

Datasheet of GLuc-ONTM Negative Control Clone (TR001) for Transcriptional Response Assay

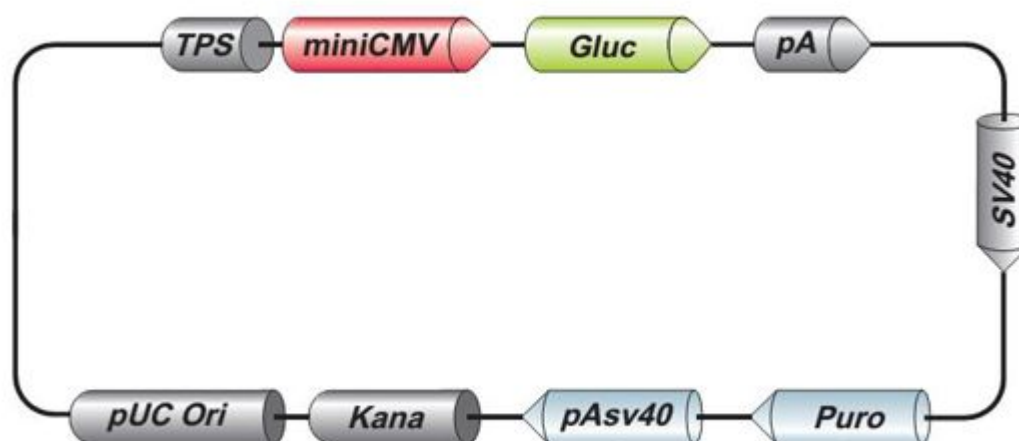


Figure 1. GLuc-ONTM Negative Control Clone (TR001) for Transcriptional Response Assay

Catalog:	TR001
Description:	Negative Control Clone TR001 contains a minimal promoter upstream of a secreted Gaussia luciferase reporter gene.
Delivery:	500ng/ μ L $\times$ 40 μ L
Storage Conditions:	-20°C
QC:	. Full sequence verification . Restriction Digestion

Restriction Enzyme Information for TR001

Table 1. Restriction Enzymes That Do Not Cleave TR001

AfeI	AgeI	AscI	BaeI	BaeI	BbvCI	BlnI	BsaAI
BstXI	BstZ17I	Bsu36I	BtrI	Ecl136II	EcoRV	FspI	NdeI
PacI	PmeI	PmlI	PpuMI	PshAI	PvuII	SacI	SanDI
SapI	SbfI	SfiI	SgrAI	SnaBI	SrfI		

Table 2. Restriction Enzymes That Cleave TR001 Once

AarI	619	AasI	4426	Acc65I	2609	AccI	1090	AccIII	2066	AcII	209	AcuI	3986	Adel	2524
AflII	343	AhdI	1055	AhII	650	Alol	780	Aor13HI	2066	Asel	3609	AsiSI	3410	Asp718I	2609
AspA2I	1876	AspEI	1055	AspI	1995	AssI	554	AsuII	458	AsuNHI	44	AvrII	1876	BamHI	740
BauI	4355	BciVI	4330	BclI	1927	BcuI	650	BfrI	343	Bful	4330	BgIII	475	BlnI	1876
BmcAI	554	BmeRI	1055	BmtI	48	Bpu10I	3427	Bpu14I	458	Bsa29I	547	BsaBI	783	BsaI	2427
Bse8I	783	BseAI	2066	BseCI	547	BseJI	783	BsePI	2408	BseRI	1872	BshVI	547	BsiWI	2009
Bso31I	2427	Bsp119I	458	Bsp13I	2066	BspDI	547	BspEI	2066	BspHI	2934	BspOI	48	BspT104I	458
BspTI	343	BssHII	2408	BssSI	4355	Bst2BI	4355	BstAFI	343	BstBI	458	BstEII	2087	BstENI	3322
BstPI	2087	Bsu15I	547	Bsul	4330	CciI	2934	CciNI	1312	Cfr9I	1897	Clal	547	CpoI	2069
CspI	2069	DraIII	2524	DrdI	4426	Dril	1055	DseDI	4426	Eam1105I	1055	Eco31I	2427	Eco57I	3986
Eco91I	2087	EcoNI	3322	EcoO65I	2087	EcoRI	483	FbaI	1927	FblI	1090	FseI	2632	HindIII	670
HpaI	2773	Kpn2I	2066	KpnI	2613	KroI	2628	Ksp22I	1927	KspAI	2773	MfeI	2782	MluI	656
MroI	2066	MroNI	2628	MspCI	343	MunI	2782	NaeI	2630	NgoMIV	2628	NheI	44	NotI	1312
NspV	458	PaeR7I	1318	PagI	2934	PasI	2002	Paul	2408	PciI	4528	PdII	2630	Pfi23II	2009
PfiFI	1995	PfoI	2428	PscI	4528	PshBI	3609	Psp1406I	209	PspEI	2087	PspLI	2009	PspXI	1318
PstI	2875	PsyI	1995	PteI	2408	RcaI	2934	RgaI	3410	RigI	2632	Rsr2I	2069	RsrII	2069
Sall	1089	Scal	554	SfaAI	3410	Sfr274I	1318	Sfal	458	Sgfl	3410	Slal	1318	SmaI	1899
Smil	665	SpeI	650	SspI	3335	StrI	1318	Swal	665	TspMI	1897	Tth111I	1995	Vha464I	343
VspI	3609	XagI	3322	XbaI	1330	XcmI	1551	XhoI	1318	XmaI	1897	XmaJI	1876	XmiI	1090
Zrml	554														

