

Datasheet of GLuc-ONTM TCF/LEF Transcriptional Response Element (TRE)
Clone (TR102)

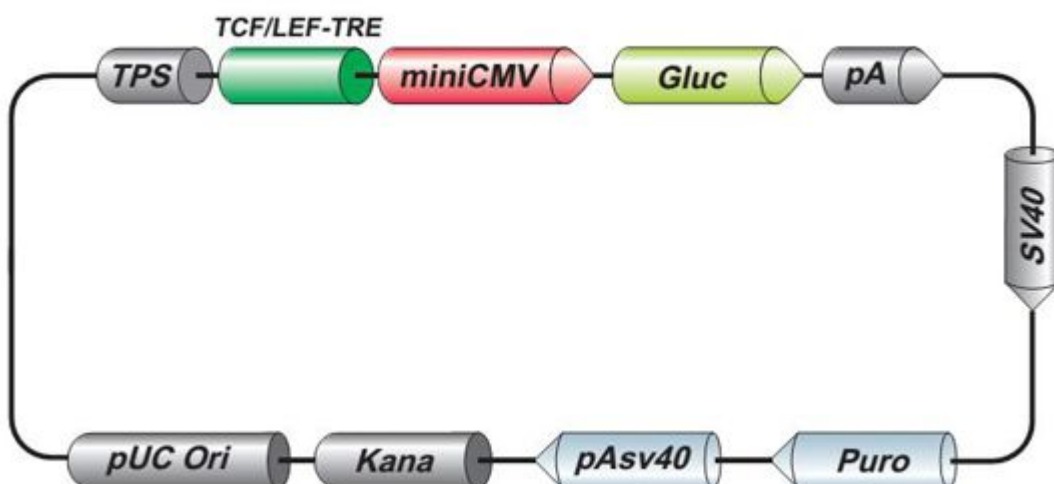


Figure 1. GLuc-ONTM TCF/LEF Transcriptional Response Element (TRE) Clone (TR102)

Catalog:

TR102

Description:

The TCF/LEF transcriptional response element (TRE) 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 × 40 μL

Storage Conditions:

-20°C

QC:

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

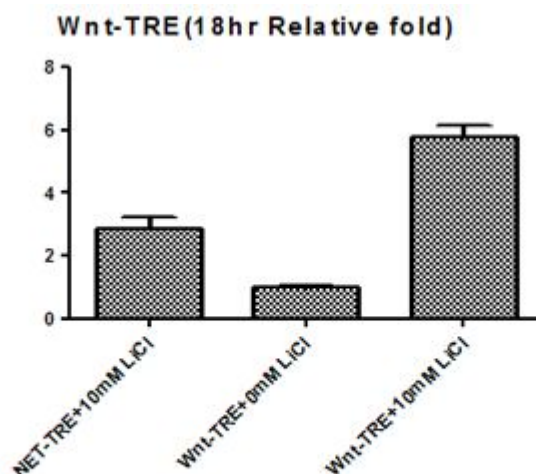


Figure 2. Validation data of GLuc-ON™ TCF/LEF Transcriptional Response Element (TRE) Clone(TR102).

Restriction Enzyme Information for TR102

Table 1. Restriction Enzymes That Do Not Cleave TR102

AfeI	AgeI	AscI	BaeI	BaeI	BbvCI	BlnI	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 TR102 Once

AarI	619	AasI	4553	AccI	1217	AccIII	2193	AcII	209	AcuI	4113	Adel	2651	AflIII	4655
AhdI	1182	Alol	907	Aor13HI	2193	Asel	3736	AsiSI	3537	AspA2I	2003	AspEI	1182	AspI	2122
AssI	554	AsuII	458	AvrII	2003	BamHI	867	Baul	4482	BciVI	4457	BclI	2054	BfuI	4457
BlnI	2003	BmcAI	554	BmeRI	1182	Bpu14I	458	Bsa29I	547	BsaBI	910	BsaI	2554	Bse8I	910
BseAI	2193	BseCI	547	BseJI	910	BsePI	2535	BseRI	1999	BshVI	547	BsiWI	2136	Bso31I	2554
Bsp119I	458	Bsp13I	2193	BspDI	547	BspEI	2193	BspHI	3061	BspT104I	458	BssHII	2535	BssSI	4482
Bst2BI	4482	BstBI	458	BstEII	2214	BstENI	3449	BstPI	2214	Bsu15I	547	BsuI	4457	CciI	3061
CciNI	1439	Cfr9I	2024	Clal	547	CpoI	2196	Cspl	2196	Dral	2961	DrallI	2651	DrdI	4553
Dril	1182	DseDI	4553	Eam1105I	1182	Ecl136II	672	Eco31I	2554	Eco53KI	672	Eco57I	4113	Eco91I	2214
EcoICRI	672	EcoNI	3449	EcoO65I	2214	EcoRI	483	FbaI	2054	FblI	1217	FseI	2759	HindIII	791
HpaI	2900	Kpn2I	2193	KroI	2755	Ksp22I	2054	KspAI	2900	MfeI	2909	MroI	2193	MroNI	2755
MunI	2909	Nael	2757	NgoMIV	2755	NotI	1439	NspV	458	Paer7I	1445	PagI	3061	PasI	2129
Paul	2535	PciI	4655	PdII	2757	Pfl23II	2136	PfIFI	2122	PfoI	2555	PscI	4655	PshBI	3736
Psp124BI	674	Psp1406I	209	PspEI	2214	PspLI	2136	PspXI	1445	PstI	3002	PsyI	2122	PteI	2535
RcaI	3061	RgaI	3537	RigI	2759	Rsr2I	2196	RsrII	2196	SacI	674	Sall	1216	Scal	554
SfaAI	3537	Sfr274I	1445	SfuI	458	Sgfl	3537	Slal	1445	SmaI	2026	Sspl	3462	SstI	674
StrI	1445	TspMI	2024	Tth111I	2122	VspI	3736	XagI	3449	XbaI	1457	XcmI	1678	XhoI	1445
XmaI	2024	XmaJI	2003	XmII	1217	Zrml	554								

