

Datasheet of GLuc-ONTM MEF2 Transcriptional Response Element (TRE)
Clone (TR109)

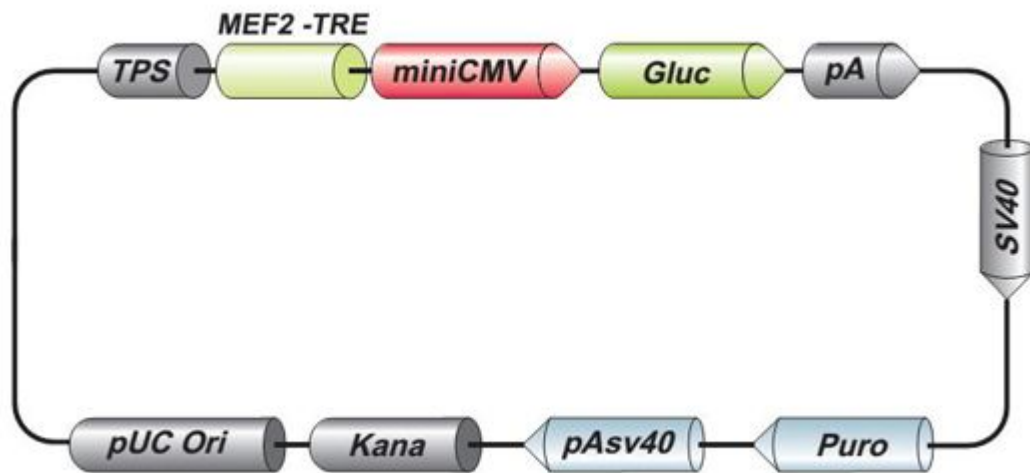


Figure 1. GLuc-ONTM MEF2 Transcriptional Response Element (TRE) Clone(TR109)

Catalog:	TR109
Description:	The MEF2 transcriptional response element (TRE) clone assays the status of the MEF2 signaling pathway through the activity of the transcription factor MEF2. The MEF2 transcriptional TRE clone contains a minimal promoter and tandem repeats of the MEF2 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

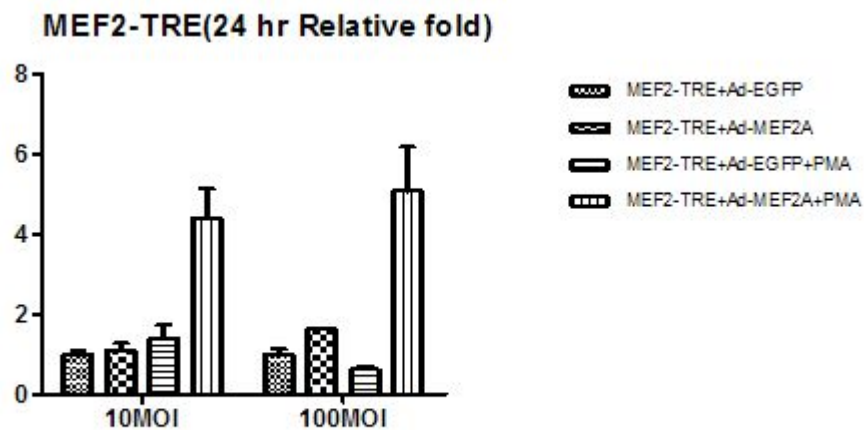


Figure 2. Validation data of GLuc-ON™ MEF2 Transcriptional Response Element (TRE) Clone(TR109).