Table 3. Restriction Enzymes That Cleave TR001 Twice

Acc36I	619	Acc36I	1225	AccB7I	479	AccB7I	3673	AfilII	656	AfilII	4528	Ajul	462	Ajul	3438
Alw44I	995	Alw44I	4214	AlwNI	967	AlwNI	4119	Apal	356	Apal	3836	ApaLI	995	ApaLI	4214
Asp700I	1014	Asp700I	1338	Ball	846	Ball	2328	BasI	479	BasI	3673	BbsI	134	BbsI	829
Bcgl	71	Bcgl	2394	BfuAI	619	BfuAI	1225	BgII	2169	BgII	2300	Bpil	134	Bpil	829
BpII	1811	BpII	1843	BpuAI	134	BpuAI	829	BpvUI	1087	BpvUI	3410	BsaXI	778	BsaXI	2331
BsmBI	2345	BsmBI	3426	Bsp120I	352	Bsp120I	3832	Bsp1407I	440	Bsp1407I	1452	Bsp19I	751	Bsp19I	1783
Bsp68I	858	Bsp68I	3067	BspMI	619	BspMI	1225	BsrGI	440	BsrGI	1452	BstAUl	440	BstAUl	1452
BstV2I	134	BstV2I	829	BtgZI	770	BtgZI	1127	BtuMI	858	BtuMI	3067	BveI	619	BveI	1225
Cail	967	Cail	4119	Cfr42I	884	Cfr42I	2167	Csil	1643	Csil	2520	DinI	2162	DinI	2293
Dral	665	Dral	2834	Drall	353	Drall	3833	Eco147I	1875	Eco147I	2284	EcoO109I	353	EcoO109I	3833
Egel	2162	Egel	2293	Ehel	2162	Ehel	2293	Esp3I	2345	Esp3I	3426	HincII	1091	HincII	2773
HindII	1091	HindII	2773	KasI	2160	KasI	2291	Kspl	884	Kspl	2167	MabI	1643	MabI	2520
MisI	846	MisI	2328	MluNI	846	MluNI	2328	Mly113I	2161	Mly113I	2292	MroXI	1014	MroXI	1338
MscI	846	MscI	2328	MslI	524	MslI	789	Msp20I	846	Msp20I	2328	MvrI	1087	MvrI	3410
NarI	2161	NarI	2292	NcoI	751	NcoI	1783	Nrul	858	Nrul	3067	PceI	1875	PceI	2284
Pdml	1014	Pdml	1338	PfiMI	479	PfiMI	3673	Ple19I	1087	Ple19I	3410	PluTI	2164	PluTI	2295
PspOMI	352	PspOMI	3832	PstNI	967	PstNI	4119	Pvul	1087	Pvul	3410	Rrul	858	Rrul	3067
Rsel	524	Rsel	789	SacII	884	SacII	2167	SexAI	1643	SexAI	2520	Sfol	2162	Sfol	2293
Sfr303I	884	Sfr303I	2167	SgrBI	884	SgrBI	2167	SmiMI	524	SmiMI	789	SseBI	1875	SseBI	2284
SspDI	2160	SspDI	2291	Stul	1875	Stul	2284	Van91I	479	Van91I	3673	VneI	995	VneI	4214
Xmnl	1014	Xmnl	1338												

Table 4. Restriction Enzymes That Cleave TR001 Three Times

AanI	427	AanI	1469	AanI	2753	AatII	66	AatII	1999	AatII	2497	Avall	2069	Avall	2179
Avall	2511	Bfml	2871	Bfml	4072	Bfml	4263	Bme18I	2069	Bme18I	2179	Bme18I	2511	Bmrl	141

Bmrl	611	Bmrl	1755	Bmul	141	Bmul	611	Bmul	1755	Bpml	933	Bpml	1152	Bpml	2451
Bse3DI	421	Bse3DI	1477	Bse3DI	2908	BseMI	421	BseMI	1477	BseMI	2908	BsrDI	421	BsrDI	1477
BsrDI	2908	Bst6I	329	Bst6I	2112	Bst6I	3223	BstAPI	1621	BstAPI	1693	BstAPI	2524	BstSFI	2871
BstSFI	4072	BstSFI	4263	Eam1104I	329	Eam1104I	2112	Eam1104I	3223	EarI	329	EarI	2112	EarI	3223
Eco47I	2069	Eco47I	2179	Eco47I	2511	Gsul	933	Gsul	1152	Gsul	2451	Psil	427	Psil	1469
Psil	2753	Sfcl	2871	Sfcl	4072	Sfcl	4263	TaqII	1971	TaqII	2595	TaqII	3664	VpaK11BI	2069
VpaK11BI	2179	VpaK11BI	2511	Zral	64	Zral	1997	Zral	2495						
