

Datasheet of GLuc-ON™ NFAT Transcriptional Response Element (TRE) Lentiviral Clone (TR201)

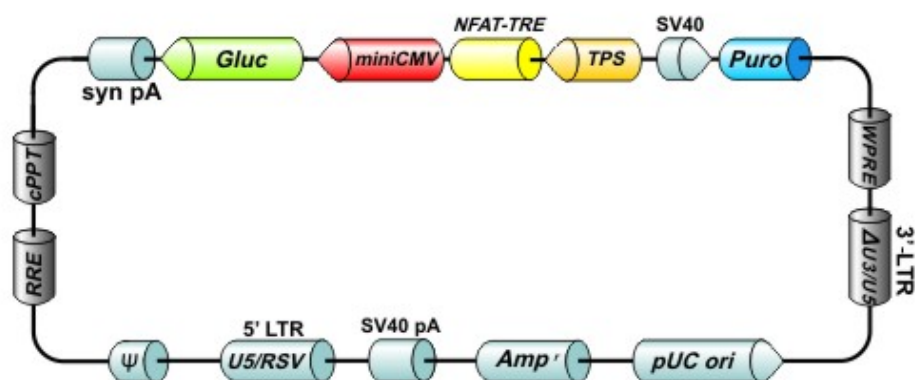


Figure 1. GLuc-ON™ NFAT Transcriptional Response Element (TRE) Lentiviral Clone (TR201)

Catalog: TR201

Description:

The NFAT transcriptional response element (TRE) lentiviral clone assays the status of the PKC/Ca⁺⁺ signaling pathway through the activity of the transcription factor NFAT. The NFAT transcriptional TRE clone contains a minimal promoter and tandem repeats of the NFAT 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: 500 ng/μL × 40 μL

Storage Conditions: -20°C

.Functional validation in common cell line

QC: .Full sequence verification

.Restriction Digestion

Validation Data

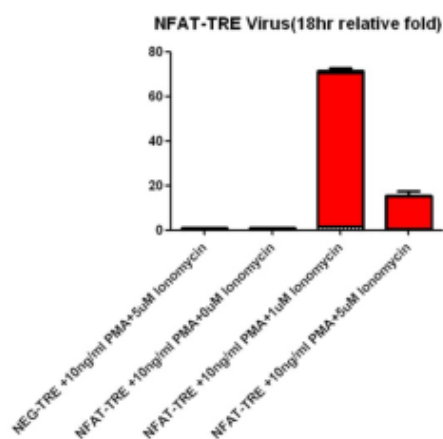


Figure 2. Validation data of GLuc-ON™ NFAT Transcriptional Response Element (TRE) Lentiviral Clone (TR201).

Restriction Enzyme Information for TR201

Table 1. Restriction Enzymes That Do Not Cleave TR201

AfeI	AgeI	AscI	BclI	BglI	BspEI	BsrGI	BstXI
BstZ17I	HpaI	MfeI	NdeI	NotI	PacI	PmeI	PshAI
SbfI	SgfI	SgrAI	SnaBI	SpeI	SrfI	SwaI	

Table 2. Restriction Enzymes That Cleave TR201 Once

Aasl	3216	Acc16l	4220	Acvl	715	Ajil	942	Alel	6556	Alol	7848	Apal	516	Aspl	1047
Asull	8300	Avill	4220	Axyl	2339	Bael	5338	BamHl	7879	BbrPl	715	BbvCl	6403	BmgBl	942
Bpu14l	8300	Bsa29l	8211	BsaAl	715	BsaBl	7840	Bse21l	2339	Bse8l	7840	BseCl	8211	BseJl	7840
Bshv1l	8211	BsiWl	1061	Bsp119l	8300	Bsp120l	512	BspDl	8211	BspT104l	8300	BstBAI	715	BstBl	8300
BstElI	1139	BstENl	6145	BstPl	1139	Bsu15l	8211	Bsu36l	2339	Btrl	942	Clal	8211	Cpol	1121
Cspl	1121	Drdl	3216	DseDl	3216	Eco32l	8532	Eco72l	715	Eco81l	2339	Eco91l	1139	EcoNl	6145
EcoO65l	1139	EcoRV	8532	Fsel	6125	Fspl	4220	Kfl	6913	Mlul	1605	Nsbl	4220	NspV	8300
Olil	6556	PaeR7l	7308	Pfl23ll	1061	PflFl	1047	PmaCl	715	Pmll	715	Ppu21l	715	PpuMl	6913
Psp5Il	6913	PspCl	715	PspEl	1139	PspLl	1061	PspOMl	512	PspPPl	6913	PspXl	7308	Pstl	2223
Psyl	1047	Rigl	6125	Rsr2l	1121	Rsrll	1121	Sfl	8106	Sfp274l	7308	Sfl	8300	SgrDl	4919
Slal	7308	Sspl	4802	Strl	7308	Tth111l	1047	Xagl	6145	Xcml	5144	Xhol	7308		