Restriction Enzyme Information for TR109

Table 1. Restriction Enzymes That Do Not Cleave TR109

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	SwI					

Table 2. Restriction Enzymes That Cleave TR109 Once

AarI	619	AasI	4594	Acc65I	663	AccI	1259	AccIII	2235	AcII	209	AcuI	4154	Adel	2693
AflII	343	AflIII	4696	AhdI	1224	Alol	949	Aor13HI	2235	AseI	3777	AsiSI	3578	Asp718I	663
AspA2I	2045	AspEI	1224	AspI	2164	AssI	554	AsuII	458	AvrII	2045	BaeI	2803	BamHI	909
BauI	4523	BciVI	4498	BclI	2096	BfrI	343	BfuI	4498	BlnI	2045	BmcAI	554	BmeRI	1224
Bpu14I	458	Bsa29I	547	BsaBI	952	BsaI	2596	Bse8I	952	BseAI	2235	BseCI	547	BseJI	952
BsePI	2577	BseRI	2041	BshVI	547	BsiWI	2178	Bso31I	2596	Bsp119I	458	Bsp13I	2235	BspDI	547
BspEI	2235	BspHI	3102	BspT104I	458	BspTI	343	BssHII	2577	BssSI	4523	Bst2BI	4523	BstAFI	343
BstBI	458	BstEII	2256	BstENI	3490	BstPI	2256	Bsu15I	547	BsuI	4498	CciI	3102	CciNI	1481
Cfr9I	2066	Clal	547	CpoI	2238	Cspl	2238	DraI	3002	DraIII	2693	DrdI	4594	Dril	1224
DseDI	4594	Eam1105I	1224	Ecl136II	672	Eco31I	2596	Eco53kl	672	Eco57I	4154	Eco91I	2256	EcoCRI	672
EcoNI	3490	EcoO65I	2256	EcoRI	483	FbaI	2096	FblI	1259	FseI	2800	HindIII	833	HpaI	2941
Kpn2I	2235	KpnI	667	KroI	2796	Ksp22I	2096	KspAI	2941	MfeI	2950	MroI	2235	MroNI	2796
MspCI	343	MunI	2950	NaeI	2798	NgoMIV	2796	NotI	1481	NspV	458	PaeR7I	1487	PagI	3102
PasI	2171	Paul	2577	PciI	4696	PdII	2798	Pfi23II	2178	PfiFI	2164	PfoI	2597	PscI	4696
PshBI	3777	Psp124BI	674	Psp1406I	209	PspEI	2256	PspLI	2178	PspXI	1487	PstI	3043	Psyl	2164
Ptel	2577	RcaI	3102	RgaI	3578	RigI	2800	Rsr2I	2238	RsrII	2238	SacI	674	Sall	1258
Scal	554	SfaAI	3578	Sfr274I	1487	SfuI	458	SgfI	3578	Slal	1487	Smal	2068	Sspl	3503
SstI	674	StrI	1487	TspMI	2066	Tth111I	2164	Vha464I	343	VspI	3777	XagI	3490	XbaI	1499
XcmI	1720	XhoI	1487	XmaI	2066	XmaJI	2045	XmiI	1259	Zrml	554				

