

DC-DON-SH02 Datasheet

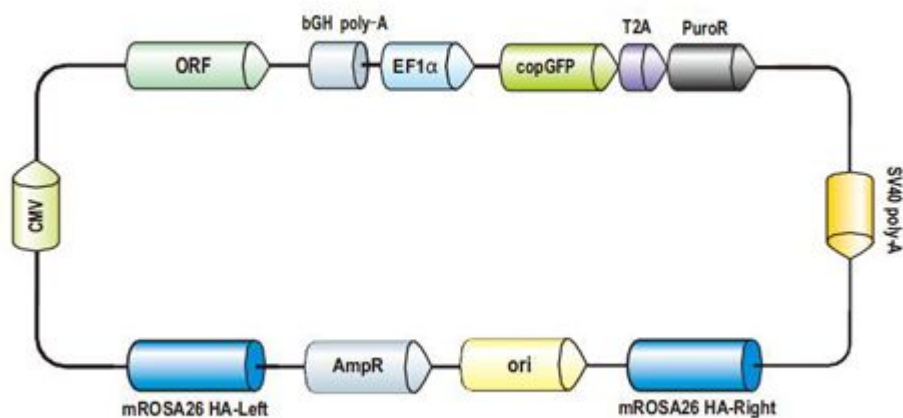
Clone Information

Catalog No.: DC-DON-SH02

Description: ROSA26 donor vector

Whole Plasmid Size: 10366 bp

Resistance: Ampicillin



Restriction Enzyme Information for DC-DON-SH02

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

Acc65I AfeI AflII AgeI AscI AvrII BaeI BaeI
 BclI ClaI KpnI NdeI NheI PacI PmeI PpuMI
 PshAI SanDI SgrAI SmaI

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

AarI	5552 AhdI	9579 AhII	1359 AIOI	289 AsiSI	4554 AspEI	9579 Aspl	5953 AsuII	2574
AxyI	4836 BbvCI	3752 BcuI	1359 BglII	4528 BmeRI	9579 Bpu14I	2574 Bse21I	4836 Bsp119I	2574
BspT104I	2574 BssNAI	3562 Bst1107I	3562 BstAPI	6482 BstBI	2574 BstEII	6045 BstPI	6045 BstSNI	2274
BstZ17I	3562 Bsu36I	4836 Cfr42I	6125 Dril	9579 Eam1105I	9579 Ecl136II	2500 Eco105I	2274 Eco32I	2183
Eco53kI	2500 Eco81I	4836 Eco91I	6045 EcoICRI	2500 EcoO65I	6045 EcoRV	2183 FseI	5444 KspI	6125
MauBI	1519 MfeI	4626 MluI	2159 MunI	4626 NspV	2574 PciI	8686 PfiFI	5953 PfoI	6386
PscI	8686 Psp124BI	2502 PspEI	6045 PspXI	2204 PstI	5953 RgaI	4554 RgiI	5444 SacI	2502
SacII	6125 SalI	4155 SbfI	7062 SdaI	7062 SfaI	4554 SfiI	1352 Sfr303I	6125 SfuI	2574
SgfI	4554 SgrBI	6125 SnaBI	2274 SpeI	1359 Sse8387I	7062 SstI	2502 Tth111I	5953 XcmI	5814

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

AasI	310 AasI	8794 AccI	3561 AccI	4156 AcvI	5283 AcvI	5415 AflIII	2159 AflIII	8686
ArsI	3294 ArsI	4931 BbrPI	5283 BbrPI	5415 BpII	736 BpII	768 BsiWI	899 BsiWI	5967
Bsp68I	5855 Bsp68I	7323 BspQI	5230 BspQI	8570 BtuMI	5855 BtuMI	7323 CpoI	1203 CpoI	6027
CspI	1203 CspI	6027 DrdI	310 DrdI	8794 DseDI	310 DseDI	8794 Eco72I	5283 Eco72I	5415
EcoRI	2171 EcoRI	3027 EcoT22I	2215 EcoT22I	7821 Fall	7745 Fall	7777 FbII	3561 FbII	4156
HindIII	4789 HindIII	7754 HpaI	1369 HpaI	6617 KspAI	1369 KspAI	6617 LguI	5230 LguI	8570
Mph1103I	2215 Mph1103I	7821 NruI	5855 NruI	7323 NsiI	2215 NsiI	7821 PaeR7I	2204 PaeR7I	7811
PciSI	5230 PciSI	8570 Pfi23II	899 Pfi23II	5967 PmaCI	5283 PmaCI	5415 PmlI	5283 PmlI	5415
PspCI	5283 PspCI	5415 PspLI	899 PspLI	5967 RruI	5855 RruI	7323 Rsr2I	1203 Rsr2I	6027
RsrII	1203 RsrII	6027 SapI	5230 SapI	8570 Sfr274I	2204 Sfr274I	7811 Slal	2204 Slal	7811
StrI	2204 StrI	7811 XbaI	2154 XbaI	2567 XhoI	2204 XhoI	7811 XmiI	3561 XmiI	4156
Zsp2I	2215 Zsp2I	7821						

