

Datasheet of GLuc-ONTM NFAT Transcriptional Response Element (TRE)
Clone (TR101)

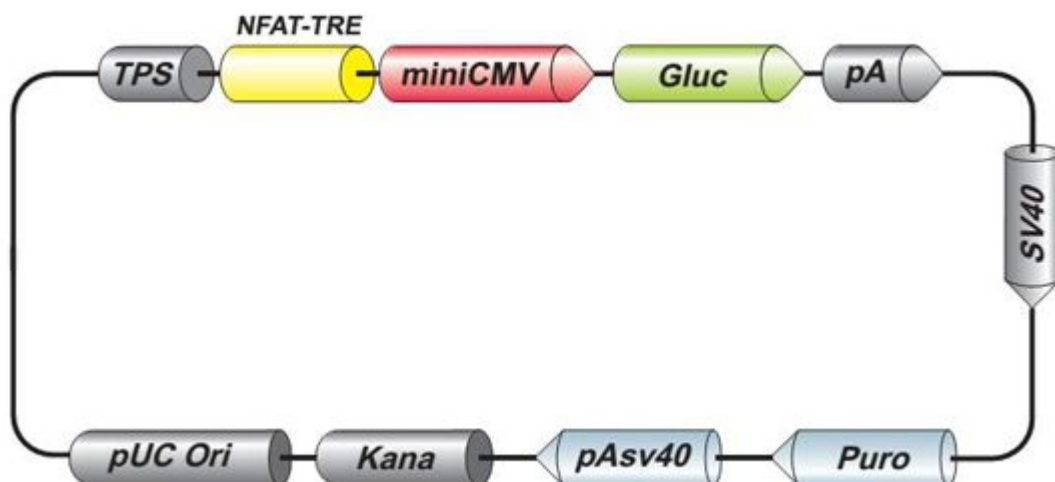


Figure 1. GLuc-ONTM NFAT Transcriptional Response Element (TRE) Clone (TR101)

Catalog:

TR101

Description:

The NFAT transcriptional response element (TRE) clone assays the status of the PKC/Ca⁺⁺ signaling pathway through the activity of the transcription factor NFAT. The NFAT transcriptional TRE clone contains a minimal promoter and tandem repeats of the NFAT 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:

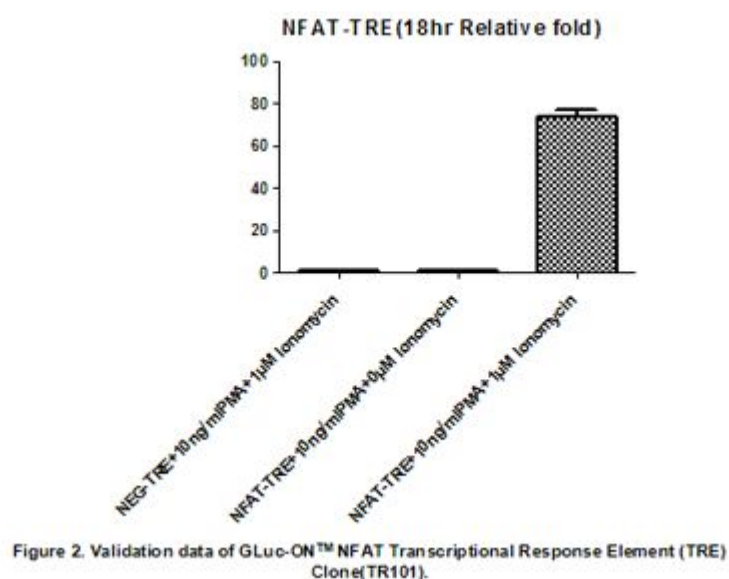
500ng/μL × 40 μL

Storage Conditions:

-20°C

QC:

