

DC-DON-SH01 Datasheet

Vector Information

Catalog No.: DC-DON-SH01

Description: AAVS1 donor vector

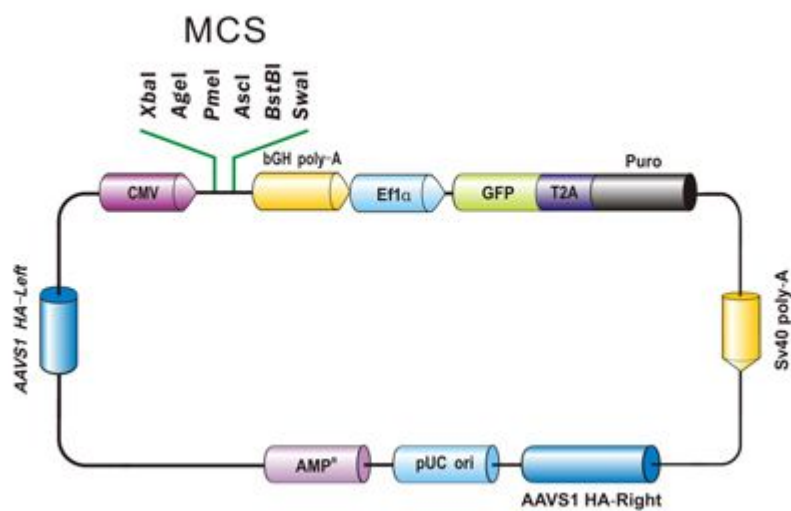
Whole Plasmid Size: 8753 bp

Resistance: Ampicillin

Suggested Sequencing Primers:

Forward: 5'-CGGTGGGAGGTCTATATAAGCAG-3'

Reverse: 5'-GACAGTGGGAGTGGCACCTTC-3'



Restriction Enzyme Information for DC-DON-SH01

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

AccIAfeI AfIII BaeI BaeI BbvCI BclI BstZ17I
 ClaI NdeI PacI SalI SbfI SgrAI

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

AarI	3911	Acc65I	1396	AgeI	2620	AhdI	7966	AhlI	1359	AjiI	4248	AjuI	1376	AscI	2635
AsiGI	2620	AsiSI	2913	Asp718I	1396	AspA2I	5930	AspEI	7966	AspI	4312	AssI	8446	AsuII	2643
AsuNHI	1984	AvrII	5930	BcuI	1359	BglII	2887	BlnI	5930	BmcAI	8446	BmeRI	7966	BmgBI	4248
BmtI	1988	BoxI	1771	Bpu14I	2643	BsaBI	2232	Bse8I	2232	BseJI	2232	BshTI	2620	Bsp119I	2643
Bsp1407I	5353	Bsp68I	4214	BspOI	1988	BspT104I	2643	BsrGI	5353	BstAPI	4841	BstAUI	5353	BstBI	2643
BstPAI	1771	BstSNI	2321	BtrI	4248	BtuMI	4214	Cfr42I	4484	CspAI	2620	DrI	7966	Eam1105I	7966
Eco105I	2321	Eco32I	2230	EcoRV	2230	FseI	3803	HindIII	3148	HpaI	4976	KfiI	6041	KpnI	1400
KspAI	4976	KspI	4484	MfeI	2985	MluI	2206	MssI	2629	MunI	2985	NheI	1984	NruI	4214
NspV	2643	PalAI	2635	PciI	7073	PfiFI	4312	PinAI	2620	PmeI	2629	PscI	7073	PshAI	1771
PspXI	2251	PsyI	4312	RgaI	2913	RigI	3803	RruI	4214	SacII	4484	Scal	8446	SfaAI	2913
SfiI	1352	Sfr303I	4484	SfuI	2643	SgfI	2913	SgrBI	4484	SgsI	2635	SmI	2650	SnaBI	2321
SpeI	1359	Swal	2650	Tth111I	4312	XbaI	2614	XmaJI	5930	Zrml	8446				

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

AanI	227	AanI	4996	Acc16I	607	Acc16I	8188	AfIII	2206	AfIII	7073	AlfI	2888	AlfI	2922
Alol	289	Alol	1921	Alw44I	7387	Alw44I	8633	AlwNI	5647	AlwNI	7489	ApaLI	7387	ApaLI	8633
ArsI	2070	ArsI	3290	Asp700I	3697	Asp700I	8565	AvIII	607	AvIII	8188	BauI	7246	BauI	8630
BlpI	2032	BlpI	3634	Bpu1102I	2032	Bpu1102I	3634	Bse3DI	8027	Bse3DI	8201	BseMI	8027	BseMI	8201
BsiWI	899	BsiWI	4326	Bsml	1417	Bsml	5057	Bsp1720I	2032	Bsp1720I	3634	BspHI	48	BspHI	7793
BsrDI	8027	BsrDI	8201	BssSI	7246	BssSI	8630	Bst2BI	7246	Bst2BI	8630	CaiI	5647	CaiI	7489

