

Datasheet of GLuc-ONTM HNF4 Transcriptional Response Element (TRE)
Clone (TR104)

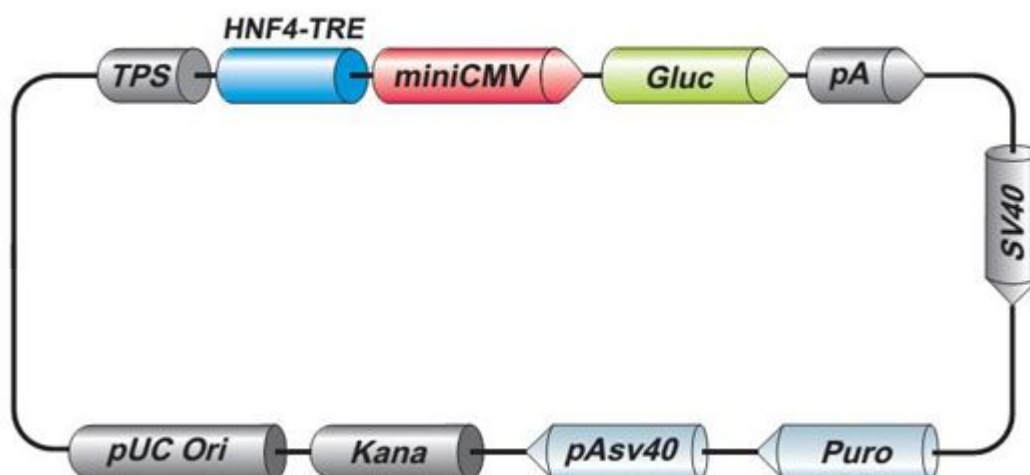


Figure 1. GLuc-ONTM HNF4 Transcriptional Response Element (TRE) Clone (TR104)

Catalog:

TR104

Description:

The HNF4 transcriptional response element (TRE) clone assays the status of the HNF4 signaling pathway through the activity of the transcription factor HNF4. The HNF4 transcriptional TRE clone contains a minimal promoter and tandem repeats of the HNF4 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

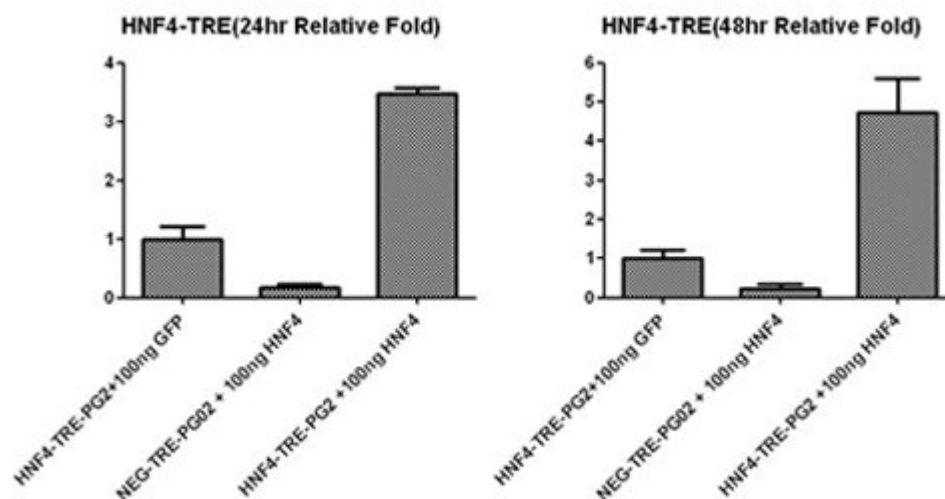


Figure 2. Validation data of GLuc-ON™ HNF4 Transcriptional Response Element (TRE) Clone(TR104).

