

Datasheet of GLuc-ONTM CREB Transcriptional Response Element (TRE)
Clone (TR100)

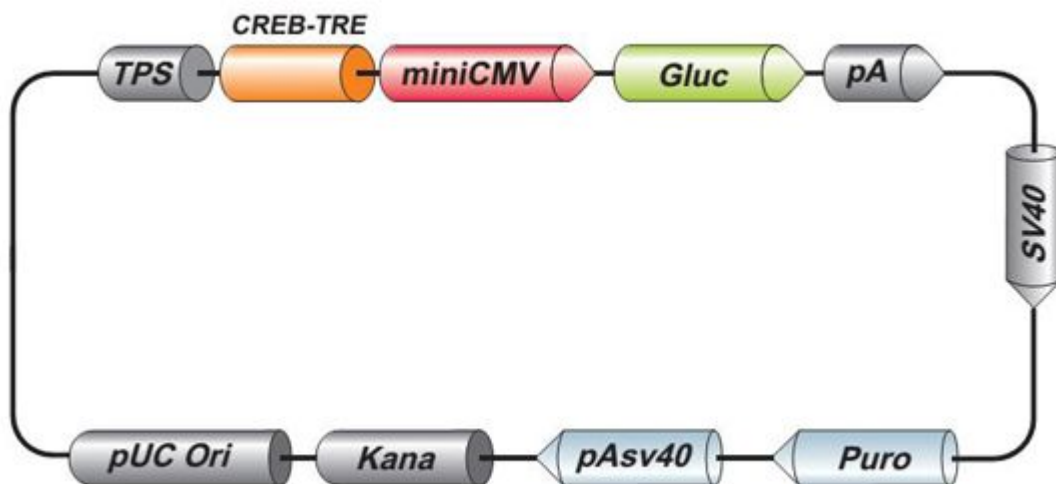


Figure 1. GLuc-ONTM CREB Transcriptional Response Element (TRE) Clone (TR100)

Catalog:

TR100

Description:

The CREB transcriptional response element (TRE) clone assays the status of the cAMP/PKA signaling pathway through the activity of the transcription factor CREB . The CREB transcriptional TRE clone contains a minimal promoter and tandem repeats of the CREB 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

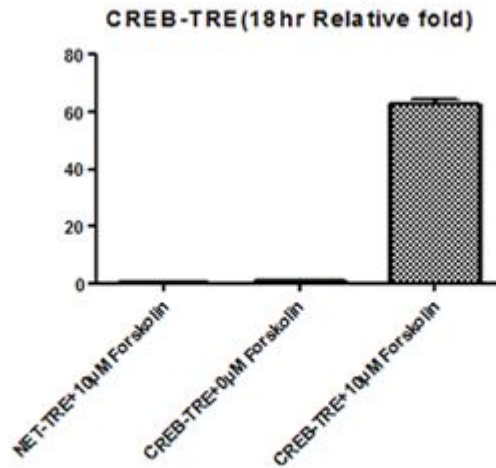


Figure 2. Validation data of GLuc-ON™ CREB Transcriptional Response Element (TRE) Clone(TR101).

Restriction Enzyme Information for TR100

Table 1. Restriction Enzymes That Do Not Cleave TR100

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

Table 2. Restriction Enzymes That Cleave TR100 Once

AarI	619	AccI	1214	AccIII	2190	AcII	209	AcuI	4110	Adel	2648	AflII	343	AflIII	4652
AhdI	1179	Alol	904	Aor13HI	2190	Asel	3733	AsiSI	3534	AspA2I	2000	AspEI	1179	AspI	2119
AssI	554	AsuII	458	AvrII	2000	BamHI	864	BauI	4479	BciVI	4454	BclI	2051	BfrI	343
BfuI	4454	BlnI	2000	BmcAI	554	BmeRI	1179	Bpu14I	458	Bsa29I	547	BsaBI	907	Bsal	2551
Bse8I	907	BseAI	2190	BseCI	547	BseJI	907	BsePI	2532	BseRI	1996	BshVI	547	BsiWI	2133
Bso31I	2551	Bsp119I	458	Bsp13I	2190	BspDI	547	BspEI	2190	BspHI	3058	BspT104I	458	BspTI	343
BssHII	2532	BssSI	4479	Bst2BI	4479	BstAFI	343	BstBI	458	BstEII	2211	BstENI	3446	BstPI	2211
Bsu15I	547	BsuI	4454	CciI	3058	CciNI	1436	Cfr9I	2021	Clal	547	Cpol	2193	Cspl	2193
DraI	2958	DraIII	2648	DriI	1179	Eam1105I	1179	Ecl136II	672	Eco31I	2551	Eco53kI	672	Eco57I	4110
Eco91I	2211	EcoICRI	672	EcoNI	3446	EcoO65I	2211	EcoRI	483	FbaI	2051	FblI	1214	FseI	2756
HindIII	788	HpaI	2897	Kpn2I	2190	KroI	2752	Ksp22I	2051	KspAI	2897	MfeI	2906	MroI	2190
MroNI	2752	MspCI	343	MunI	2906	NaeI	2754	NgoMIV	2752	NotI	1436	NspV	458	Paer7I	1442
PagI	3058	PasI	2126	Paul	2532	PciI	4652	PdII	2754	Pfl23II	2133	PflFI	2119	PfoI	2552
PscI	4652	PshBI	3733	Psp124BI	674	Psp1406I	209	PspEI	2211	PspLI	2133	PspXI	1442	PstI	2999
PsyI	2119	PteI	2532	PvuII	701	RcaI	3058	RgaI	3534	RglI	2756	Rsr2I	2193	RsrII	2193
SacI	674	Sall	1213	Scal	554	SfaAI	3534	Sfr274I	1442	SfuI	458	Sgfl	3534	Slal	1442
Smal	2023	SspI	3459	SstI	674	StrI	1442	TspMI	2021	Tth111I	2119	Vha464I	343	VspI	3733
XagI	3446	XbaI	1454	XcmI	1675	XhoI	1442	XmaI	2021	XmaJI	2000	XmiI	1214	Zrml	554

