

TN-ROSA26-R Datasheet

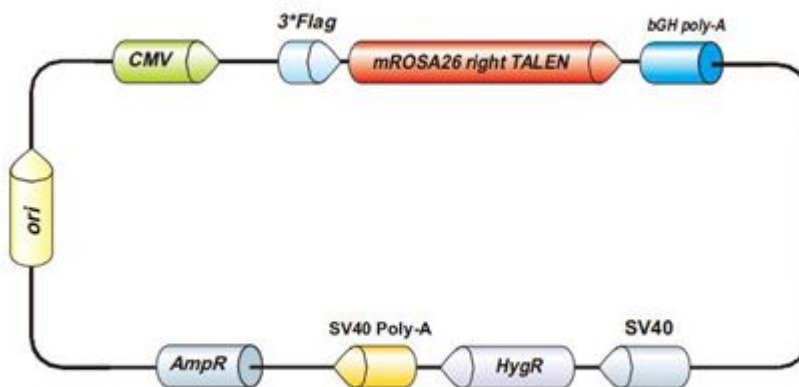
Clone Information

Catalog No.: TN-ROSA26-R

Description: Right TALEN clone for mouse ROSA26 safe harbor site

Whole Plasmid Size: 7914 bp

Resistance: Hygromycin



Restriction Enzyme Information for TN-ROSA26-R

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

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-R Once

Acc16I	7171 AflII	961 AhiI	1473 ApaI	974 ArsI	6375 AsiSI	5911 AspA2I	5483 AviII	7171
AvrII	5483 BcuI	1473 BfrI	961 BlnI	5483 BoxI	5574 Bpu10I	4076 BsePI	32 BsmBI	5615
BsmI	6782 Bsp120I	970 BspTI	961 BssHII	32 BssNAI	6834 Bst1107I	6834 Bst6I	4327 BstAFI	961
BstPAI	5574 BstXI	4682 BstZ17I	6834 Cfr42I	6327 Cfr9I	5504 CpoI	5955 CsiI	5250 CspCI	1850
CspI	5955 DniI	1093 Eam1104I	4327 EarI	4327 Ecl136II	2040 Eco147I	5482 Eco32I	1314 Eco53kI	2040
EcoICRI	2040 EcoRI	5790 EcoRV	1314 EgeI	1093 EheI	1093 Esp3I	5615 FspI	7171 KasI	1091
KroI	6645 KspI	6327 MabI	5250 MauBI	32 MluI	2603 Mly113I	1092 MroNI	6645 MspCI	961
Mva1269I	6782 NaeI	6647 NarI	1092 NgoMIV	6645 NsbI	7171 PaeI	976 PaeI	32 PceI	5482
PctI	6782 PdiI	6647 PfoI	6637 PluTI	1095 PshAI	5574 Psp124BI	2042 PspOMI	970 PteI	32
PvuII	2538 RgaI	5911 Rsr2I	5955 RsrII	5955 SacI	2042 SacII	6327 SexAI	5250 SfaI	5911
SfoI	1093 Sfr303I	6327 SgfI	5911 SgrBI	6327 Smal	5506 SpeI	1473 SseBI	5482 SspDI	1091
SstI	2042 StuI	5482 TspMI	5504 Vha464I	961 XcmI	4682 XmaI	5504 XmaJI	5483	

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

AcvI	5543 AcvI	6575 Adel	5835 Adel	6128 AhdI	956 AhdI	6949 AjuI	4814 AjuI	6260
AlwNI	321 AlwNI	2556 Asp700I	5155 Asp700I	7548 AspEI	956 AspEI	6949 Aspl	5607 Aspl	6051
AssI	6517 AssI	7429 BbrPI	5543 BbrPI	6575 BcgI	1425 BcgI	7454 BmcAI	6517 BmcAI	7429
BmeRI	956 BmeRI	6949 Bmrl	1763 Bmrl	5362 Bmul	1763 Bmul	5362 Bse3DI	2619 Bse3DI	4803
BseMI	2619 BseMI	4803 Bsp68I	2566 Bsp68I	7873 BsrDI	2619 BsrDI	4803 BstSNI	1814 BstSNI	4383
BtuMI	2566 BtuMI	7873 Cail	321 Cail	2556 DraI	1160 DraI	7526 DrallI	5835 DrallI	6128

Dril	956	Dril	6949	Eam1105I	956	Eam1105I	6949	Eco105I	1814	Eco105I	4383	Eco72I	5543	Eco72I	6575
EcoT22I	5233	EcoT22I	5305	FauNDI	1708	FauNDI	5997	Mph1103I	5233	Mph1103I	5305	MroXI	5155	MroXI	7548
NdeI	1708	NdeI	5997	Nrul	2566	Nrul	7873	Nsil	5233	Nsil	5305	PaeI	5231	PaeI	5303
PciI	730	PciI	1326	Pdml	5155	Pdml	7548	PfiFI	5607	PfiFI	6051	PmaCI	5543	PmaCI	6575
PmlI	5543	PmlI	6575	Psci	730	Psci	1326	PspCI	5543	PspCI	6575	PstI	2379	PstI	5884
PstNI	321	PstNI	2556	Psyl	5607	Psyl	6051	Rrul	2566	Rrul	7873	Scal	6517	Scal	7429
SnaBI	1814	SnaBI	4383	SphI	5231	SphI	5303	Tth111I	5607	Tth111I	6051	XmnI	5155	XmnI	7548
Zrml	6517	Zrml	7429	Zsp2I	5233	Zsp2I	5305								

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

AanI	1134	AanI	5085	AanI	6721	AasI	628	AasI	6051	AasI	6432	AccI	4632	AccI	7175
AccI	7548	AflIII	730	AflIII	1326	AflIII	2603	BclI	4182	BclI	4844	BclI	5534	BpvUI	2450
BpvUI	5911	BpvUI	7319	BsaXI	4369	BsaXI	4583	BsaXI	5595	Bse118I	5885	Bse118I	6645	Bse118I	7029
BseRI	2591	BseRI	4360	BseRI	5479	BseX3I	5752	BseX3I	5917	BseX3I	6487	BspHI	2	BspHI	2148
BspHI	7784	BsrFI	5885	BsrFI	6645	BsrFI	7029	BssAI	5885	BssAI	6645	BssAI	7029	BstZI	5752
BstZI	5917	BstZI	6487	CciI	2	CciI	2148	CciI	7784	Cfr10I	5885	Cfr10I	6645	Cfr10I	7029
DrdI	628	DrdI	6051	DrdI	6432	DseDI	628	DseDI	6051	DseDI	6432	EagI	5752	EagI	5917
EagI	6487	EclXI	5752	EclXI	5917	EclXI	6487	Eco52I	5752	Eco52I	5917	Eco52I	6487	FbaI	4182
FbaI	4844	FbaI	5534	Ksp22I	4182	Ksp22I	4844	Ksp22I	5534	MvrI	2450	MvrI	5911	MvrI	7319
PagI	2	PagI	2148	PagI	7784	Ple19I	2450	Ple19I	5911	Ple19I	7319	Psil	1134	Psil	5085
Psil	6721	Psp1406I	4632	Psp1406I	7175	Psp1406I	7548	PvuI	2450	PvuI	5911	PvuI	7319	RcaI	2
RcaI	2148	RcaI	7784	Sall	2295	Sall	4570	Sall	6839	SspI	1083	SspI	1244	SspI	7753
