

TN-ROSA26-L Datasheet

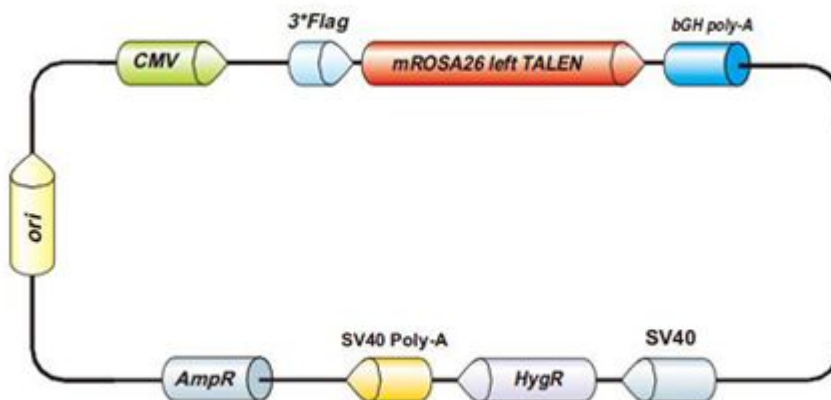
Clone Information

Catalog No.: TN-ROSA26-L

Description: Left TALEN clone for mouse ROSA26 safe harbor site

Whole Plasmid Size: 8016 bp

Resistance: Hygromycin



Restriction Enzyme Information for TN-ROSA26-L

Table 1. Restriction Enzymes That Do Not Cleave TN-ROSA26-L

Acc65I AgeI AscI BaeI BaeI BamHI BbsI BbvCI
 BglII BlnI BsaI BsaBI BsgI BsiWI BsrGI BstBI
 BstEII Bsu36I BtrI ClaI EcoNI FseI HindIII HpaI
 KpnI MscI NheI NotI PacI PflMI PmeI PpuMI
 SanDI SapI SbfI SfiI SgrAI SrfI SwaI TliI
 XbaI XhoI

Table 2. Restriction Enzymes That Cleave TN-ROSA26-L Once

Acc16I	7273 AflII	961 AhiI	1473 ApaI	974 ArsI	6477 AsiSI	6013 AspA2I	5585 AviII	7273
AvrII	5585 BcuI	1473 BfrI	961 BlnI	5585 BoxI	5676 Bpu10I	4178 BsePI	32 BsmBI	5717
BsmI	6884 Bsp120I	970 BspTI	961 BssHII	32 BssNAI	6936 Bst1107I	6936 Bst6I	4429 BstAFI	961
BstPAI	5676 BstXI	4784 BstZ17I	6936 Cfr42I	6429 Cfr9I	5606 CpoI	6057 CsiI	5352 CspCI	1850
CspI	6057 DniI	1093 Eam1104I	4429 EarI	4429 Ecl136II	2040 Eco147I	5584 Eco32I	1314 Eco53kI	2040
EcoICRI	2040 EcoRI	5892 EcoRV	1314 EgeI	1093 EheI	1093 Esp3I	5717 FspI	7273 KasI	1091
KroI	6747 KspI	6429 Mabi	5352 MauBI	32 MluI	2603 Mly113I	1092 MroNI	6747 MspCI	961
Mva1269I	6884 NaeI	6749 NarI	1092 NgoMIV	6747 NsbI	7273 Pasi	976 Paul	32 PciI	5584
PctI	6884 PdiI	6749 PfoI	6739 PluTI	1095 PshAI	5676 Psp124BI	2042 PspOMI	970 PteI	32
PvuII	2538 RgaI	6013 Rsr2I	6057 RsrII	6057 SacI	2042 SacII	6429 SexAI	5352 SfaI	6013
SfoI	1093 Sfr303I	6429 SgfI	6013 SgrBI	6429 Smal	5608 SpeI	1473 SseBI	5584 SspDI	1091
SstI	2042 StuI	5584 TspMI	5606 Vha464I	961 XcmI	4784 XmaI	5606 XmaJI	5585	

