

Datasheet of GLuc-ONTM JNK Transcriptional Response Element (TRE)
Clone (TR105)

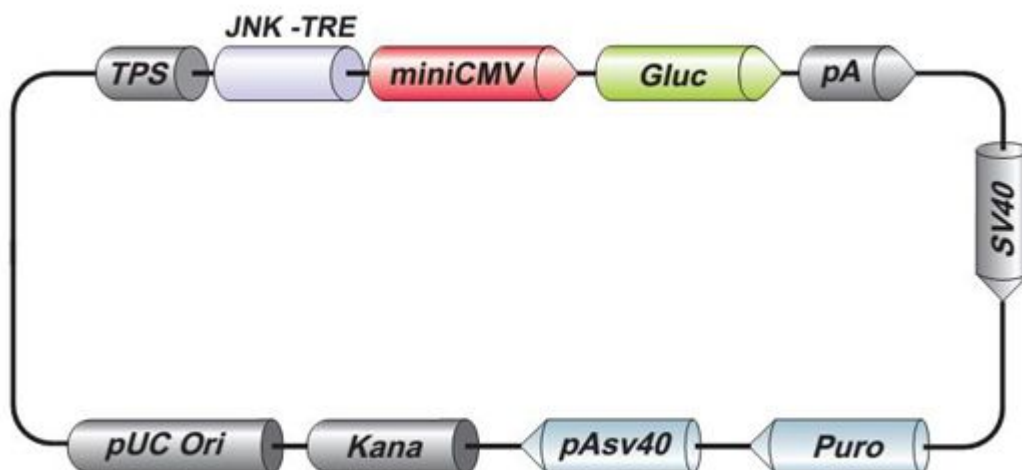


Figure 1. GLuc-ONTM JNK Transcriptional Response Element (TRE) Clone (TR105)

Catalog:

TR105

Description:

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

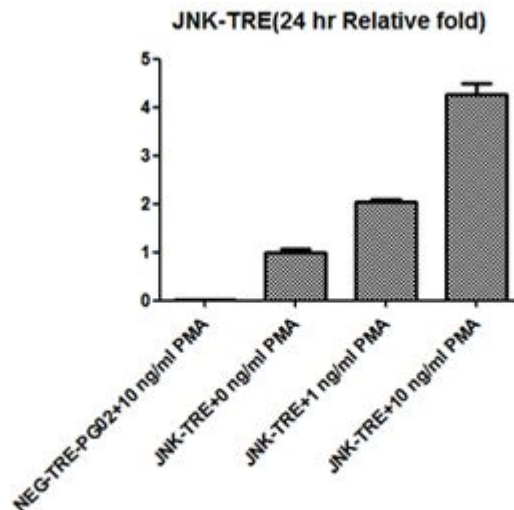


Figure 2. Validation data of GLuc-ON™ JNK Transcriptional Response Element (TRE) Clone(TR105).

Restriction Enzyme Information for TR105

Table 1. Restriction Enzymes That Do Not Cleave TR105

AfeI	AgeI	AscI	BaeI	BaeI	BbvCI	BlpI	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 TR105 Once

AarI	619	AccI	1205	AccIII	2181	AcII	209	AcuI	4101	Adel	2639	AflII	343	AflIII	4643
Alol	895	Aor13HI	2181	Asel	3724	AsiSI	3525	AspA2I	1991	AspI	2110	AssI	554	AsuII	458
AvrII	1991	BamHI	855	Baul	4470	BciVI	4445	BclI	2042	BfrI	343	Bful	4445	BlnI	1991
BmcAI	554	Bpu14I	458	Bsa29I	547	BsaBI	898	Bsal	2542	Bse8I	898	BseAI	2181	BseCI	547
BseJI	898	BsePI	2523	BseRI	1987	BshVI	547	BsiWI	2124	Bso31I	2542	Bsp119I	458	Bsp13I	2181
BspDI	547	BspEI	2181	BspT104I	458	BspTI	343	BssHII	2523	BssSI	4470	Bst2BI	4470	BstAFI	343
BstBI	458	BstEII	2202	BstENI	3437	BstPI	2202	Bsu15I	547	Bsul	4445	CciNI	1427	Cfr9I	2012
Clal	547	CpoI	2184	Cspl	2184	Dral	2949	DraIII	2639	Ecl136II	672	Eco31I	2542	Eco53kl	672
Eco57I	4101	Eco91I	2202	EcoICRI	672	EcoNI	3437	EcoO65I	2202	EcoRI	483	FbaI	2042	FblI	1205
FseI	2747	HindIII	779	HpaI	2888	Kpn2I	2181	KroI	2743	Ksp22I	2042	KspAI	2888	MfeI	2897
MroI	2181	MroNI	2743	MspCI	343	MunI	2897	NaeI	2745	NgoMIV	2743	NotI	1427	NspV	458
PaeR7I	1433	PasI	2117	Paul	2523	PciI	4643	PdII	2745	Pfl23II	2124	PfIFI	2110	PfoI	2543
PscI	4643	PshBI	3724	Psp124BI	674	Psp1406I	209	PspEI	2202	PspLI	2124	PspXI	1433	PstI	2990
PsyI	2110	PteI	2523	PvuII	736	RgaI	3525	RigI	2747	Rsr2I	2184	RsrII	2184	SacI	674
Sall	1204	Scal	554	SfaAI	3525	Sfr274I	1433	SfuI	458	Sgfl	3525	Slal	1433	Smal	2014
Sspl	3450	SstI	674	StrI	1433	TspMI	2012	Tth111I	2110	Vha464I	343	VspI	3724	XagI	3437
XbaI	1445	XcmI	1666	XhoI	1433	XmaI	2012	XmaJI	1991	XmiI	1205	Zrml	554		

