

DC-RFP-SH01 Datasheet

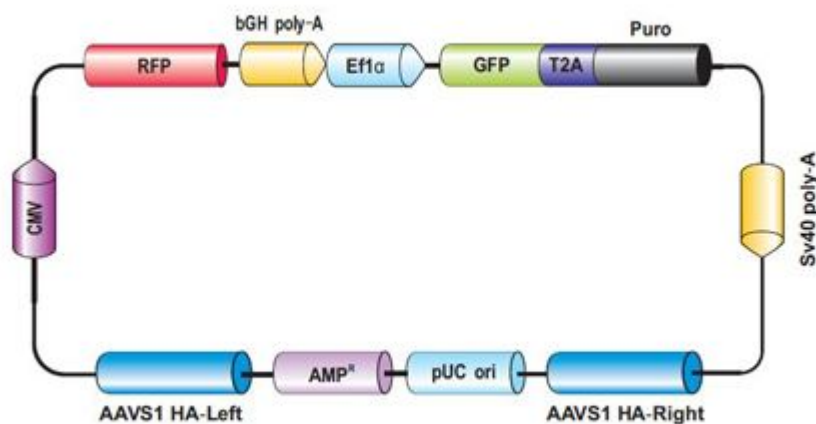
Clone Information

Catalog No.: DC-RFP-SH01

Description: AAVS1 RFP control

Whole Plasmid Size: 9430 bp

Resistance: Ampicillin



Restriction Enzyme Information for DC-RFP-SH01

Table 1. Restriction Enzymes That Do Not Cleave DC-RFP-SH01

AccI Afel AfII Agel AscI BaeI BaeI BbvCI
 BclI BstZ17I ClaI PacI PmeI SalI SbfI SgrAI

Table 2. Restriction Enzymes That Cleave DC-RFP-SH01 Once

AarI	4588	Acc65I	1396	AhII	1359	Ajil	4925	Ajul	1376	AsiSI	3590	Asp718I	1396	AspA2I	6607
AspI	4989	AssI	9123	AsuII	3320	AsuNHI	1984	AvrII	6607	BcuI	1359	BgIII	3564	BlnI	6607
BmcAI	9123	BmgBI	4925	BmtI	1988	BoxI	1771	Bpu14I	3320	BsaBI	2232	Bse8I	2232	BseJI	2232
Bsp119I	3320	Bsp1407I	6030	Bsp68I	4891	BspOI	1988	BspT104I	3320	BsrGI	6030	BstAPI	5518	BstAUI	6030
BstBI	3320	BstPAI	1771	BstSNI	2321	BtrI	4925	BtuMI	4891	Cfr42I	5161	Eco105I	2321	FauNDI	2669
FseI	4480	HindIII	3825	HpaI	5653	KfiI	6718	KpnI	1400	KspAI	5653	KspI	5161	MfeI	3662
MluI	2206	MunI	3662	NdeI	2669	NheI	1984	NruI	4891	NspV	3320	PciI	7750	PfiFI	4989
PscI	7750	PshAI	1771	PspXI	2251	PsrI	2962	PsyI	4989	RgaI	3590	RigI	4480	RruI	4891
SacII	5161	Scal	9123	SfaAI	3590	SfiI	1352	Sfr303I	5161	SfuI	3320	SgfI	3590	SgrBI	5161
SmiI	3327	SnaBI	2321	SpeI	1359	Swal	3327	Tth111I	4989	XbaI	2614	XmaJI	6607	Zrml	9123

Table 3. Restriction Enzymes That Cleave DC-RFP-SH01 Twice

AanI	227	AanI	5673	Acc16I	607	Acc16I	8865	AflIII	2206	AflIII	7750	AhdI	2799	AhdI	8643
AlfI	3565	AlfI	3599	Alol	289	Alol	1921	Alw44I	8064	Alw44I	9310	AlwNI	6324	AlwNI	8166
ApaLI	8064	ApaLI	9310	ArsI	2070	ArsI	3967	AspEI	2799	AspEI	8643	AvII	607	AvII	8865
BauI	7923	BauI	9307	BlpI	2032	BlpI	4311	BmeRI	2799	BmeRI	8643	Bpu1102I	2032	Bpu1102I	4311
BsiWI	899	BsiWI	5003	Bsml	1417	Bsml	5734	Bsp1720I	2032	Bsp1720I	4311	BspHI	48	BspHI	8470
BssSI	7923	BssSI	9307	Bst2BI	7923	Bst2BI	9307	Cail	6324	Cail	8166	Ccil	48	Ccil	8470
CciNI	2245	CciNI	3571	CelII	2032	CelII	4311	Cpol	1203	Cpol	5063	Cspl	1203	Cspl	5063
Dril	2799	Dril	8643	Eam1105I	2799	Eam1105I	8643	Ecl136II	2031	Ecl136II	2547	Eco32I	2230	Eco32I	2949
Eco53kI	2031	Eco53kI	2547	EcoICRI	2031	EcoICRI	2547	EcoRI	1384	EcoRI	2218	EcoRV	2230	EcoRV	2949
EcoT22I	2262	EcoT22I	6885	Fall	6188	Fall	6220	FspI	607	FspI	8865	HincII	5653	HincII	5915
HindII	5653	HindII	5915	Mph1103I	2262	Mph1103I	6885	Mva1269I	1417	Mva1269I	5734	NotI	2245	NotI	3571