Table 3. Restriction Enzymes That Cleave TR100 Twice

AasI	694	AasI	4550	Acc36I	619	Acc36I	1349	Acc65I	663	Acc65I	2733	AccB7I	479	AccB7I	3797
AjuI	462	AjuI	3562	Alw44I	1119	Alw44I	4338	AlwNI	1091	AlwNI	4243	Apal	356	Apal	3960
ApaLI	1119	ApaLI	4338	Asp700I	1138	Asp700I	1462	Asp718I	663	Asp718I	2733	AsuNHI	44	AsuNHI	676
Ball	970	Ball	2452	BasI	479	BasI	3797	BbsI	134	BbsI	953	Bcgl	71	Bcgl	2518
BfuAI	619	BfuAI	1349	BglII	475	BglII	769	BmtI	48	BmtI	680	Bpil	134	Bpil	953
BpII	1935	BpII	1967	Bpu10I	668	Bpu10I	3551	BpuAI	134	BpuAI	953	BpvUI	1211	BpvUI	3534
BsaXI	902	BsaXI	2455	BsmBI	2469	BsmBI	3550	Bsp120I	352	Bsp120I	3956	Bsp1407I	440	Bsp1407I	1576
Bsp68I	982	Bsp68I	3191	BspMI	619	BspMI	1349	BspOI	48	BspOI	680	BsrGI	440	BsrGI	1576
BstAUI	440	BstAUI	1576	BstV2I	134	BstV2I	953	BtgZI	894	BtgZI	1251	BtuMI	982	BtuMI	3191
BveI	619	BveI	1349	Cail	1091	Cail	4243	Cfr42I	1008	Cfr42I	2291	Csil	1767	Csil	2644
DinI	2286	DinI	2417	Drall	353	Drall	3957	DrdI	694	DrdI	4550	DseDI	694	DseDI	4550
Eco147I	1999	Eco147I	2408	EcoO109I	353	EcoO109I	3957	EgeI	2286	EgeI	2417	Ehel	2286	Ehel	2417
Esp3I	2469	Esp3I	3550	HincII	1215	HincII	2897	HindII	1215	HindII	2897	KasI	2284	KasI	2415
KpnI	667	KpnI	2737	KspI	1008	KspI	2291	MabI	1767	MabI	2644	MisI	970	MisI	2452
MluNI	970	MluNI	2452	Mly113I	2285	Mly113I	2416	MroXI	1138	MroXI	1462	MscI	970	MscI	2452
MslI	524	MslI	913	Msp20I	970	Msp20I	2452	MvrI	1211	MvrI	3534	NarI	2285	NarI	2416
NheI	44	NheI	676	NruI	982	NruI	3191	PceI	1999	PceI	2408	Pdml	1138	Pdml	1462
PfiMI	479	PfiMI	3797	Ple19I	1211	Ple19I	3534	PluTI	2288	PluTI	2419	PspOMI	352	PspOMI	3956
PstNI	1091	PstNI	4243	PvuI	1211	PvuI	3534	RruI	982	RruI	3191	RseI	524	RseI	913
SacII	1008	SacII	2291	SexAI	1767	SexAI	2644	SfiI	657	SfiI	782	SfoI	2286	SfoI	2417
Sfr303I	1008	Sfr303I	2291	SgrBI	1008	SgrBI	2291	SmiMI	524	SmiMI	913	SseBI	1999	SseBI	2408
SspDI	2284	SspDI	2415	StuI	1999	StuI	2408	Van91I	479	Van91I	3797	VneI	1119	VneI	4338
XmnI	1138	XmnI	1462												

Table 4. Restriction Enzymes That Cleave TR100 Three Times

AanI	427	AanI	1593	AanI	2877	Avall	2193	Avall	2303	Avall	2635	Bfml	2995	Bfml	4196
Bfml	4387	Bme18I	2193	Bme18I	2303	Bme18I	2635	Bmrl	141	Bmrl	611	Bmrl	1879	Bmul	141
Bmul	611	Bmul	1879	Bpml	1057	Bpml	1276	Bpml	2575	Bse3DI	421	Bse3DI	1601	Bse3DI	3032
BseMI	421	BseMI	1601	BseMI	3032	Bsp19I	712	Bsp19I	875	Bsp19I	1907	BsrDI	421	BsrDI	1601
BsrDI	3032	Bst6I	329	Bst6I	2236	Bst6I	3347	BstAPI	1745	BstAPI	1817	BstAPI	2648	BstSFI	2995
BstSFI	4196	BstSFI	4387	Eam1104I	329	Eam1104I	2236	Eam1104I	3347	EarI	329	EarI	2236	EarI	3347
Eco47I	2193	Eco47I	2303	Eco47I	2635	Gsul	1057	Gsul	1276	Gsul	2575	NcoI	712	NcoI	875
NcoI	1907	Psil	427	Psil	1593	Psil	2877	Sfcl	2995	Sfcl	4196	Sfcl	4387	TaqII	2095
TaqII	2719	TaqII	3788	VpaK11BI	2193	VpaK11BI	2303	VpaK11BI	2635						
