

DC-RFP-SH02 Datasheet

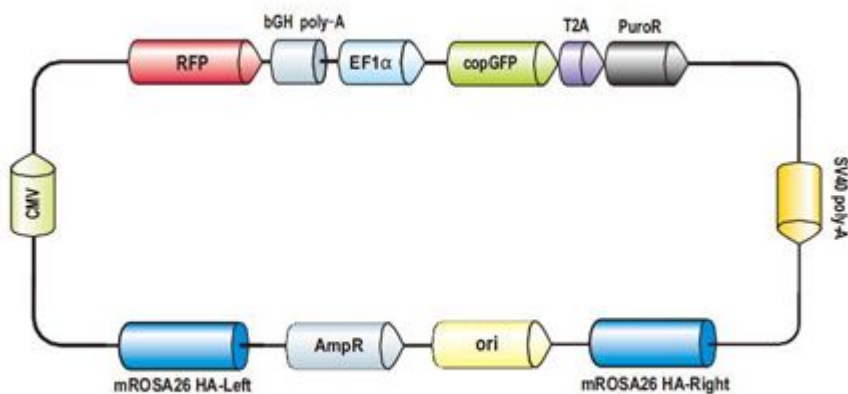
Clone Information

Catalog No.: DC-RFP-SH02

Description: ROSA26 RFP control

Whole Plasmid Size: 9355 bp

Resistance: Ampicillin



Restriction Enzyme Information for DC-RFP-SH02

Table 1. Restriction Enzymes That Do Not Cleave DC-RFP-SH02

AccI AfelAflII Agel AscI AvrII BaeI BaeI
 BbvCI BclI BstZ17I ClaI NheI PacI PmeI PpuMI
 PshAI SalI SanDI SgrAI

Table 2. Restriction Enzymes That Cleave DC-RFP-SH02 Once

AarI	4579	Acc65I	1396	AhlI	1359	AjuI	1376	Alol	289	ArsI	3958	AsiSI	3581	Asp718I	1396
AspI	4980	AsuII	3311	AxyI	3863	BamHI	3324	BcuI	1359	BglII	3555	Bpu10I	3957	Bpu14I	3311
BsaBI	2223	Bse21I	3863	Bse8I	2223	BseJI	2223	Bsp119I	3311	Bsp1407I	6021	BspT104I	3311	BsrGI	6021
BstAPI	5509	BstAUI	6021	BstBI	3311	BstEI	5072	BstPI	5072	BstSNI	2312	Bsu36I	3863	Cfr42I	5152
Ecl136II	2538	Eco105I	2312	Eco53kI	2538	Eco81I	3863	Eco91I	5072	EcoICRI	2538	EcoO65I	5072	FauNDI	2660
FseI	4471	KpnI	1400	KspI	5152	MauBI	1556	MfeI	3653	MluI	2197	MunI	3653	NdeI	2660
NspV	3311	PciI	7713	PfiFI	4980	PfoI	5413	PscI	7713	Psp124BI	2540	PspEI	5072	PspXI	2242
Psrl	2953	PsyI	4980	RgaI	3581	RigI	4471	SacI	2540	SacII	5152	SbfI	6089	SdaI	6089
SfaAI	3581	SfiI	1352	Sfr303I	5152	SfuI	3311	SgfI	3581	SgrBI	5152	SmiI	3318	SnaBI	2312
SpeI	1359	Sse8387I	6089	SstI	2540	Swal	3318	Tth111I	4980	XcmI	4841				

Table 3. Restriction Enzymes That Cleave DC-RFP-SH02 Twice

AasI	310	AasI	7821	Acc36I	4306	Acc36I	4579	AccB7I	1455	AccB7I	4313	AcvI	4310	AcvI	4442
AflIII	2197	AflIII	7713	AhdI	2790	AhdI	8606	AjiI	1492	AjiI	4916	AlfI	3556	AlfI	3590
Alw44I	8027	Alw44I	9273	AlwNI	1770	AlwNI	8129	ApaLI	8027	ApaLI	9273	AspEI	2790	AspEI	8606
AssI	6399	AssI	9086	Ball	5313	Ball	7092	BasI	1455	BasI	4313	BbrPI	4310	BbrPI	4442
BfuAI	4306	BfuAI	4579	BmcAI	6399	BmcAI	9086	BmeRI	2790	BmeRI	8606	BmgBI	1492	BmgBI	4916
BsiWI	899	BsiWI	4994	BsmBI	2561	BsmBI	5330	BsmI	5725	BsmI	6809	Bsp68I	4882	Bsp68I	6350
BspMI	4306	BspMI	4579	BtrI	1492	BtrI	4916	BtuMI	4882	BtuMI	6350	BveI	4306	BveI	4579
CaiI	1770	CaiI	8129	Cfr9I	1187	Cfr9I	1251	CpoI	1203	CpoI	5054	Cspl	1203	Cspl	5054
DrdI	310	DrdI	7821	Dril	2790	Dril	8606	DseDI	310	DseDI	7821	Eam1105I	2790	Eam1105I	8606
Eco32I	2221	Eco32I	2940	Eco72I	4310	Eco72I	4442	EcoRI	1384	EcoRI	2209	EcoRV	2221	EcoRV	2940
EcoT22I	2253	EcoT22I	6848	Esp3I	2561	Esp3I	5330	Fall	6772	Fall	6804	HindIII	3816	HindIII	6781

