

Datasheet of GLuc-ON™ PR Transcriptional Response Element (TRE) Clone (TR113)

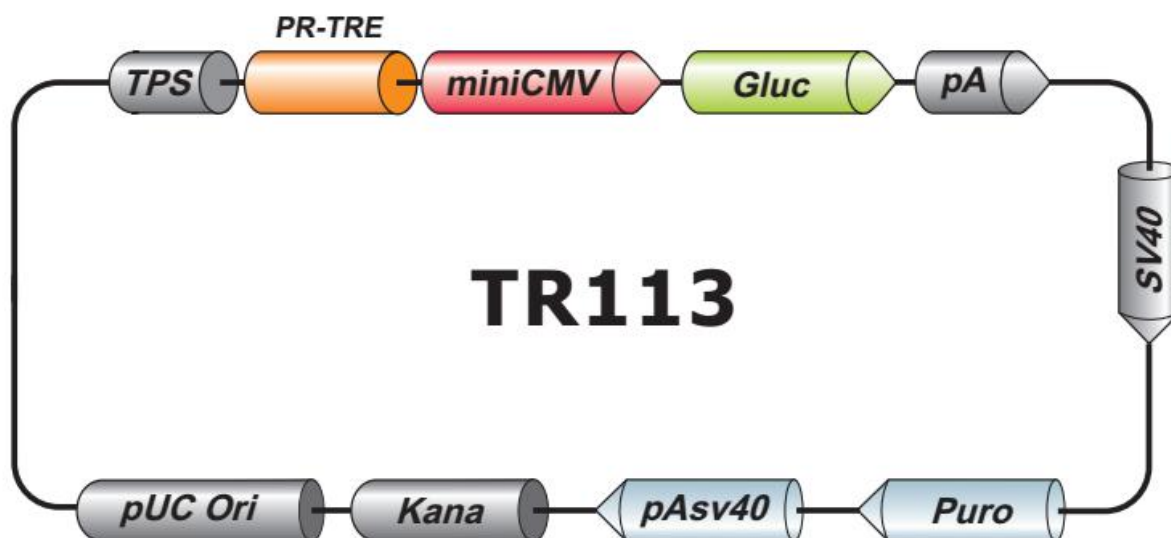


Figure 1. GLuc-ON™ PR Transcriptional Response Element (TRE) Clone (TR113)

Cat. No	TR113
Description	GeneCopoeia 's GLuc-ON™ PR transcriptional response element (TRE) clone enables analysis of the activity of signal transduction mediated by progesterone receptors (PR). Progesterone receptors are transcription factors that, in response to stimulation by progesterone, lead to proliferation of breast tissue. Aberrant function and/or expression of PRs is an important factor in development of breast cancer. The PR transcriptional TRE clone contains a minimal promoter and tandem repeats of the PR 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 × 40μL
Storage Conditions	-20°C
QC	• Functional validation in common cell line • Full sequence verification • Restriction Digestion

Validation Data

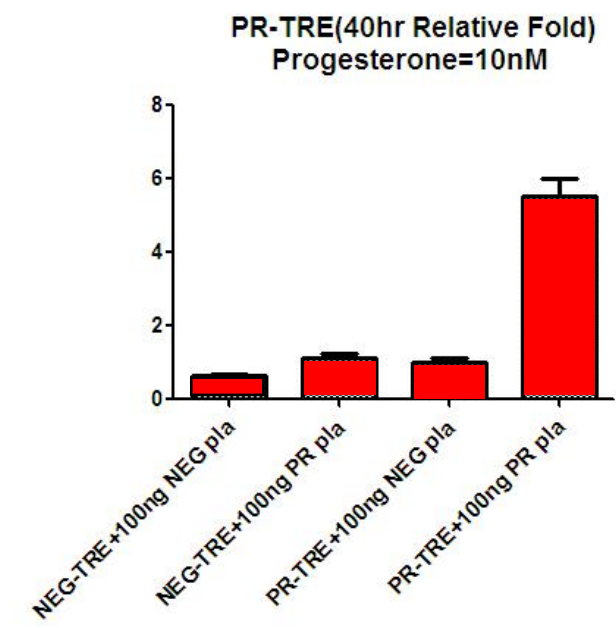


Figure 2. Activation of the GLuc-ON™ PR TRE by exogenous expression of PR. 293T cells were co-transfected with PR TRE clone + PR expression clone, PR TRE clone + negative expression clone, negative control clone + PR expression clone, or negative TRE clone + negative expression clone. Transfected cells were incubated for 40 hours. A Gaussia Luciferase assay was performed, and transcriptional response activity values are expressed as luminance fold activation.