Table 3. Restriction Enzymes That Cleave TR105 Twice

AasI	729	AasI	4541	Acc36I	619	Acc36I	1340	Acc65I	663	Acc65I	2724	AccB7I	479	AccB7I	3788
AhdI	744	AhdI	1170	AjuI	462	AjuI	3553	Alw44I	1110	Alw44I	4329	AlwNI	1082	AlwNI	4234
Apal	356	Apal	3951	ApalI	1110	ApalI	4329	Asp700I	1129	Asp700I	1453	Asp718I	663	Asp718I	2724
AspEI	744	AspEI	1170	AsuNHI	44	AsuNHI	676	Ball	961	Ball	2443	BasI	479	BasI	3788
BbsI	134	BbsI	944	Bcgl	71	Bcgl	2509	BfuAI	619	BfuAI	1340	BglII	475	BglII	760
BmeRI	744	BmeRI	1170	BmtI	48	BmtI	680	Bpil	134	Bpil	944	BpII	1926	BpII	1958
Bpu10I	668	Bpu10I	3542	BpuAI	134	BpuAI	944	BpvUI	1202	BpvUI	3525	BsaXI	893	BsaXI	2446
BsmBI	2460	BsmBI	3541	Bsp120I	352	Bsp120I	3947	Bsp1407I	440	Bsp1407I	1567	Bsp19I	866	Bsp19I	1898
Bsp68I	973	Bsp68I	3182	BspHI	726	BspHI	3049	BspMI	619	BspMI	1340	BspOI	48	BspOI	680
BsrGI	440	BsrGI	1567	BstAUI	440	BstAUI	1567	BstV2I	134	BstV2I	944	BtgZI	885	BtgZI	1242
BtuMI	973	BtuMI	3182	Bvel	619	Bvel	1340	Cail	1082	Cail	4234	Ccil	726	Ccil	3049
Cfr42I	999	Cfr42I	2282	Csil	1758	Csil	2635	DinI	2277	DinI	2408	Drall	353	Drall	3948
DrdI	729	DrdI	4541	Dril	744	Dril	1170	DseDI	729	DseDI	4541	Eam1105I	744	Eam1105I	1170
Eco147I	1990	Eco147I	2399	EcoO109I	353	EcoO109I	3948	Egel	2277	Egel	2408	Ehel	2277	Ehel	2408
Esp3I	2460	Esp3I	3541	HincII	1206	HincII	2888	HindII	1206	HindII	2888	KasI	2275	KasI	2406
KpnI	667	KpnI	2728	KspI	999	KspI	2282	MabI	1758	MabI	2635	MisI	961	MisI	2443
MluNI	961	MluNI	2443	Mly113I	2276	Mly113I	2407	MroXI	1129	MroXI	1453	MscI	961	MscI	2443
MslI	524	MslI	904	Msp20I	961	Msp20I	2443	MvrI	1202	MvrI	3525	NarI	2276	NarI	2407
NcoI	866	NcoI	1898	NheI	44	NheI	676	NruI	973	NruI	3182	PagI	726	PagI	3049
PceI	1990	PceI	2399	Pdml	1129	Pdml	1453	PfIM	479	PfIM	3788	Ple19I	1202	Ple19I	3525
PluTI	2279	PluTI	2410	PspOMI	352	PspOMI	3947	PstNI	1082	PstNI	4234	PvuI	1202	PvuI	3525
RcaI	726	RcaI	3049	RruI	973	RruI	3182	RseI	524	RseI	904	SacI	999	SacI	2282
SexAI	1758	SexAI	2635	SfiI	657	SfiI	773	SfoI	2277	SfoI	2408	Sfr303I	999	Sfr303I	2282
SgrBI	999	SgrBI	2282	SmiMI	524	SmiMI	904	SseBI	1990	SseBI	2399	SspDI	2275	SspDI	2406
StuI	1990	StuI	2399	Van91I	479	Van91I	3788	VneI	1110	VneI	4329	XmnI	1129	XmnI	1453

Table 4. Restriction Enzymes That Cleave TR105 Three Times

AanI	427	AanI	1584	AanI	2868	AatII	66	AatII	2114	AatII	2612	Avall	2184	Avall	2294
Avall	2626	BfmI	2986	BfmI	4187	BfmI	4378	Bme18I	2184	Bme18I	2294	Bme18I	2626	Bmrl	141
Bmrl	611	Bmrl	1870	Bmul	141	Bmul	611	Bmul	1870	Bpml	1048	Bpml	1267	Bpml	2566
Bse3DI	421	Bse3DI	1592	Bse3DI	3023	BseMI	421	BseMI	1592	BseMI	3023	BsrDI	421	BsrDI	1592
BsrDI	3023	Bst6I	329	Bst6I	2227	Bst6I	3338	BstAPI	1736	BstAPI	1808	BstAPI	2639	BstSFI	2986
BstSFI	4187	BstSFI	4378	Eam1104I	329	Eam1104I	2227	Eam1104I	3338	EarI	329	EarI	2227	EarI	3338
Eco47I	2184	Eco47I	2294	Eco47I	2626	Gsul	1048	Gsul	1267	Gsul	2566	Psil	427	Psil	1584
Psil	2868	Sfcl	2986	Sfcl	4187	Sfcl	4378	TaqII	2086	TaqII	2710	TaqII	3779	VpaK11BI	2184
VpaK11BI	2294	VpaK11BI	2626	Zral	64	Zral	2112	Zral	2610						
