

Datasheet of GLuc-ONTM Oct4 Transcriptional Response Element (TRE)
Clone (TR103)

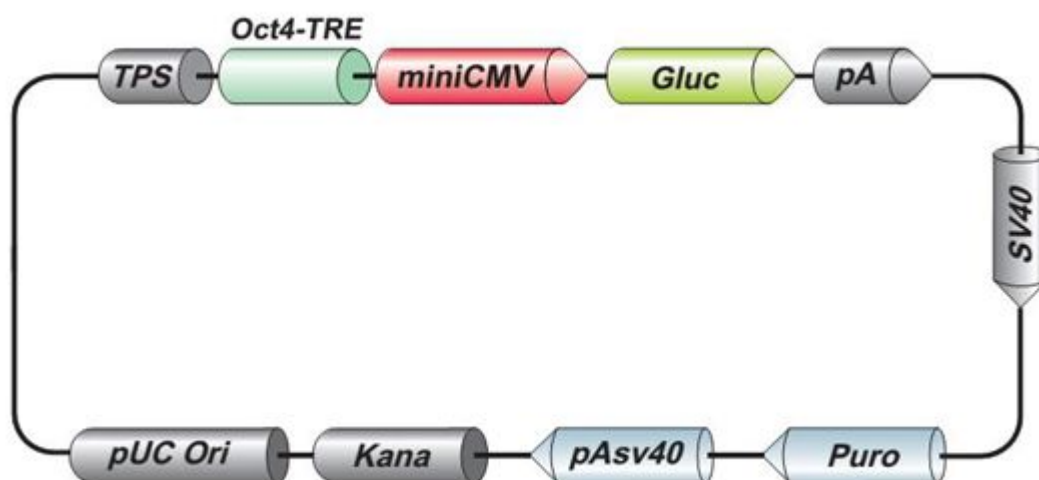


Figure 1. GLuc-ONTM Oct4 Transcriptional Response Element (TRE) Clone (TR103)

Catalog:

TR103

Description:

The Oct4 transcriptional response element (TRE) clone assays the status of the Oct4 signaling pathway through the activity of the transcription factor Oct4. The Oct4 transcriptional TRE clone contains a minimal promoter and tandem repeats of the Oct4 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

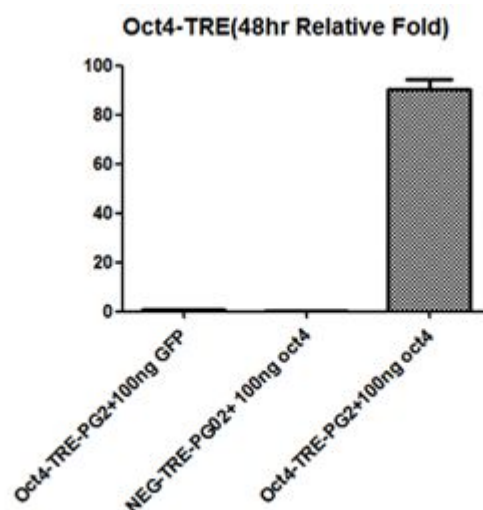


Figure 2. Validation data of GLuc-ON™ Oct4 Transcriptional Response Element (TRE) Clone(TR103).

Restriction Enzyme Information for TR103

Table 1. Restriction Enzymes That Do Not Cleave TR103

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 TR103 Once

AarI	619	AasI	4523	AccI	1187	AccIII	2163	AclI	209	AcuI	4083	Adel	2621	AflII	343
AflIII	4625	AhdI	1152	Alol	877	Aor13HI	2163	AseI	3706	AsiSI	3507	AspA2I	1973	AspEI	1152
AspI	2092	AssI	554	AsuII	458	AvrII	1973	BamHI	837	BauI	4452	BciVI	4427	BclI	2024
BfrI	343	BfuI	4427	BlnI	1973	BmcAI	554	BmeRI	1152	Bpu14I	458	Bsa29I	547	BsaBI	880
BsaI	2524	Bse8I	880	BseAI	2163	BseCI	547	BseJI	880	BsePI	2505	BseRI	1969	BshVI	547
BsiWI	2106	Bso31I	2524	Bsp119I	458	Bsp13I	2163	BspDI	547	BspEI	2163	BspHI	3031	BspT104I	458
BspTI	343	BssHII	2505	BssSI	4452	Bst2BI	4452	BstAFI	343	BstBI	458	BstEII	2184	BstENI	3419
BstPI	2184	Bsu15I	547	Bsul	4427	CciI	3031	CciNI	1409	Cfr9I	1994	ClaI	547	CpoI	2166
CsPI	2166	DraI	2931	DraIII	2621	DrdI	4523	DriI	1152	DseDI	4523	Eam1105I	1152	Ecl136II	672
Eco31I	2524	Eco53kI	672	Eco57I	4083	Eco91I	2184	EcoCRI	672	EcoNI	3419	EcoO65I	2184	EcoRI	483
FbaI	2024	FblI	1187	FseI	2729	HindIII	761	HpaI	2870	Kpn2I	2163	KroI	2725	Ksp22I	2024
KspAI	2870	MfeI	2879	MroI	2163	MroNI	2725	MspCI	343	MunI	2879	NaeI	2727	NgoMIV	2725
NotI	1409	NspV	458	PaeR7I	1415	PagI	3031	PasI	2099	Paul	2505	PciI	4625	PdII	2727
Pfi23II	2106	PfiFI	2092	PfoI	2525	PscI	4625	PshBI	3706	Psp124BI	674	Psp1406I	209	PspEI	2184
PspLI	2106	PspXI	1415	PstI	2972	PsyI	2092	PteI	2505	RcaI	3031	RgaI	3507	RigI	2729
Rsr2I	2166	RsrII	2166	SacI	674	Sall	1186	Scal	554	SfaAI	3507	Sfr274I	1415	SfuI	458
SgfI	3507	Slal	1415	SmaI	1996	SspI	3432	SstI	674	StrI	1415	TspMI	1994	Tth111I	2092
Vha464I	343	VspI	3706	XagI	3419	XbaI	1427	XcmI	1648	XhoI	1415	XmaI	1994	XmaJI	1973
XmiI	1187	Zrml	554												

