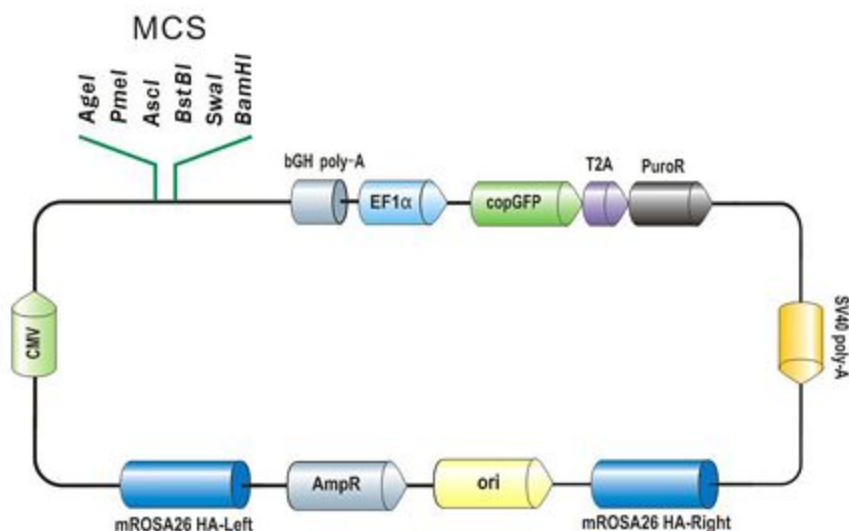


DC-DON-SH02 Datasheet

Clone Information

Catalog No.: DC-DON-SH02
Description: ROSA26 donor vector
Whole Plasmid Size: 8715 bp
Resistance: Ampicillin

Vector Information for DC-DON-SH02



Restriction Enzyme Information for DC-DON-SH02

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

AccI Afel AfII AvrII BaeI BaeII BbvCI BclI
 BstZ17I ClaI NdeI NheI PaeI PpuMI PshAI SalI
 SanDI SgrAI

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

AarI	3901	Acc65I	1396	AgeI	2610	AhdI	7928	AhlI	1359	AjuI	1376	AlaI	289	ArS	3280
AscI	2625	AsiGI	2610	AsiSI	2903	Asp718I	1396	AspEI	7928	AspI	4302	AsuII	2633	AxyI	3185
BamHI	2646	BcuI	1359	BglII	2877	BmeRI	7928	Bpu10I	3279	Bpu14I	2633	BsaBI	2222	Bse21I	3185
Bse8I	2222	BseJI	2222	BshTI	2610	Bsp119I	2633	Bsp1407I	5343	BspT104I	2633	BsrGI	5343	BstAPI	4831
BstAU	5343	BstBI	2633	BstEI	4394	BstPI	4394	BstSNI	2311	Bsu36I	3185	Cfr42I	4474	CspAI	2610
DrI	7928	Eam1105I	7928	Ecl136II	2537	Eco105I	2311	Eco32I	2220	Eco53kI	2537	Eco81I	3185	Eco91I	4394
EcoICRI	2537	EcoO65I	4394	EcoRV	2220	FseI	3793	KpnI	1400	KspI	4474	MauBI	1556	MfeI	2975
MluI	2196	MssI	2619	MunI	2975	NspV	2633	PalAI	2625	PciI	7035	PfiFI	4302	PfoI	4735
PinAI	2610	PmeI	2619	PscI	7035	Psp124BI	2539	PspEI	4394	PspXI	2241	PsyI	4302	RgaI	2903
RigI	3793	SacI	2539	SacII	4474	SbfI	5411	SdaI	5411	SfaAI	2903	SfiI	1352	Sfr303I	4474
SfuI	2633	SgfI	2903	SgrBI	4474	SgsI	2625	SmiI	2640	SnaBI	2311	SpeI	1359	Sse8387I	5411
SstI	2539	Swal	2640	Tth111I	4302	XcmI	4163								