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



Restriction Enzyme Information for TR101

Table 1. Restriction Enzymes That Do Not Cleave TR101

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 TR101 Once

AarI	619	AasI	4563	AccI	1227	AccIII	2203	AcII	209	AcuI	4123	Adel	2661	AflII	343
AflIII	4665	AhdI	1192	Alol	917	Aor13HI	2203	Asel	3746	AsiSI	3547	AspA2I	2013	AspEI	1192
AspI	2132	AssI	554	AsuII	458	AvrII	2013	BamHI	877	BauI	4492	BciVI	4467	BclI	2064
BfrI	343	BfuI	4467	BlnI	2013	BmcAI	554	BmeRI	1192	Bpu14I	458	Bsa29I	547	BsaBI	920
BsaI	2564	Bse8I	920	BseAI	2203	BseCI	547	BseJI	920	BsePI	2545	BseRI	2009	BshVI	547
BsiWI	2146	Bso31I	2564	Bsp119I	458	Bsp13I	2203	BspDI	547	BspEI	2203	BspHI	3071	BspT104I	458
BspTI	343	BssHII	2545	BssSI	4492	Bst2BI	4492	BstAFI	343	BstBI	458	BstEII	2224	BstENI	3459
BstPI	2224	Bsu15I	547	Bsul	4467	CciI	3071	CciNI	1449	Cfr9I	2034	Clal	547	Cpol	2206
CsPI	2206	DraI	2971	DraIII	2661	DrdI	4563	Dril	1192	DseDI	4563	Eam1105I	1192	Ecl136II	672
Eco31I	2564	Eco53kI	672	Eco57I	4123	Eco91I	2224	EcoICRI	672	EcoNI	3459	EcoO65I	2224	FbaI	2064
FblI	1227	FseI	2769	HindIII	842	HpaI	2910	Kpn2I	2203	KroI	2765	Ksp22I	2064	KspAI	2910
MfeI	2919	MroI	2203	MroNI	2765	MspCI	343	MunI	2919	NaeI	2767	NgoMIV	2765	NotI	1449
NspV	458	PaeR7I	1455	PagI	3071	PasI	2139	Paul	2545	PciI	4665	PdII	2767	Pfl23II	2146
PfiFI	2132	PfoI	2565	PscI	4665	PshBI	3746	Psp124BI	674	Psp1406I	209	PspEI	2224	PspLI	2146
PspXI	1455	PstI	3012	PsyI	2132	Ptel	2545	RcaI	3071	RgaI	3547	RigI	2769	Rsr2I	2206
RsrII	2206	SacI	674	Sall	1226	Scal	554	SfaAI	3547	SfiI	657	Sfr274I	1455	SfuI	458
SgfI	3547	SlaI	1455	SmaI	2036	SspI	3472	SstI	674	StrI	1455	TspMI	2034	Tth111I	2132
Vha464I	343	VspI	3746	XagI	3459	XcmI	1688	XhoI	1455	XmaI	2034	XmaJI	2013	XmiI	1227
Zrml	554														