Table 3. Restriction Enzymes That Cleave TR103 Twice

Acc36I	619	Acc36I	1322	Acc65I	663	Acc65I	2706	AccB7I	479	AccB7I	3770	Ajul	462	Ajul	3535
Alw44I	1092	Alw44I	4311	AlwNI	1064	AlwNI	4216	Apal	356	Apal	3933	ApaLI	1092	ApaLI	4311
Asp700I	1111	Asp700I	1435	Asp718I	663	Asp718I	2706	AsuNHI	44	AsuNHI	676	Ball	943	Ball	2425
BasI	479	BasI	3770	BbsI	134	BbsI	926	Bcgl	71	Bcgl	2491	BfuAI	619	BfuAI	1322
BglII	475	BglII	742	BmtI	48	BmtI	680	Bpil	134	Bpil	926	BpII	1908	BpII	1940
Bpu10I	668	Bpu10I	3524	BpuAI	134	BpuAI	926	BpvUI	1184	BpvUI	3507	BsaXI	875	BsaXI	2428
BsmBI	2442	BsmBI	3523	Bsp120I	352	Bsp120I	3929	Bsp1407I	440	Bsp1407I	1549	Bsp19I	848	Bsp19I	1880
Bsp68I	955	Bsp68I	3164	BspMI	619	BspMI	1322	BspOI	48	BspOI	680	BsrGI	440	BsrGI	1549
BstAUI	440	BstAUI	1549	BstV2I	134	BstV2I	926	BtgZI	867	BtgZI	1224	BtuMI	955	BtuMI	3164
BveI	619	BveI	1322	Cail	1064	Cail	4216	Cfr42I	981	Cfr42I	2264	Csil	1740	Csil	2617
DinI	2259	DinI	2390	Drall	353	Drall	3930	Eco147I	1972	Eco147I	2381	EcoO109I	353	EcoO109I	3930
EgeI	2259	EgeI	2390	Ehel	2259	Ehel	2390	Esp3I	2442	Esp3I	3523	HincII	1188	HincII	2870
HindII	1188	HindII	2870	KasI	2257	KasI	2388	KpnI	667	KpnI	2710	KspI	981	KspI	2264
MabI	1740	MabI	2617	MlsI	943	MlsI	2425	MluNI	943	MluNI	2425	Mly113I	2258	Mly113I	2389
MroXI	1111	MroXI	1435	MscI	943	MscI	2425	MslI	524	MslI	886	Msp20I	943	Msp20I	2425
MvrI	1184	MvrI	3507	NarI	2258	NarI	2389	NcoI	848	NcoI	1880	NheI	44	NheI	676
NruI	955	NruI	3164	PceI	1972	PceI	2381	Pdml	1111	Pdml	1435	PfiMI	479	PfiMI	3770
Ple19I	1184	Ple19I	3507	PluTI	2261	PluTI	2392	PspOMI	352	PspOMI	3929	PstNI	1064	PstNI	4216
PvuI	1184	PvuI	3507	RruI	955	RruI	3164	RseI	524	RseI	886	SacII	981	SacII	2264
SexAI	1740	SexAI	2617	SfiI	657	SfiI	755	SfoI	2259	SfoI	2390	Sfr303I	981	Sfr303I	2264
SgrBI	981	SgrBI	2264	SmiMI	524	SmiMI	886	SseBI	1972	SseBI	2381	SspDI	2257	SspDI	2388
StuI	1972	StuI	2381	Van91I	479	Van91I	3770	VneI	1092	VneI	4311	XmnI	1111	XmnI	1435

Table 4. Restriction Enzymes That Cleave TR103 Three Times

AanI	427	AanI	1566	AanI	2850	AatII	66	AatII	2096	AatII	2594	Avall	2166	Avall	2276
Avall	2608	BfmI	2968	BfmI	4169	BfmI	4360	Bme18I	2166	Bme18I	2276	Bme18I	2608	Bmrl	141
Bmrl	611	Bmrl	1852	Bmul	141	Bmul	611	Bmul	1852	Bpml	1030	Bpml	1249	Bpml	2548
Bse3DI	421	Bse3DI	1574	Bse3DI	3005	BseMI	421	BseMI	1574	BseMI	3005	BsrDI	421	BsrDI	1574
BsrDI	3005	Bst6I	329	Bst6I	2209	Bst6I	3320	BstAPI	1718	BstAPI	1790	BstAPI	2621	BstSFI	2968
BstSFI	4169	BstSFI	4360	Eam1104I	329	Eam1104I	2209	Eam1104I	3320	EarI	329	EarI	2209	EarI	3320
Eco47I	2166	Eco47I	2276	Eco47I	2608	Gsul	1030	Gsul	1249	Gsul	2548	Psil	427	Psil	1566
Psil	2850	Sfcl	2968	Sfcl	4169	Sfcl	4360	TaqII	2068	TaqII	2692	TaqII	3761	VpaK11BI	2166
VpaK11BI	2276	VpaK11BI	2608	Zral	64	Zral	2094	Zral	2592						