Restriction Enzyme Information for TR113

Table 1. Restriction Enzymes That Do Not Cleave TR113

AfeI	AgeI	AscI	BaeI	BaeI	BbvCI	BlpI	BsaAI
BstXI	BstZ17I	Bsu36I	BtrI	EcoRV	FspI	MluI	NdeI
PacI	PmeI	PmlI	PpuMI	PshAI	PvuII	SanDI	SapI
SbfI	SgrAI	S	naBI	SpeI	SrfI	Swal	

Table 2. Restriction Enzymes That Cleave TR113 Once

AarI	619 AasI	4553 AccI	1217 AccIII	2193 AclI	209 AcuI	4113 Adel	2651 AflII	343
AflIII	4655 AhdI	1182 Aor13HI	2193 AseI	3736 AsiSI	3537 AspA2I	2003 AspEI	1182 Aspl	2122
AssI	554 AsuII	458 AvrII	2003 BamHI	867 Baul	4482 BciVI	4457 BclI	2054 BfrI	343
BfuI	4457 BlnI	2003 BmcAI	554 BmeRI	1182 Bpu14I	458 Bsa29I	547 BsaBI	910 Bsal	2554
Bse8I	910 BseAI	2193 BseCI	547 BseJI	910 BsePI	2535 BseRI	1999 BshVI	547 BsiWI	2136
Bso31I	2554 Bsp119I	458 Bsp13I	2193 BspDI	547 BspEI	2193 BspHI	3061 BspT104I	458 BspTI	343
BssHII	2535 BssSI	4482 Bst2BI	4482 BstAFI	343 BstBI	458 BstEII	2214 BstENI	3449 BstPI	2214
Bsu15I	547 Bsul	4457 CciI	3061 CciNI	1439 Cfr9I	2024 Clal	547 Cpol	2196 Cspl	2196
DraI	2961 DraIII	2651 DrdI	4553 Dril	1182 DseDI	4553 Eam1105I	1182 Ecl136II	672 Eco31I	2554
Eco53kl	672 Eco57I	4113 Eco91I	2214 EcoICRI	672 EcoNI	3449 EcoO65I	2214 EcoRI	483 Fbal	2054
FblI	1217 FseI	2759 HindIII	791 HpaI	2900 Kpn2I	2193 Krol	2755 Ksp22I	2054 KspAI	2900
MfeI	2909 MroI	2193 MroNI	2755 MspCI	343 MunI	2909 NaeI	2757 NgoMIV	2755 NotI	1439
NspV	458 Paer7I	1445 PagI	3061 PasI	2129 Paul	2535 PciI	4655 PdiI	2757 Pfl23II	2136
PfIFI	2122 PfoI	2555 Psci	4655 PshBI	3736 Psp124BI	674 Psp1406I	209 PspEI	2214 PspLI	2136
PspXI	1445 PstI	3002 Psyl	2122 PteI	2535 Rcal	3061 Rgal	3537 Rigl	2759 Rsr2I	2196
RsrII	2196 SacI	674 Sall	1216 Scal	554 SfaAI	3537 Sfr274I	1445 Sful	458 Sgfl	3537
Slal	1445 Smal	2026 Sspl	3462 SstI	674 StrI	1445 TspMI	2024 Tth111I	2122 Vha464I	343
Vspl	3736 XagI	3449 XcmI	1678 XhoI	1445 XmaI	2024 XmaJI	2003 Xmil	1217 Zrml	554