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

AatII	2381	AatII	5957	AatII	6455	Acc36I	4142	Acc36I	5279	Acc36I	5552	AccB7I	1418	AccB7I	3259
AccB7I	5286	Adel	355	Adel	1418	Adel	6482	Ajil	1455	Ajil	3932	Ajil	5889	Alw44I	3927
Alw44I	9000	Alw44I	10246	AlwNI	1733	AlwNI	3806	AlwNI	9102	Apal	1611	Apal	5689	Apal	6982
Apal	3927	Apal	9000	Apal	10246	Asp700I	2114	Asp700I	5338	Asp700I	10178	AssI	3444	AssI	7372
AssI	10059	BamHI	2777	BamHI	3480	BamHI	4297	BasI	1418	BasI	3259	BasI	5286	BbsI	4492
BbsI	5863	BbsI	7005	Bcgl	6352	Bcgl	7802	Bcgl	10084	BfuAI	4142	BfuAI	5279	BfuAI	5552
BlpI	1799	BlpI	1861	BlpI	5275	BmcAI	3444	BmcAI	7372	BmcAI	10059	BmgBI	1455	BmgBI	3932
BmgBI	5889	Bpil	4492	Bpil	5863	Bpil	7005	Bpu10I	2800	Bpu10I	3752	Bpu10I	4930	Bpu1102I	1799
Bpu1102I	1861	Bpu1102I	5275	BpuAI	4492	BpuAI	5863	BpuAI	7005	BpvUI	628	BpvUI	4554	BpvUI	9949
BsaBI	2185	BsaBI	3911	BsaBI	4291	Bse3DI	3046	Bse3DI	9640	Bse3DI	9814	Bse8I	2185	Bse8I	3911
Bse8I	4291	BseJI	2185	BseJI	3911	BseJI	4291	BseMI	3046	BseMI	9640	BseMI	9814	BseRI	1698
BseRI	5728	BseRI	5908	BsmBI	2523	BsmBI	3251	BsmBI	6303	Bsp120I	1607	Bsp120I	5685	Bsp120I	6978
Bsp1720I	1799	Bsp1720I	1861	Bsp1720I	5275	BspHI	48	BspHI	7528	BspHI	9406	BspMI	4142	BspMI	5279
BspMI	5552	BsrDI	3046	BsrDI	9640	BsrDI	9814	BstENI	4644	BstENI	4835	BstENI	6973	BstV2I	4492
BstV2I	5863	BstV2I	7005	Btrl	1455	Btrl	3932	Btrl	5889	Bvel	4142	Bvel	5279	Bvel	5552
Cail	1733	Cail	3806	Cail	9102	Ccil	48	Ccil	7528	Ccil	9406	CellI	1799	CellI	1861
CellI	5275	Cfr9I	1187	Cfr9I	1251	Cfr9I	3896	Csil	6478	Csil	6881	Csil	7037	CspCI	149
CspCI	2310	CspCI	4873	DrallI	355	DrallI	1418	DrallI	6482	Eco147I	1346	Eco147I	1355	Eco147I	6242
EcoNI	4644	EcoNI	4835	EcoNI	6973	Esp3I	2523	Esp3I	3251	Esp3I	6303	MabI	6478	MabI	6881
MabI	7037	MroXI	2114	MroXI	5338	MroXI	10178	Mvrl	628	Mvrl	4554	Mvrl	9949	Pael	2213
Pael	4503	Pael	7823	Pagl	48	Pagl	7528	Pagl	9406	Pcel	1346	Pcel	1355	Pcel	6242
Pdml	2114	Pdml	5338	Pdml	10178	PflMI	1418	PflMI	3259	PflMI	5286	Ple19I	628	Ple19I	4554
Ple19I	9949	PspOMI	1607	PspOMI	5685	PspOMI	6978	PstNI	1733	PstNI	3806	PstNI	9102	PvuI	628
PvuI	4554	PvuI	9949	Rcal	48	Rcal	7528	Rcal	9406	Scal	3444	Scal	7372	Scal	10059
SexAI	6478	SexAI	6881	SexAI	7037	Smal	1189	Smal	1253	Smal	3898	SphI	2213	SphI	4503
SphI	7823	SseBI	1346	SseBI	1355	SseBI	6242	Stul	1346	Stul	1355	Stul	6242	TspMI	1187
TspMI	1251	TspMI	3896	Van91I	1418	Van91I	3259	Van91I	5286	Vnel	3927	Vnel	9000	Vnel	10246
Xagl	4644	Xagl	4835	Xagl	6973	Xmal	1187	Xmal	1251	Xmal	3896	Xmnl	2114	Xmnl	5338
Xmnl	10178	Zral	2379	Zral	5955	Zral	6453	Zrml	3444	Zrml	7372	Zrml	10059		