Table 3. Restriction Enzymes That Cleave TR109 Twice

Acc36I	619	Acc36I	1394	AccB7I	479	AccB7I	3841	Ajul	462	Ajul	3606	Alw44I	1164	Alw44I	4382
AlwNI	1136	AlwNI	4287	Apal	356	Apal	4004	ApaLI	1164	ApaLI	4382	Asp700I	1183	Asp700I	1507
AsuNHI	44	AsuNHI	676	Ball	1015	Ball	2497	BasI	479	BasI	3841	BbsI	134	BbsI	998
Bcgl	71	Bcgl	2563	BfuAI	619	BfuAI	1394	BgIII	475	BgIII	814	BmtI	48	BmtI	680
Bpil	134	Bpil	998	BpII	1980	BpII	2012	Bpu10I	668	Bpu10I	3595	BpuAI	134	BpuAI	998
BpvUI	1256	BpvUI	3578	BsaXI	947	BsaXI	2500	BsmBI	2514	BsmBI	3594	Bsp120I	352	Bsp120I	4000
Bsp1407I	440	Bsp1407I	1621	Bsp19I	920	Bsp19I	1952	Bsp68I	1027	Bsp68I	3235	BspMI	619	BspMI	1394
BspOI	48	BspOI	680	BsrGI	440	BsrGI	1621	BstAUI	440	BstAUI	1621	BstV2I	134	BstV2I	998
BtgZI	939	BtgZI	1296	BtuMI	1027	BtuMI	3235	BveI	619	BveI	1394	Cail	1136	Cail	4287
Cfr42I	1053	Cfr42I	2336	Csil	1812	Csil	2689	DinI	2331	DinI	2462	Drall	353	Drall	4001
Eco147I	2044	Eco147I	2453	EcoO109I	353	EcoO109I	4001	EgeI	2331	EgeI	2462	EheI	2331	EheI	2462
Esp3I	2514	Esp3I	3594	HincII	1260	HincII	2941	HindII	1260	HindII	2941	KasI	2329	KasI	2460
KspI	1053	KspI	2336	MabI	1812	MabI	2689	MisI	1015	MisI	2497	MluNI	1015	MluNI	2497
Mly113I	2330	Mly113I	2461	MroXI	1183	MroXI	1507	MscI	1015	MscI	2497	MslI	524	MslI	958
Msp20I	1015	Msp20I	2497	Mvrl	1256	Mvrl	3578	NarI	2330	NarI	2461	NcoI	920	NcoI	1952
NheI	44	NheI	676	NruI	1027	NruI	3235	PceI	2044	PceI	2453	Pdml	1183	Pdml	1507
PfiMI	479	PfiMI	3841	Ple19I	1256	Ple19I	3578	PluTI	2333	PluTI	2464	PspOMI	352	PspOMI	4000
PstNI	1136	PstNI	4287	PvuI	1256	PvuI	3578	RruI	1027	RruI	3235	RseI	524	RseI	958
SacII	1053	SacII	2336	SexAI	1812	SexAI	2689	SfiI	657	SfiI	827	SfoI	2331	SfoI	2462
Sfr303I	1053	Sfr303I	2336	SgrBI	1053	SgrBI	2336	SmiMI	524	SmiMI	958	SseBI	2044	SseBI	2453
SspDI	2329	SspDI	2460	StuI	2044	StuI	2453	Van91I	479	Van91I	3841	VneI	1164	VneI	4382
XmnI	1183	XmnI	1507												

Table 4. Restriction Enzymes That Cleave TR109 Three Times

AanI	427	AanI	1638	AanI	2921	AatII	66	AatII	2168	AatII	2666	Avall	2238	Avall	2348
Avall	2680	Bfml	3039	Bfml	4240	Bfml	4431	Bme18I	2238	Bme18I	2348	Bme18I	2680	Bmrl	141
Bmrl	611	Bmrl	1924	Bmul	141	Bmul	611	Bmul	1924	Bpml	1102	Bpml	1321	Bpml	2620
Bse3DI	421	Bse3DI	1646	Bse3DI	3076	BseMI	421	BseMI	1646	BseMI	3076	BsrDI	421	BsrDI	1646
BsrDI	3076	Bst6I	329	Bst6I	2281	Bst6I	3391	BstAPI	1790	BstAPI	1862	BstAPI	2693	BstSFI	3039
BstSFI	4240	BstSFI	4431	Eam1104I	329	Eam1104I	2281	Eam1104I	3391	EarI	329	EarI	2281	EarI	3391
Eco47I	2238	Eco47I	2348	Eco47I	2680	Gsul	1102	Gsul	1321	Gsul	2620	Psil	427	Psil	1638
Psil	2921	Sfcl	3039	Sfcl	4240	Sfcl	4431	TaqII	2140	TaqII	2764	TaqII	3832	VpaK11BI	2238
VpaK11BI	2348	VpaK11BI	2680	Zral	64	Zral	2166	Zral	2664						