Restriction Enzyme Information for TR104

Table 1. Restriction Enzymes That Do Not Cleave TR104

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 TR104 Once

AarI	619	AasI	4535	AccI	1199	AccIII	2175	AclI	209	AcuI	4095	Adel	2633	AflII	343
AflIII	4637	AhdI	1164	Alol	889	Aor13HI	2175	Asel	3718	AsiSI	3519	AspA2I	1985	AspEI	1164
AspI	2104	AssI	554	AsuII	458	AvrII	1985	BamHI	849	BauI	4464	BciVI	4439	BclI	2036
BfrI	343	BfuI	4439	BlnI	1985	BmcAI	554	BmeRI	1164	Bpu14I	458	Bsa29I	547	BsaBI	892
BsaI	2536	Bse8I	892	BseAI	2175	BseCI	547	BseJI	892	BsePI	2517	BseRI	1981	BshVI	547
BsiWI	2118	Bso31I	2536	Bsp119I	458	Bsp13I	2175	BspDI	547	BspEI	2175	BspHI	3043	BspT104I	458
BspTI	343	BssHII	2517	BssSI	4464	Bst2BI	4464	BstAFI	343	BstBI	458	BstEII	2196	BstENI	3431
BstPI	2196	Bsu15I	547	Bsul	4439	Ccil	3043	CciNI	1421	Cfr9I	2006	Clal	547	CpoI	2178
Cspl	2178	Dral	2943	DrallI	2633	DrdI	4535	Dril	1164	DseDI	4535	Eam1105I	1164	Ecl136II	672
Eco31I	2536	Eco53kI	672	Eco57I	4095	Eco91I	2196	EcoICRI	672	EcoNI	3431	EcoO65I	2196	EcoRI	483
FbaI	2036	FblI	1199	FseI	2741	HindIII	773	HpaI	2882	Kpn2I	2175	KroI	2737	Ksp22I	2036
KspAI	2882	MfeI	2891	MroI	2175	MroNI	2737	MspCI	343	MunI	2891	NaeI	2739	NgoMIV	2737
NotI	1421	NspV	458	PaeR7I	1427	PagI	3043	PasI	2111	Paul	2517	Pcil	4637	PdII	2739
Pfi23II	2118	PfiFI	2104	PfoI	2537	PscI	4637	PshBI	3718	Psp124BI	674	Psp1406I	209	PspEI	2196
PspLI	2118	PspXI	1427	PstI	2984	PsyI	2104	Ptel	2517	Rcal	3043	Rgal	3519	RigI	2741
Rsr2I	2178	RsrII	2178	SacI	674	Sall	1198	Scal	554	SfaAI	3519	Sfr274I	1427	SfuI	458
SgfI	3519	Slal	1427	SmaI	2008	SspI	3444	SstI	674	StrI	1427	TspMI	2006	Tth111I	2104
Vha464I	343	Vspl	3718	XagI	3431	XbaI	1439	XcmI	1660	XhoI	1427	XmaI	2006	XmaJI	1985
XmiI	1199	Zrml	554												