Table 3. Restriction Enzymes That Cleave TR102 Twice

Acc36I	619	Acc36I	1352	Acc65I	663	Acc65I	2736	AccB7I	479	AccB7I	3800	AfIII	343	AfIII	768
AjuI	462	AjuI	3565	Alw44I	1122	Alw44I	4341	AlwNI	1094	AlwNI	4246	ApaLI	1122	ApaLI	4341
Asp700I	1141	Asp700I	1465	Asp718I	663	Asp718I	2736	AsuNHI	44	AsuNHI	676	Ball	973	Ball	2455
BasI	479	BasI	3800	BbsI	134	BbsI	956	BcgI	71	BcgI	2521	BfRI	343	BfRI	768
BfuAI	619	BfuAI	1352	BglII	475	BglII	772	BmtI	48	BmtI	680	BpII	134	BpII	956
BpII	1938	BpII	1970	Bpu10I	668	Bpu10I	3554	BpuAI	134	BpuAI	956	BpvUI	1214	BpvUI	3537
BsaXI	905	BsaXI	2458	BsmBI	2472	BsmBI	3553	Bsp1407I	440	Bsp1407I	1579	Bsp19I	878	Bsp19I	1910
Bsp68I	985	Bsp68I	3194	BspMI	619	BspMI	1352	BspOI	48	BspOI	680	BspTI	343	BspTI	768
BsrGI	440	BsrGI	1579	BstAFI	343	BstAFI	768	BstAUI	440	BstAUI	1579	BstV2I	134	BstV2I	956
BtgZI	897	BtgZI	1254	BtuMI	985	BtuMI	3194	BveI	619	BveI	1352	CaiI	1094	CaiI	4246
Cfr42I	1011	Cfr42I	2294	CsiI	1770	CsiI	2647	DinI	2289	DinI	2420	Eco147I	2002	Eco147I	2411
EgeI	2289	EgeI	2420	EheI	2289	EheI	2420	Esp3I	2472	Esp3I	3553	HincII	1218	HincII	2900
HindII	1218	HindII	2900	KasI	2287	KasI	2418	KpnI	667	KpnI	2740	KspI	1011	KspI	2294
MabI	1770	MabI	2647	MisI	973	MisI	2455	MluNI	973	MluNI	2455	Mly113I	2288	Mly113I	2419
MroXI	1141	MroXI	1465	MscI	973	MscI	2455	MslI	524	MslI	916	Msp20I	973	Msp20I	2455
MspCI	343	MspCI	768	MvrI	1214	MvrI	3537	NarI	2288	NarI	2419	NcoI	878	NcoI	1910
NheI	44	NheI	676	NruI	985	NruI	3194	PceI	2002	PceI	2411	PdmI	1141	PdmI	1465
PfIMI	479	PfIMI	3800	Ple19I	1214	Ple19I	3537	PluTI	2291	PluTI	2422	PstNI	1094	PstNI	4246
PvuI	1214	PvuI	3537	RruI	985	RruI	3194	RseI	524	RseI	916	SacII	1011	SacII	2294
SexAI	1770	SexAI	2647	SfiI	657	SfiI	785	SfoI	2289	SfoI	2420	Sfr303I	1011	Sfr303I	2294
SgrBI	1011	SgrBI	2294	SmiMI	524	SmiMI	916	SseBI	2002	SseBI	2411	SspDI	2287	SspDI	2418
StuI	2002	StuI	2411	Van91I	479	Van91I	3800	Vha464I	343	Vha464I	768	VneI	1122	VneI	4341
XmnI	1141	XmnI	1465												

Table 4. Restriction Enzymes That Cleave TR102 Three Times

AanI	427	AanI	1596	AanI	2880	AatII	66	AatII	2126	AatII	2624	Apal	356	Apal	727
Apal	3963	Avall	2196	Avall	2306	Avall	2638	Bfml	2998	Bfml	4199	Bfml	4390	Bme18I	2196
Bme18I	2306	Bme18I	2638	Bmrl	141	Bmrl	611	Bmrl	1882	Bmul	141	Bmul	611	Bmul	1882
Bpml	1060	Bpml	1279	Bpml	2578	Bse3DI	421	Bse3DI	1604	Bse3DI	3035	BseMI	421	BseMI	1604
BseMI	3035	Bsp120I	352	Bsp120I	723	Bsp120I	3959	BsrDI	421	BsrDI	1604	BsrDI	3035	Bst6I	329
Bst6I	2239	Bst6I	3350	BstAPI	1748	BstAPI	1820	BstAPI	2651	BstSFI	2998	BstSFI	4199	BstSFI	4390
Eam1104I	329	Eam1104I	2239	Eam1104I	3350	EarI	329	EarI	2239	EarI	3350	Eco47I	2196	Eco47I	2306
Eco47I	2638	Gsul	1060	Gsul	1279	Gsul	2578	Psil	427	Psil	1596	Psil	2880	PspOMI	352
PspOMI	723	PspOMI	3959	Sfcl	2998	Sfcl	4199	Sfcl	4390	TaqII	2098	TaqII	2722	TaqII	3791
VpaK11BI	2196	VpaK11BI	2306	VpaK11BI	2638	Zral	64	Zral	2124	Zral	2622				