Nsbl	607	Nsbl	8865	Nsil	2262	Nsil	6885	PaeR7I	2251	PaeR7I	6875	PagI	48	PagI	8470
PctI	1417	PctI	5734	Pfi23II	899	Pfi23II	5003	PfoI	5422	PfoI	6163	PpuMI	1646	PpuMI	6718
Psil	227	Psil	5673	Psp124BI	2033	Psp124BI	2549	Psp5II	1646	Psp5II	6718	PspLI	899	PspLI	5003
PspPPI	1646	PspPPI	6718	PstNI	6324	PstNI	8166	RcaI	48	RcaI	8470	Rsr2I	1203	Rsr2I	5063
RsrII	1203	RsrII	5063	SacI	2033	SacI	2549	Sfr274I	2251	Sfr274I	6875	Slal	2251	Slal	6875
SstI	2033	SstI	2549	StrI	2251	StrI	6875	VneI	8064	VneI	9310	XcmI	4850	XcmI	6538
XhoI	2251	XhoI	6875	Zsp2I	2262	Zsp2I	6885								

Table 4. Restriction Enzymes That Cleave DC-RFP-SH01 Three Times

AasI	310	AasI	6343	AasI	7858	AatII	2428	AatII	4993	AatII	5491	AccB7I	4322	AccB7I	6164
AccB7I	6625	AccI	3784	AccI	8869	AccI	9242	AcvI	1611	AcvI	4319	AcvI	4451	ApaI	2062
ApaI	4725	ApaI	6018	Asp700I	2974	Asp700I	4374	Asp700I	9242	AxyI	3872	AxyI	6073	AxyI	6730
Ball	5322	Ball	6297	Ball	7129	BasI	4322	BasI	6164	BasI	6625	BbrPI	1611	BbrPI	4319
BbrPI	4451	Bpu10I	3966	Bpu10I	6325	Bpu10I	6786	BpvUI	628	BpvUI	3590	BpvUI	9013	Bse21I	3872
Bse21I	6073	Bse21I	6730	Bse3DI	3122	Bse3DI	8704	Bse3DI	8878	BseMI	3122	BseMI	8704	BseMI	8878
BsePI	747	BsePI	5402	BsePI	7389	BsmBI	1855	BsmBI	2570	BsmBI	5339	Bsp120I	2058	Bsp120I	4721
Bsp120I	6014	BsrDI	3122	BsrDI	8704	BsrDI	8878	BssHII	747	BssHII	5402	BssHII	7389	BstEII	1468
BstEII	1923	BstEII	5081	BstPI	1468	BstPI	1923	BstPI	5081	BstXI	1380	BstXI	2240	BstXI	6025
Bsu36I	3872	Bsu36I	6073	Bsu36I	6730	Cfr9I	1187	Cfr9I	1251	Cfr9I	2055	Csil	5514	Csil	5917
Csil	6618	CspCI	149	CspCI	2357	CspCI	3909	DrdI	310	DrdI	6343	DrdI	7858	DseDI	310
DseDI	6343	DseDI	7858	Eco72I	1611	Eco72I	4319	Eco72I	4451	Eco81I	3872	Eco81I	6073	Eco81I	6730
Eco91I	1468	Eco91I	1923	Eco91I	5081	EcoO65I	1468	EcoO65I	1923	EcoO65I	5081	Esp3I	1855	Esp3I	2570
Esp3I	5339	MabI	5514	MabI	5917	MabI	6618	MisI	5322	MisI	6297	MisI	7129	MluNI	5322
MluNI	6297	MluNI	7129	MroXI	2974	MroXI	4374	MroXI	9242	MscI	5322	MscI	6297	MscI	7129
Msp20I	5322	Msp20I	6297	Msp20I	7129	MvrI	628	MvrI	3590	MvrI	9013	PaeI	2260	PaeI	3539
PaeI	6887	Paul	747	Paul	5402	Paul	7389	Pdml	2974	Pdml	4374	Pdml	9242	PfiMI	4322
PfiMI	6164	PfiMI	6625	Ple19I	628	Ple19I	3590	Ple19I	9013	PmaCI	1611	PmaCI	4319	PmaCI	4451
PmlI	1611	PmlI	4319	PmlI	4451	Psp1406I	3784	Psp1406I	8869	Psp1406I	9242	PspCI	1611	PspCI	4319
PspCI	4451	PspEI	1468	PspEI	1923	PspEI	5081	PspOMI	2058	PspOMI	4721	PspOMI	6014	Ptel	747
Ptel	5402	Ptel	7389	PvuI	628	PvuI	3590	PvuI	9013	SexAI	5514	SexAI	5917	SexAI	6618
SmaI	1189	SmaI	1253	SmaI	2057	SphI	2260	SphI	3539	SphI	6887	Sspl	17	Sspl	147
Sspl	5864	TspMI	1187	TspMI	1251	TspMI	2055	Van91I	4322	Van91I	6164	Van91I	6625	XmaI	1187
XmaI	1251	XmaI	2055	Xmnl	2974	Xmnl	4374	Xmnl	9242	Zral	2426	Zral	4991	Zral	5489