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

AasI	310	AasI	7143	Acc36I	3628	Acc36I	3901	AccB7I	1455	AccB7I	3635	AcvI	3632	AcvI	3764
AflIII	2196	AflIII	7035	AjiI	1492	AjiI	4238	AlfI	2878	AlfI	2912	Alw44I	7349	Alw44I	8595
AlwNI	1770	AlwNI	7451	ApaLI	7349	ApaLI	8595	AssI	5721	AssI	8408	Ball	4635	Ball	6414
BasI	1455	BasI	3635	BbrPI	3632	BbrPI	3764	BfuAI	3628	BfuAI	3901	BmcAI	5721	BmcAI	8408
BmgBI	1492	BmgBI	4238	BpII	736	BpII	768	Bse3DI	7989	Bse3DI	8163	BseMI	7989	BseMI	8163
BsiWI	899	BsiWI	4316	BsmBI	2560	BsmBI	4652	BsmI	5047	BsmI	6131	Bsp68I	4204	Bsp68I	5672
BspMI	3628	BspMI	3901	BspQI	3579	BspQI	6919	BsrDI	7989	BsrDI	8163	BtrI	1492	BtrI	4238
BtuMI	4204	BtuMI	5672	Bvel	3628	Bvel	3901	CaiI	1770	CaiI	7451	Cfr9I	1187	Cfr9I	1251
CpoI	1203	CpoI	4376	Cspl	1203	Cspl	4376	DrdI	310	DrdI	7143	DseDI	310	DseDI	7143
Eco72I	3632	Eco72I	3764	EcoRI	1384	EcoRI	2208	EcoT22I	2252	EcoT22I	6170	Esp3I	2560	Esp3I	4652
FalI	6094	FalI	6126	HindIII	3138	HindIII	6103	HpaI	1406	HpaI	4966	KspAI	1406	KspAI	4966
LguI	3579	LguI	6919	MisI	4635	MisI	6414	MluNI	4635	MluNI	6414	Mph1103I	2252	Mph1103I	6170
MscI	4635	MscI	6414	Msp20I	4635	Msp20I	6414	Mva1269I	5047	Mva1269I	6131	NruI	4204	NruI	5672
NsiI	2252	NsiI	6170	PaeR7I	2241	PaeR7I	6160	PciSI	3579	PciSI	6919	PctI	5047	PctI	6131
Pfi23II	899	Pfi23II	4316	PfiMI	1455	PfiMI	3635	PmaCI	3632	PmaCI	3764	PmII	3632	PmII	3764
PspCI	3632	PspCI	3764	PspLI	899	PspLI	4316	PstNI	1770	PstNI	7451	RruI	4204	RruI	5672
Rsr2I	1203	Rsr2I	4376	RsrII	1203	RsrII	4376	SapI	3579	SapI	6919	Scal	5721	Scal	8408
Sfr274I	2241	Sfr274I	6160	Slal	2241	Slal	6160	Smal	1189	Smal	1253	StrI	2241	StrI	6160
TspGWI8368	18368	TspGWI8368	18368	TspMI	1187	TspMI	1251	Van91I	1455	Van91I	3635	VneI	7349	VneI	8595

XbaI	2191 XbaI	2604 XhoI	2241 XhoI	6160 XmaI	1187 XmaI	1251 Zrml	5721 Zrml	8408
Zsp2I	2252 Zsp2I	6170						

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

AatII	2418 AatII	4306 AatII	4804 AclI	3097 AclI	8154 AclI	8527 Adel	355 Adel	1455
Adel	4831 Apal	1648 Apal	4038 Apal	5331 Asp700I	2151 Asp700I	3687 Asp700I	8527 BbsI	2841
BbsI	4212 BbsI	5354 Bcgl	4701 Bcgl	6151 Bcgl	8433 Blpl	1836 Blpl	1898 Blpl	3624
Bpil	2841 Bpil	4212 Bpil	5354 Bpu1102I	1836 Bpu1102I	1898 Bpu1102I	3624 BpuAI	2841 BpuAI	4212
BpuAI	5354 BpvUI	628 BpvUI	2903 BpvUI	8298 BseRI	1735 BseRI	4077 BseRI	4257 Bsp120I	1644
Bsp120I	4034 Bsp120I	5327 Bsp1720I	1836 Bsp1720I	1898 Bsp1720I	3624 BspHI	48 BspHI	5877 BspHI	7755
BstENI	2993 BstENI	3184 BstENI	5322 BstV2I	2841 BstV2I	4212 BstV2I	5354 CciI	48 CciI	5877
CciI	7755 CciNI	1565 CciNI	2235 CciNI	2884 CellI	1836 CellI	1898 CellI	3624 Csil	4827
Csil	5230 Csil	5386 CspCI	149 CspCI	2347 CspCI	3222 DralII	355 DralII	1455 DralII	4831
Eco147I	1346 Eco147I	1355 Eco147I	4591 EcoNI	2993 EcoNI	3184 EcoNI	5322 HincII	1406 HincII	4966
HincII	5228 HindII	1406 HindII	4966 HindII	5228 Mabl	4827 Mabl	5230 Mabl	5386 MroXI	2151
MroXI	3687 MroXI	8527 Mvrl	628 Mvrl	2903 Mvrl	8298 NotI	1565 NotI	2235 NotI	2884
PaeI	2250 PaeI	2852 PaeI	6172 PagI	48 PagI	5877 PagI	7755 PasI	1138 PasI	2134
PasI	4309 Pcel	1346 Pcel	1355 Pcel	4591 PdmI	2151 PdmI	3687 PdmI	8527 Ple19I	628
Ple19I	2903 Ple19I	8298 Psp1406I	3097 Psp1406I	8154 Psp1406I	8527 PspOMI	1644 PspOMI	4034 PspOMI	5327
PvuI	628 PvuI	2903 PvuI	8298 PvuII	657 PvuII	3134 PvuII	6859 Rcal	48 Rcal	5877
Rcal	7755 SexAI	4827 SexAI	5230 SexAI	5386 SphI	2250 SphI	2852 SphI	6172 SseBI	1346
SseBI	1355 SseBI	4591 Sspl	17 Sspl	147 Sspl	5177 Stul	1346 Stul	1355 Stul	4591
XagI	2993 XagI	3184 XagI	5322 XmnI	2151 XmnI	3687 XmnI	8527 Zral	2416 Zral	4304
Zral	4802							