Ple19I	1298	Ple19I	3620	PluTI	2375	PluTI	2506	PspOMI	352	PspOMI	4042
PstNI	1178	PstNI	4329	PvuI	1298	PvuI	3620	Rrul	1069	Rrul	3277	RseI	524	RseI	1000
SacII	1095	SacII	2378	SexAI	1854	SexAI	2731	SfiI	657	SfiI	869	SfoI	2373	SfoI	2504
Sfr303I	1095	Sfr303I	2378	SgrBI	1095	SgrBI	2378	SmiMI	524	SmiMI	1000	SseBI	2086	SseBI	2495
SspDI	2371	SspDI	2502	StuI	2086	StuI	2495	Van91I	479	Van91I	3883	VneI	1206	VneI	4424
XmnI	1225	XmnI	1549												

Table 4. Restriction Enzymes That Cleave TR110 Three Times

AanI	427	AanI	1680	AanI	2963	AatII	66	AatII	2210	AatII	2708	Avall	2280	Avall	2390
Avall	2722	Bfml	3081	Bfml	4282	Bfml	4473	Bme18I	2280	Bme18I	2390	Bme18I	2722	Bmrl	141
Bmrl	611	Bmrl	1966	Bmul	141	Bmul	611	Bmul	1966	Bpml	1144	Bpml	1363	Bpml	2662
Bse3DI	421	Bse3DI	1688	Bse3DI	3118	BseMI	421	BseMI	1688	BseMI	3118	BsrDI	421	BsrDI	1688
BsrDI	3118	Bst6I	329	Bst6I	2323	Bst6I	3433	BstAPI	1832	BstAPI	1904	BstAPI	2735	BstSFI	3081
BstSFI	4282	BstSFI	4473	Eam1104I	329	Eam1104I	2323	Eam1104I	3433	EarI	329	EarI	2323	EarI	3433
Eco47I	2280	Eco47I	2390	Eco47I	2722	Gsul	1144	Gsul	1363	Gsul	2662	Psil	427	Psil	1680
Psil	2963	Sfcl	3081	Sfcl	4282	Sfcl	4473	TaqII	2182	TaqII	2806	TaqII	3874	VpaK11BI	2280
VpaK11BI	2390	VpaK11BI	2722	Zral	64	Zral	2208	Zral	2706						