HpaI	1406 HpaI	5644 KspAI	1406 KspAI	5644 MlsI	5313 MlsI	7092 MluNI	5313 MluNI	7092
Mph1103I	2253 Mph1103I	6848 MscI	5313 MscI	7092 Msp20I	5313 Msp20I	7092 Mva1269I	5725 Mva1269I	6809
NruI	4882 NruI	6350 NsiI	2253 NsiI	6848 PaeR7I	2242 PaeR7I	6838 PctI	5725 PctI	6809
Pfi23II	899 Pfi23II	4994 PfiMI	1455 PfiMI	4313 PmaCI	4310 PmaCI	4442 PmlI	4310 PmlI	4442
PspCI	4310 PspCI	4442 PspLI	899 PspLI	4994 PstNI	1770 PstNI	8129 RruI	4882 RruI	6350
Rsr2I	1203 Rsr2I	5054 RsrII	1203 RsrII	5054 Scal	6399 Scal	9086 Sfr274I	2242 Sfr274I	6838
Slal	2242 Slal	6838 Smal	1189 Smal	1253 StrI	2242 StrI	6838 TspMI	1187 TspMI	1251
Van91I	1455 Van91I	4313 VneI	8027 VneI	9273 XbaI	2191 XbaI	2605 XhoI	2242 XhoI	6838
XmaI	1187 XmaI	1251 Zrml	6399 Zrml	9086 Zsp2I	2253 Zsp2I	6848		

Table 4. Restriction Enzymes That Cleave DC-RFP-SH02 Three Times

AatII	2419 AatII	4984 AatII	5482 AclI	3775 AclI	8832 AclI	9205 ApaI	1648 ApaI	4716
ApaI	6009 BbsI	3519 BbsI	4890 BbsI	6032 BplI	1836 BplI	1898 BplI	4302 BpII	3519
BpII	4890 BpII	6032 Bpu1102I	1836 Bpu1102I	1898 Bpu1102I	4302 BpuAI	3519 BpuAI	4890 BpuAI	6032
BpvUI	628 BpvUI	3581 BpvUI	8976 Bse3DI	3113 Bse3DI	8667 Bse3DI	8841 BseMI	3113 BseMI	8667
BseMI	8841 Bsp120I	1644 Bsp120I	4712 Bsp120I	6005 Bsp1720I	1836 Bsp1720I	1898 Bsp1720I	4302 BspHI	48
BspHI	6555 BspHI	8433 BspQI	2908 BspQI	4257 BspQI	7597 BsrDI	3113 BsrDI	8667 BsrDI	8841
BstV2I	3519 BstV2I	4890 BstV2I	6032 CciI	48 CciI	6555 CciI	8433 CciNI	1565 CciNI	2236
CciNI	3562 CeiII	1836 CeiII	1898 CeiII	4302 CsiI	5505 CsiI	5908 CsiI	6064 CspCI	149
CspCI	2348 CspCI	3900 Eco147I	1346 Eco147I	1355 Eco147I	5269 HincII	1406 HincII	5644 HincII	5906
HindII	1406 HindII	5644 HindII	5906 LguI	2908 LguI	4257 LguI	7597 MabI	5505 MabI	5908
MabI	6064 Mvrl	628 Mvrl	3581 Mvrl	8976 NotI	1565 NotI	2236 NotI	3562 PaeI	2251
PaeI	3530 PaeI	6850 PagI	48 PagI	6555 PagI	8433 Pasi	1138 Pasi	2134 Pasi	4987
PceI	1346 PceI	1355 PceI	5269 PciSI	2908 PciSI	4257 PciSI	7597 Ple19I	628 Ple19I	3581
Ple19I	8976 Psp1406I	3775 Psp1406I	8832 Psp1406I	9205 PspOMI	1644 PspOMI	4712 PspOMI	6005 Pvul	628
Pvul	3581 Pvul	8976 Rcal	48 Rcal	6555 Rcal	8433 SapI	2908 SapI	4257 SapI	7597
SexAI	5505 SexAI	5908 SexAI	6064 SphI	2251 SphI	3530 SphI	6850 SseBI	1346 SseBI	1355
SseBI	5269 Sspl	17 Sspl	147 Sspl	5855 Stul	1346 Stul	1355 Stul	5269 TspGWI	2940
TspGWI	9046 TspGWI	9388 Zral	2417 Zral	4982 Zral	5480			