Table 3. Restriction Enzymes That Cleave TR104 Twice

Acc36I	619	Acc36I	1334	Acc65I	663	Acc65I	2718	AccB7I	479	AccB7I	3782	Ajul	462	Ajul	3547
Alw44I	1104	Alw44I	4323	AlwNI	1076	AlwNI	4228	Apal	356	Apal	3945	ApaLI	1104	ApaLI	4323
Asp700I	1123	Asp700I	1447	Asp718I	663	Asp718I	2718	AsuNHI	44	AsuNHI	676	Ball	955	Ball	2437
BasI	479	BasI	3782	BbsI	134	BbsI	938	Bcgl	71	Bcgl	2503	BfuAI	619	BfuAI	1334
BglII	475	BglII	754	BmtI	48	BmtI	680	Bpil	134	Bpil	938	BpII	1920	BpII	1952
Bpu10I	668	Bpu10I	3536	BpuAI	134	BpuAI	938	BpvUI	1196	BpvUI	3519	BsaXI	887	BsaXI	2440
BsmBI	2454	BsmBI	3535	Bsp120I	352	Bsp120I	3941	Bsp1407I	440	Bsp1407I	1561	Bsp19I	860	Bsp19I	1892
Bsp68I	967	Bsp68I	3176	BspMI	619	BspMI	1334	BspOI	48	BspOI	680	BsrGI	440	BsrGI	1561
BstAUI	440	BstAUI	1561	BstV2I	134	BstV2I	938	BtgZI	879	BtgZI	1236	BtuMI	967	BtuMI	3176
BveI	619	BveI	1334	Cail	1076	Cail	4228	Cfr42I	993	Cfr42I	2276	Csil	1752	Csil	2629
DinI	2271	DinI	2402	Drall	353	Drall	3942	Eco147I	1984	Eco147I	2393	EcoO109I	353	EcoO109I	3942
EgeI	2271	EgeI	2402	EheI	2271	EheI	2402	Esp3I	2454	Esp3I	3535	HincII	1200	HincII	2882
HindII	1200	HindII	2882	KasI	2269	KasI	2400	KpnI	667	KpnI	2722	KspI	993	KspI	2276
MabI	1752	MabI	2629	MlsI	955	MlsI	2437	MluNI	955	MluNI	2437	Mly113I	2270	Mly113I	2401
MroXI	1123	MroXI	1447	MscI	955	MscI	2437	MslI	524	MslI	898	Msp20I	955	Msp20I	2437
MvrI	1196	MvrI	3519	NarI	2270	NarI	2401	NcoI	860	NcoI	1892	NheI	44	NheI	676
NruI	967	NruI	3176	PceI	1984	PceI	2393	Pdml	1123	Pdml	1447	PfiMI	479	PfiMI	3782
Ple19I	1196	Ple19I	3519	PluTI	2273	PluTI	2404	PspOMI	352	PspOMI	3941	PstNI	1076	PstNI	4228
PvuI	1196	PvuI	3519	RruI	967	RruI	3176	RseI	524	RseI	898	SacII	993	SacII	2276
SexAI	1752	SexAI	2629	SfiI	657	SfiI	767	SfoI	2271	SfoI	2402	Sfr303I	993	Sfr303I	2276
SgrBI	993	SgrBI	2276	SmiMI	524	SmiMI	898	SseBI	1984	SseBI	2393	SspDI	2269	SspDI	2400
StuI	1984	StuI	2393	Van91I	479	Van91I	3782	VneI	1104	VneI	4323	XmnI	1123	XmnI	1447

Table 4. Restriction Enzymes That Cleave TR104 Three Times

AanI	427	AanI	1578	AanI	2862	AatII	66	AatII	2108	AatII	2606	Avall	2178	Avall	2288
Avall	2620	BfmI	2980	BfmI	4181	BfmI	4372	Bme18I	2178	Bme18I	2288	Bme18I	2620	Bmrl	141
Bmrl	611	Bmrl	1864	Bmul	141	Bmul	611	Bmul	1864	Bpml	1042	Bpml	1261	Bpml	2560
Bse3DI	421	Bse3DI	1586	Bse3DI	3017	BseMI	421	BseMI	1586	BseMI	3017	BsrDI	421	BsrDI	1586
BsrDI	3017	Bst6I	329	Bst6I	2221	Bst6I	3332	BstAPI	1730	BstAPI	1802	BstAPI	2633	BstSFI	2980
BstSFI	4181	BstSFI	4372	Eam1104I	329	Eam1104I	2221	Eam1104I	3332	EarI	329	EarI	2221	EarI	3332
Eco47I	2178	Eco47I	2288	Eco47I	2620	Gsul	1042	Gsul	1261	Gsul	2560	Psil	427	Psil	1578
Psil	2862	Sfcl	2980	Sfcl	4181	Sfcl	4372	TaqII	2080	TaqII	2704	TaqII	3773	VpaK11BI	2178
VpaK11BI	2288	VpaK11BI	2620	Zral	64	Zral	2106	Zral	2604						