Table 3. Restriction Enzymes That Cleave TR201 Twice

Aanl	2684	Aanl	4957	Aarl	738	Aarl	8137	Adel	762	Adel	1576	Aflll	2532	Aflll	5492
Ahdl	3998	Ahdl	7569	AspA2l	324	AspA2l	550	AspEl	3998	AspEl	7569	Assl	4478	Assl	8206
AsuNHI	2720	AsuNHI	8080	Avrll	324	Avrll	550	Ball	1380	Ball	7777	Baul	3281	Baul	4662
Bcgl	1446	Bcgl	4503	Bfrl	2532	Bfrl	5492	Blnl	324	Blnl	550	BmcAl	4478	BmcAl	8206
BmeRl	3998	BmeRl	7569	Bmtl	2724	Bmtl	8084	Bpu10l	6403	Bpu10l	8089	BsePl	1460	BsePl	5686
BsmBl	986	BsmBl	1397	Bsp68l	5808	Bsp68l	7765	BspHl	3825	BspHl	4833	BspOl	2724	BspOl	8084
BspQl	2992	BspQl	6077	BspTl	2532	BspTl	5492	BssHll	1460	BssHll	5686	BssSl	3281	BssSl	4662
Bst2Bl	3281	Bst2Bl	4662	BstAFI	2532	BstAFI	5492	BtuMl	5808	BtuMl	7765	Ccil	3825	Ccil	4833
Cfr9l	345	Cfr9l	1054	Csil	91	Csil	1572	Dralll	762	Dralll	1576	Dril	3998	Dril	7569
Eam1105l	3998	Eam1105l	7569	Eco147l	323	Eco147l	1336	EcoRl	7949	EcoRl	8273	EcoT22l	74	EcoT22l	146
Esp3l	986	Esp3l	1397	Krol	2130	Krol	6121	Lgul	2992	Lgul	6077	Mabl	91	Mabl	1572
Misl	1380	Misl	7777	MluNl	1380	MluNl	7777	Mph1103l	74	Mph1103l	146	MroNl	2130	MroNl	6121
MscI	1380	MscI	7777	Msp20l	1380	Msp20l	7777	MspCl	2532	MspCl	5492	Nael	2132	Nael	6123
NgoMIV	2130	NgoMIV	6121	Nhel	2720	Nhel	8080	Nrul	5808	Nrul	7765	Nsil	74	Nsil	146
Pagl	3825	Pagl	4833	Paul	1460	Paul	5686	Pcel	323	Pcel	1336	Pcil	889	Pcil	3108
PciSl	2992	PciSl	6077	Pdil	2132	Pdil	6123	Pfol	1480	Pfol	1891	PscI	889	PscI	3108
Psil	2684	Psil	4957	Ptel	1460	Ptel	5686	Pvull	2376	Pvull	2932	Rcal	3825	Rcal	4833
Rrul	5808	Rrul	7765	Sapl	2992	Sapl	6077	Scal	4478	Scal	8206	SexAl	91	SexAl	1572
Smal	347	Smal	1056	SseBl	323	SseBl	1336	Stul	323	Stul	1336	TspMl	345	TspMl	1054
Vha464l	2532	Vha464l	5492	Xbal	7974	Xbal	7981	Xmal	345	Xmal	1054	XmaJl	324	XmaJl	550
Zrml	4478	Zrml	8206	Zsp2l	74	Zsp2l	146								

Table 4. Restriction Enzymes That Cleave TR201 Three Times

[illegible]