

MCP-ROSA26-CG01 Datasheet

Clone Information

Catalog No.: MCP-ROSA26-CG01

Description: All-in-one sgRNA/Cas9 expression clone for mouse ROSA26 safe harbor site

Target site: GACTCCAGTCTTTCTAGAAGA

Whole Plasmid Size: 10442 bp

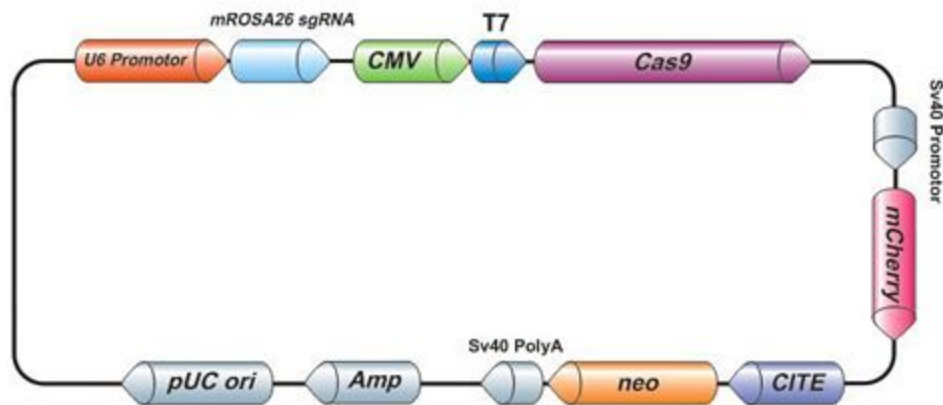
Resistance: Ampicillin

Suggested Sequencing Primers:

Forward: 5'-TTCTTGGGTAGTTTGCAG-3'

Reverse: 5'-CCTATTGGCGTTACTATG-3'

Vector Information



Restriction Enzyme Information for MCP-ROSA26-CG01

Table 1. Restriction Enzymes That Do Not Cleave MCP-ROSA26-CG01

AflIII BsaBI BsiWI BsmBI BstBI Bsu36I ClaI EcoRV
 FseI HpaI NheI PacI PmeI SfiI SgfI SrfI
 SwaI

Table 2. Restriction Enzymes That Cleave MCP-ROSA26-CG01 Once

AanI	8059	AccIII	5299	AhlI	537	Aor13HI	5299	AscI	5748	AspI	7360	BaeI	1206	BamHI	333
BbvCI	6308	BcuI	537	BoxI	343	BpvUI	8730	BseAI	5299	Bsp13I	5299	Bsp68I	496	BspEI	5299
BstPAI	343	BstSNI	878	BtuMI	496	CciNI	6482	Cfr9I	5712	CpoI	7758	CsiI	5458	CspCI	914
CspI	7758	Ecl136II	1104	Eco105I	878	Eco53kI	1104	EcoICRI	1104	EcoRI	1	KfiI	4566	Kpn2I	5299
MabI	5458	MfeI	449	MluI	516	MreI	4837	MroI	5299	MunI	449	MvrI	8730	NotI	6482
NruI	496	PaeR7I	5362	PalAI	5748	PfiFI	7360	Ple19I	8730	PshAI	343	PsiI	8059	Psp124BI	1106
PspXI	5362	PsyI	7360	PvuI	8730	RruI	496	Rsr2I	7758	RsrII	7758	SacI	1106	Sall	8177
SexAI	5458	Sfr274I	5362	Sgsl	5748	Slal	5362	SmaI	5714	SnaBI	878	SpeI	537	SspI	8294
SstI	1106	StrI	5362	TspMI	5712	Tth111I	7360	XhoI	5362	XmaI	5712				