Table 3. Restriction Enzymes That Cleave TR101 Twice

Acc36I	619	Acc36I	1362	Acc65I	663	Acc65I	2746	AccB7I	479	AccB7I	3810	Alw44I	1132	Alw44I	4351
AlwNI	1104	AlwNI	4256	Apal	356	Apal	3973	ApalI	1132	ApalI	4351	Asp700I	1151	Asp700I	1475
Asp718I	663	Asp718I	2746	AsuNHI	44	AsuNHI	676	Ball	983	Ball	2465	BasI	479	BasI	3810
BbsI	134	BbsI	966	Bcgl	71	Bcgl	2531	BfuAI	619	BfuAI	1362	BglII	475	BglII	772
BmtI	48	BmtI	680	Bpil	134	Bpil	966	BpII	1948	BpII	1980	Bpu10I	668	Bpu10I	3564
BpuAI	134	BpuAI	966	BpvUI	1224	BpvUI	3547	BsaXI	915	BsaXI	2468	BsmBI	2482	BsmBI	3563
Bsp120I	352	Bsp120I	3969	Bsp1407I	440	Bsp1407I	1589	Bsp19I	888	Bsp19I	1920	Bsp68I	995	Bsp68I	3204
BspMI	619	BspMI	1362	BspOI	48	BspOI	680	BsrGI	440	BsrGI	1589	BstAUI	440	BstAUI	1589
BstV2I	134	BstV2I	966	BtgZI	907	BtgZI	1264	BtuMI	995	BtuMI	3204	BveI	619	BveI	1362
Cail	1104	Cail	4256	Cfr42I	1021	Cfr42I	2304	Csil	1780	Csil	2657	DinI	2299	DinI	2430
Drall	353	Drall	3970	Eco147I	2012	Eco147I	2421	EcoO109I	353	EcoO109I	3970	EcoRI	483	EcoRI	807
Egel	2299	Egel	2430	Ehel	2299	Ehel	2430	Esp3I	2482	Esp3I	3563	HincII	1228	HincII	2910
HindII	1228	HindII	2910	KasI	2297	KasI	2428	KpnI	667	KpnI	2750	KspI	1021	KspI	2304
MabI	1780	MabI	2657	MisI	983	MisI	2465	MluNI	983	MluNI	2465	Mly113I	2298	Mly113I	2429
MroXI	1151	MroXI	1475	MscI	983	MscI	2465	MslI	524	MslI	926	Msp20I	983	Msp20I	2465
Mvrl	1224	Mvrl	3547	NarI	2298	NarI	2429	NcoI	888	NcoI	1920	NheI	44	NheI	676
Nrul	995	Nrul	3204	PceI	2012	PceI	2421	Pdml	1151	Pdml	1475	PfiMI	479	PfiMI	3810
Ple19I	1224	Ple19I	3547	PluTI	2301	PluTI	2432	PspOMI	352	PspOMI	3969	PstNI	1104	PstNI	4256
Pvul	1224	Pvul	3547	Rrul	995	Rrul	3204	RseI	524	RseI	926	SacII	1021	SacII	2304
SexAI	1780	SexAI	2657	SfoI	2299	SfoI	2430	Sfr303I	1021	Sfr303I	2304	SgrBI	1021	SgrBI	2304
SmiMI	524	SmiMI	926	SseBI	2012	SseBI	2421	SspDI	2297	SspDI	2428	StuI	2012	StuI	2421
Van91I	479	Van91I	3810	VneI	1132	VneI	4351	XmnI	1151	XmnI	1475				

Table 4. Restriction Enzymes That Cleave TR101 Three Times

AanI	427	AanI	1606	AanI	2890	AatII	66	AatII	2136	AatII	2634	AjuI	462	AjuI	799
AjuI	3575	Avall	2206	Avall	2316	Avall	2648	Bfml	3008	Bfml	4209	Bfml	4400	BglI	657
BglI	2306	BglI	2437	Bme18I	2206	Bme18I	2316	Bme18I	2648	Bmrl	141	Bmrl	611	Bmrl	1892
Bmul	141	Bmul	611	Bmul	1892	Bpml	1070	Bpml	1289	Bpml	2588	Bse3DI	421	Bse3DI	1614
Bse3DI	3045	BseMI	421	BseMI	1614	BseMI	3045	BsrDI	421	BsrDI	1614	BsrDI	3045	Bst6I	329
Bst6I	2249	Bst6I	3360	BstAPI	1758	BstAPI	1830	BstAPI	2661	BstSFI	3008	BstSFI	4209	BstSFI	4400
Eam1104I	329	Eam1104I	2249	Eam1104I	3360	EarI	329	EarI	2249	EarI	3360	Eco47I	2206	Eco47I	2316
Eco47I	2648	Gsul	1070	Gsul	1289	Gsul	2588	Psil	427	Psil	1606	Psil	2890	Sfcl	3008
Sfcl	4209	Sfcl	4400	TaqII	2108	TaqII	2732	TaqII	3801	VpaK11BI	2206	VpaK11BI	2316	VpaK11BI	2648
XbaI	775	XbaI	782	XbaI	1467	Zral	64	Zral	2134	Zral	2632				