Table 3. Restriction Enzymes That Cleave TN-ROSA26-L Twice

AcvI	5645 AcvI	6677 Adel	5937 Adel	6230 AhdI	956 AhdI	7051 Ajul	4916 Ajul	6362
AlwNI	321 AlwNI	2556 Asp700I	5257 Asp700I	7650 AspEI	956 AspEI	7051 Aspl	5709 Aspl	6153
AssI	6619 AssI	7531 BbrPI	5645 BbrPI	6677 BcgI	1425 BcgI	7556 BmcAI	6619 BmcAI	7531
BmeRI	956 BmeRI	7051 Bmri	1763 Bmri	5464 Bmul	1763 Bmul	5464 Bse3DI	2619 Bse3DI	4905
BseMI	2619 BseMI	4905 Bsp68I	2566 Bsp68I	7975 BsrDI	2619 BsrDI	4905 BstSNI	1814 BstSNI	4485
BtuMI	2566 BtuMI	7975 Cail	321 Cail	2556 DraI	1160 DraI	7628 DrallI	5937 DrallI	6230

Dril	956	Dril	7051	Eam1105I	956	Eam1105I	7051	Eco105I	1814	Eco105I	4485	Eco72I	5645	Eco72I	6677
EcoT22I	5335	EcoT22I	5407	FauNDI	1708	FauNDI	6099	Mph1103I	5335	Mph1103I	5407	MroXI	5257	MroXI	7650
NdeI	1708	NdeI	6099	Nrul	2566	Nrul	7975	Nsil	5335	Nsil	5407	PaeI	5333	PaeI	5405
PciI	730	PciI	1326	Pdml	5257	Pdml	7650	PfiFI	5709	PfiFI	6153	PmaCI	5645	PmaCI	6677
PmlI	5645	PmlI	6677	Psci	730	Psci	1326	PspCI	5645	PspCI	6677	PstI	2379	PstI	5986
PstNI	321	PstNI	2556	Psyl	5709	Psyl	6153	Rrul	2566	Rrul	7975	Scal	6619	Scal	7531
SnaBI	1814	SnaBI	4485	SphI	5333	SphI	5405	Tth111I	5709	Tth111I	6153	XmnI	5257	XmnI	7650
Zrml	6619	Zrml	7531	Zsp2I	5335	Zsp2I	5407								

Table 4. Restriction Enzymes That Cleave TN-ROSA26-L Three Times

AanI	1134	AanI	5187	AanI	6823	AasI	628	AasI	6153	AasI	6534	AccI	4734	AccI	7277
AccI	7650	AflIII	730	AflIII	1326	AflIII	2603	BclI	4284	BclI	4946	BclI	5636	BpvUI	2450
BpvUI	6013	BpvUI	7421	BsaXI	4471	BsaXI	4685	BsaXI	5697	Bse118I	5987	Bse118I	6747	Bse118I	7131
BseRI	2591	BseRI	4462	BseRI	5581	BseX3I	5854	BseX3I	6019	BseX3I	6589	BspHI	2	BspHI	2148
BspHI	7886	BsrFI	5987	BsrFI	6747	BsrFI	7131	BssAI	5987	BssAI	6747	BssAI	7131	BstZI	5854
BstZI	6019	BstZI	6589	CciI	2	CciI	2148	CciI	7886	Cfr10I	5987	Cfr10I	6747	Cfr10I	7131
DrdI	628	DrdI	6153	DrdI	6534	DseDI	628	DseDI	6153	DseDI	6534	EagI	5854	EagI	6019
EagI	6589	EclXI	5854	EclXI	6019	EclXI	6589	Eco52I	5854	Eco52I	6019	Eco52I	6589	FbaI	4284
FbaI	4946	FbaI	5636	Ksp22I	4284	Ksp22I	4946	Ksp22I	5636	MvrI	2450	MvrI	6013	MvrI	7421
PagI	2	PagI	2148	PagI	7886	Ple19I	2450	Ple19I	6013	Ple19I	7421	Psil	1134	Psil	5187
Psil	6823	Psp1406I	4734	Psp1406I	7277	Psp1406I	7650	PvuI	2450	PvuI	6013	PvuI	7421	RcaI	2
RcaI	2148	RcaI	7886	Sall	2295	Sall	4672	Sall	6941	SspI	1083	SspI	1244	SspI	7855