Table 3. Restriction Enzymes That Cleave MCP-ROSA26-CG01 Twice

AarI	1607 AarI	6835 Acc65I	1222 Acc65I	6937 AcsI	1 AcsI	8090 AfeI	2158 AfeI	5191
AgeI	5245 AgeI	5757 AhdI	5964 AhdI	9099 Ajul	3279 Ajul	4314 AleI	5770 AleI	6412
Aor51HI	2158 Aor51HI	5191 Apal	2548 Apal	6613 Apol	1 Apol	8090 AsiGI	5245 AsiGI	5757
Asp718I	1222 Asp718I	6937 AspA2I	5691 AspA2I	6647 AspEI	5964 AspEI	9099 AvrII	5691 AvrII	6647
BciVI	7461 BciVI	9789 BfuI	7461 BfuI	9789 BglII	1439 BglII	2234 BlnI	5691 BlnI	6647
BmeRI	5964 BmeRI	9099 BshTI	5245 BshTI	5757 Bsp120I	2544 Bsp120I	6609 Bsp1407I	2747 Bsp1407I	6470
BspHI	4313 BspHI	9267 BsrGI	2747 BsrGI	6470 BssNAI	5328 BssNAI	8172 Bst1107I	5328 Bst1107I	8172
BstAU1	2747 BstAU1	6470 BstZ17I	5328 BstZ17I	8172 Bsul	7461 Bsul	9789 CciI	4313 CciI	9267
CspAI	5245 CspAI	5757 Dril	5964 Dril	9099 Eam1105I	5964 Eam1105I	9099 Eco47III	2158 Eco47III	5191
EcoT22I	5441 EcoT22I	5513 FauNDI	268 FauNDI	772 KpnI	1226 KpnI	6941 Mph1103I	5441 Mph1103I	5513
NdeI	268 NdeI	772 NsiI	5441 NsiI	5513 OiiI	5770 OiiI	6412 PagI	4313 PagI	9267
PinAI	5245 PinAI	5757 PspOMI	2544 PspOMI	6609 Rcal	4313 Rcal	9267 SbfI	1436 SbfI	6125
SdaI	1436 SdaI	6125 SgrAI	4837 SgrAI	6451 Sse8387I	1436 Sse8387I	6125 XapI	1 XapI	8090
XbaI	351 XbaI	441 XmaJI	5691 XmaJI	6647 Zsp2I	5441 Zsp2I	5513		

Table 4. Restriction Enzymes That Cleave MCP-ROSA26-CG01 Three Times

AasI	6310 AasI	7269 AasI	9885 AccI	5327 AccI	8171 AccI	8178 AclI	6519 AclI	8497
AclI	8870 AcvI	40 AcvI	3978 AcvI	6812 Asp700I	4321 Asp700I	6708 Asp700I	8499 BbrPI	40
BbrPI	3978 BbrPI	6812 BsmI	6642 BsmI	6681 BsmI	8120 BstAPI	2724 BstAPI	5436 BstAPI	5508
BstEII	2862 BstEII	5925 BstEII	6096 BstPI	2862 BstPI	5925 BstPI	6096 BstXI	1233 BstXI	1803
BstXI	3978 Cfr42I	1683 Cfr42I	2559 Cfr42I	3780 DrdI	6310 DrdI	7269 DrdI	9885 DseDI	6310
DseDI	7269 DseDI	9885 Eco147I	5690 Eco147I	5995 Eco147I	6217 Eco72I	40 Eco72I	3978 Eco72I	6812
Eco91I	2862 Eco91I	5925 Eco91I	6096 EcoO65I	2862 EcoO65I	5925 EcoO65I	6096 FbII	5327 FbII	8171
FbII	8178 HincII	522 HincII	6369 HincII	8179 HindII	522 HindII	6369 HindII	8179 KspI	1683
KspI	2559 KspI	3780 MroXI	4321 MroXI	6708 MroXI	8499 Mva1269I	6642 Mva1269I	6681 Mva1269I	8120
PceI	5690 PceI	5995 PceI	6217 PciI	3650 PciI	6986 PciI	9987 PctI	6642 PctI	6681
PctI	8120 PdmI	4321 PdmI	6708 PdmI	8499 PmaCI	40 PmaCI	3978 PmaCI	6812 PmII	40
PmII	3978 PmII	6812 PscI	3650 PscI	6986 PscI	9987 Psp1406I	6519 Psp1406I	8497 Psp1406I	8870
PspCI	40 PspCI	3978 PspCI	6812 PspEI	2862 PspEI	5925 PspEI	6096 PsrI	3997 PsrI	4812
PsrI	4879 SacII	1683 SacII	2559 SacII	3780 Sfr303I	1683 Sfr303I	2559 Sfr303I	3780 SgrBI	1683
SgrBI	2559 SgrBI	3780 SseBI	5690 SseBI	5995 SseBI	6217 Stul	5690 Stul	5995 Stul	6217
XmI	5327 XmI	8171 XmI	8178 XmnI	4321 XmnI	6708 XmnI	8499		