Table 3. Restriction Enzymes That Cleave TR113 Twice

Acc36I	619 Acc36I	1352 Acc65I	663 Acc65I	2736 AccB7I	479 AccB7I	3800 Ajul	462 Ajul3	565
Alw44I	1122 Alw44I	4341 AlwNI	1094 AlwNI	4246 Apal	356 Apal	3963 ApaLI	1122 ApaLI	4341
Asp700I	1141 Asp700I	1465 Asp718I	663 Asp718I	2736 AsuNHI	44 AsuNHI	676 Ball	973 Ball	2455
BasI	479 BasI	3800 BbsI	134 BbsI	956 Bcgl	71 Bcgl	2521 BfuAI	619 BfuAI	1352
BglII	475 BglII	772 BmtI	48 BmtI	680 Bpil	134 Bpil	956 BpII	1938 BpII	1970
Bpu10I	668 Bpu10I	3554 BpuAI	134 BpuAI	956 BpvUI	1214 BpvUI	3537 BsaXI	905 BsaXI	2458
BsmBI	2472 BsmBI	3553 Bsp120I	352 Bsp120I	3959 Bsp1407I	440 Bsp1407I	1579 Bsp19I	878 Bsp19I	1910
Bsp68I	985 Bsp68I	3194 BspMI	619 BspMI	1352 BspOI	48 BspOI	680 BsrGI	440 BsrGI	1579
BstAUI	440 BstAUI	1579 BstV2I	134 BstV2I	956 BtgZI	897 BtgZI	1254 BtuMI	985 BtuMI	3194
BveI	619 BveI	1352 Cail	1094 Cail	4246 Cfr42I	1011 Cfr42I	2294 Csil	1770 Csil	2647
DinI	2289 DinI	2420 Drall	353 Drall	3960 Eco147I	2002 Eco147I	2411 EcoO109I	353 EcoO109I	3960
EgeI	2289 EgeI	2420 Ehel	2289 Ehel	2420 Esp3I	2472 Esp3I	3553 HincII	1218 HincII	2900
HindII	1218 HindII	2900 KasI	2287 KasI	2418 KpnI	667 KpnI	2740 KspI	1011 KspI	2294
MabI	1770 MabI	2647 MlsI	973 MlsI	2455 MluNI	973 MluNI	2455 Mly113I	2288 Mly113I	2419
MroXI	1141 MroXI	1465 MscI	973 MscI	2455 MslI	524 MslI	916 Msp20I	973 Msp20I	2455
MvrI	1214 MvrI	3537 NarI	2288 NarI	2419 NcoI	878 NcoI	1910 NheI	44 NheI	676
NruI	985 NruI	3194 PceI	2002 PceI	2411 PdmI	1141 PdmI	1465 PflMI	479 PflMI	3800
Ple19I	1214 Ple19I	3537 PluTI	2291 PluTI	2422 PspOMI	352 PspOMI	3959 PstNI	1094 PstNI	4246
PvuI	1214 PvuI	3537 RruI	985 RruI	3194 RseI	524 RseI	916 SacII	1011 SacII	2294
SexAI	1770 SexAI	2647 SfiI	657 SfiI	785 SfoI	2289 SfoI	2420 Sfr303I	1011 Sfr303I	2294
SgrBI	1011 SgrBI	2294 SmiMI	524 SmiMI	916 SseBI	2002 SseBI	2411 SspDI	2287 SspDI	2418
StuI	2002 StuI	2411 Van91I	479 Van91I	3800 VneI	1122 VneI	4341 XbaI	769 XbaI	1457
XmnI	1141 XmnI	1465						

Table 4. Restriction Enzymes That Cleave TR113 Three times

AanI	427 AanI	1596 AanI	2880 AatII	66 AatII	2126 AatII	2624 Avall	2196 Avall	2306
Avall	2638 Bfml	2998 Bfml	4199 Bfml	4390 Bme18I	2196 Bme18I	2306 Bme18I	2638 Bmrl	141
Bmrl	611 Bmrl	1882 Bmul	141 Bmul	611 Bmul	1882 Bpml	1060 Bpml	1279 Bpml	2578
Bse3DI	421 Bse3DI	1604 Bse3DI	3035 BseMI	421 BseMI	1604 BseMI	3035 BsrDI	421 BsrDI	1604
BsrDI	3035 Bst6I	329 Bst6I	2239 Bst6I	3350 BstAPI	1748 BstAPI	1820 BstAPI	2651 BstSFI	2998
BstSFI	4199 BstSFI	4390 Eam1104I	329 Eam1104I	2239 Eam1104I	3350 Earl	329 Earl	2239 Earl	3350
Eco47I	2196 Eco47I	2306 Eco47I	2638 Gsul	1060 Gsul	1279 Gsul	2578 Psil	427 Psil	1596
Psil	2880 Sfcl	2998 Sfcl	4199 Sfcl	4390 TaqII	2098 TaqII	2722 TaqII	3791 VpaK11BI	2196
VpaK11BI	2306 VpaK11BI	2638 Zral	64 Zral	2124 Zral	2622			