CciI	48	CciI	7793 CciNI	2245 CciNI	2894 CeliI	2032 CeliI	3634 Cpol	1203 Cpol	4386
Cspl	1203 Cspl	4386 Ecl136II	2031 Ecl136II	2547 Eco53kl	2031 Eco53kl	2547 EcoICRI	2031 EcoICRI	2547	
EcoRI	1384 EcoRI	2218 EcoT22I	2262 EcoT22I	6208 Fall	5511 Fall	5543 Fspl	607 Fspl	8188	
HincII	4976 HincII	5238 HindII	4976 HindII	5238 Mph1103I	2262 Mph1103I	6208 MroXI	3697 MroXI	8565	
Mva1269I	1417 Mva1269I	5057 NotI	2245 NotI	2894 Nsbl	607 Nsbl	8188 Nsil	2262 Nsil	6208	
PaeR7I	2251 PaeR7I	6198 PagI	48 PagI	7793 Pctl	1417 Pctl	5057 Pdml	3697 Pdml	8565	
Pfl23II	899 Pfl23II	4326 PfoI	4745 PfoI	5486 PpuMI	1646 PpuMI	6041 Psil	227 Psil	4996	
Psp124BI	2033 Psp124BI	2549 Psp5II	1646 Psp5II	6041 PspLI	899 PspLI	4326 PspPPI	1646 PspPPI	6041	
PstNI	5647 PstNI	7489 Rcal	48 Rcal	7793 Rsr2I	1203 Rsr2I	4386 RsrII	1203 RsrII	4386	
SacI	2033 SacI	2549 Sfr274I	2251 Sfr274I	6198 Slal	2251 Slal	6198 SstI	2033 SstI	2549	
StrI	2251 StrI	6198 Vnel	7387 Vnel	8633 Xcml	4173 Xcml	5861 Xhol	2251 Xhol	6198	
XmnI	3697 XmnI	8565 Zsp2I	2262 Zsp2I	6208					

AasI	310 AasI	5666 AasI	7181 AatII	2428 AatII	4316 AatII	4814 AccB7I	3645 AccB7I	5487
AccB7I	5948 AcII	3107 AcII	8192 AcII	8565 AcvI	1611 AcvI	3642 AcvI	3774 Adel	355
Adel	1611 Adel	4841 Apal	2062 Apal	4048 Apal	5341 AxyI	3195 AxyI	5396 AxyI	6053
Ball	4645 Ball	5620 Ball	6452 BasI	3645 BasI	5487 BasI	5948 BbrPI	1611 BbrPI	3642
BbrPI	3774 Bcgl	4711 Bcgl	6189 Bcgl	8471 Bpu10I	3289 Bpu10I	5648 Bpu10I	6109 BpvUI	628
BpvUI	2913 BpvUI	8336 Bse21I	3195 Bse21I	5396 Bse21I	6053 BsmBI	1855 BsmBI	2570 BsmBI	4662
Bsp120I	2058 Bsp120I	4044 Bsp120I	5337 BstEI	1468 BstEI	1923 BstEI	4404 BstPI	1468 BstPI	1923
BstPI	4404 BstXI	1380 BstXI	2240 BstXI	5348 Bsu36I	3195 Bsu36I	5396 Bsu36I	6053 Cfr9I	1187
Cfr9I	1251 Cfr9I	2055 CsII	4837 CsII	5240 CsII	5941 CspCI	149 CspCI	2357 CspCI	3232
DraIII	355 DraIII	1611 DraIII	4841 DrdI	310 DrdI	5666 DrdI	7181 DseDI	310 DseDI	5666
DseDI	7181 Eco72I	1611 Eco72I	3642 Eco72I	3774 Eco81I	3195 Eco81I	5396 Eco81I	6053 Eco91I	1468
Eco91I	1923 Eco91I	4404 EcoO65I	1468 EcoO65I	1923 EcoO65I	4404 Esp3I	1855 Esp3I	2570 Esp3I	4662
MabI	4837 MabI	5240 MabI	5941 MIsI	4645 MIsI	5620 MIsI	6452 MluNI	4645 MluNI	5620
MluNI	6452 MscI	4645 MscI	5620 MscI	6452 Msp20I	4645 Msp20I	5620 Msp20I	6452 MvrI	628
MvrI	2913 MvrI	8336 PaeI	2260 PaeI	2862 PaeI	6210 PfIM	3645 PfIM	5487 PfIM	5948
Ple19I	628 Ple19I	2913 Ple19I	8336 PmaCI	1611 PmaCI	3642 PmaCI	3774 PmII	1611 PmII	3642
PmII	3774 Psp1406I	3107 Psp1406I	8192 Psp1406I	8565 PspCI	1611 PspCI	3642 PspCI	3774 PspEI	1468
PspEI	1923 PspEI	4404 PspOMI	2058 PspOMI	4044 PspOMI	5337 Pvul	628 Pvul	2913 Pvul	8336
PvuII	657 Pvull	3144 Pvull	6897 SexAI	4837 SexAI	5240 SexAI	5941 Smal	1189 Smal	1253
Smal	2057 SphI	2260 SphI	2862 SphI	6210 Sspl	17 Sspl	147 Sspl	5187 TspMI	1187
TspMI	1251 TspMI	2055 Van91I	3645 Van91I	5487 Van91I	5948 XmaI	1187 XmaI	1251 XmaI	2055
Zral	2426 Zral	4314 Zral	4812					
