

HCP-AAVS1-CG02 Datasheet

Clone Information

Catalog No.: HCP-AAVS1-CG02

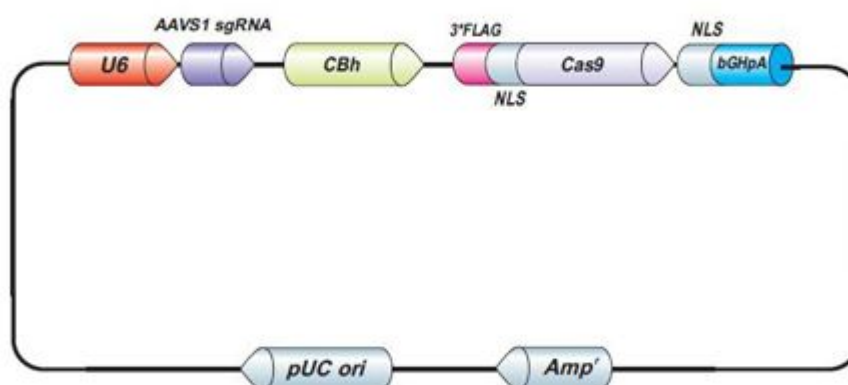
Description: All-in-one sgRNA/Cas9 expression clone for human AAVS1 safe harbor site

Target site: GGGGCCACTAGGGACAGGAT

Whole Plasmid Size: 8508 bp

Resistance: Ampicillin

Vector Information



Restriction Enzyme Information for HCP-AAVS1-CG02

Table 1. Restriction Enzymes That Do Not Cleave HCP-AAVS1-CG02

AccI	AfeI	AflII	AscI	AvrII	BaeI	BaeI	BamHI
BbsI	BbvCI	BpII	BsaI	BsiWI	BsrGI	BstBI	BstEII
BstZ17I	Bsu36I	Clal	HincII	HindIII	HpaI	MfeI	MluI
NheI	NruI	NsiI	PacI	PmeI	Ppu10I	PshAI	RsrII
Sall	SapI	SexAI	SfiI	SgfI	SgrAI	SpeI	StuI
Swal	TliI	Tth111I	XhoI				

Table 2. Restriction Enzymes That Cleave HCP-AAVS1-CG02 Once

AanI	5994	AarI	869	Acc16I	7042	Acc65I	86	AcvI	3809	AflIII	8153	AgeI	891	ApaI	2379
ArsI	659	AsiGI	891	Asp718I	86	AviII	7042	BbrPI	3809	BglII	1270	BpvUI	6896	BsaBI	4518
Bse8I	4518	BseJI	4518	BseRI	764	BshTI	891	BsmI	4660	Bsp120I	2375	BstSNI	353	CciNI	5414
Cfr42I	7200	CspAI	891	CspCI	3790	DinI	5564	Eco105I	353	Eco32I	2873	Eco72I	3809	EcoRI	5175
EcoRV	2873	EgeI	5564	EheI	5564	FseI	5152	FspI	7042	KasI	5562	KfiI	4397	KpnI	90
KspI	7200	Mly113I	5563	Mva1269I	4660	MvrI	6896	NarI	5563	NotI	5414	NsbI	7042	PasI	2789
PciI	8153	PctI	4660	PinAI	891	Ple19I	6896	PluTI	5566	PmaCI	3809	PmlI	3809	PscI	8153
PsiI	5994	PspCI	3809	PspOMI	2375	PvuI	6896	RigI	5152	SacII	7200	SbfI	5559	SdaI	5559
SfoI	5564	Sfr303I	7200	SgrBI	7200	SnaBI	353	Sse8387I	5559	SspDI	5562	XbaI	80		

Table 3. Restriction Enzymes That Cleave HCP-AAVS1-CG02 Twice

AccB7I	1319	AccB7I	3500	AccIII	3916	AccIII	4189	Adel	3419	Adel	5869	AjiI	1559	AjiI	4070
AjuI	3110	AjuI	4145	AlfI	5536	AlfI	5570	Aor13HI	3916	Aor13HI	4189	BarI	2585	BarI	4433
BasI	1319	BasI	3500	Bcgl	3171	Bcgl	6761	BmgBI	1559	BmgBI	4070	Bpml	874	Bpml	7196
BseAI	3916	BseAI	4189	BseX3I	4643	BseX3I	5414	BsmBI	6228	BsmBI	6270	Bsp13I	3916	Bsp13I	4189
Bsp19I	373	Bsp19I	901	BspEI	3916	BspEI	4189	BstAPI	2045	BstAPI	2555	BstENI	1527	BstENI	1794
BstXI	1064	BstXI	1634	BstZI	4643	BstZI	5414	Btrl	1559	Btrl	4070	BtsI	6846	BtsI	6866
Cfr9I	5506	Cfr9I	5517	DrallI	3419	DrallI	5869	EagI	4643	EagI	5414	Ecl136II	3257	Ecl136II	5186

EclXI	4643 EclXI	5414 Eco52I	4643 Eco52I	5414 Eco53kl	3257 Eco53kl	5186 EcoICRI	3257 EcoICRI	5186
EcoNI	1527 EcoNI	1794 Esp3I	6228 Esp3I	6270 FauNDI	247 FauNDI	8337 Gsul	874 Gsul	7196
Kpn2I	3916 Kpn2I	4189 MroI	3916 MroI	4189 NcoI	373 NcoI	901 NdeI	247 NdeI	8337
PaeI	4286 PaeI	5405 PfiMI	1319 PfiMI	3500 Psp124BI	3259 Psp124BI	5188 SacI	3259 SacI	5188
SmaI	5508 SmaI	5519 SphI	4286 SphI	5405 SspI	6074 SspI	6460 SstI	3259 SstI	5188
TspMI	5506 TspMI	5517 Van91I	1319 Van91I	3500 XagI	1527 XagI	1794 XcmI	4559 XcmI	5042
XmaI	5506 XmaI	5517						

Table 4. Restriction Enzymes That Cleave HCP-AAVS1-CG02 Three Times

AasI	5913 AasI	6182 AasI	8051 AclI	6079 AclI	6663 AclI	7036 AhdI	3038 AhdI	5492
AhdI	7265 Alw44I	6096 Alw44I	6593 Alw44I	7839 ApaLI	6096 ApaLI	6593 ApaLI	7839 AseI	838
AseI	7090 AseI	8229 Asp700I	3924 Asp700I	4152 Asp700I	6665 AspEI	3038 AspEI	5492 AspEI	7265
BciVI	1011 BciVI	6428 BciVI	7955 BfuI	1011 BfuI	6428 BfuI	7955 BmeRI	3038 BmeRI	5492
BmeRI	7265 BmrI	302 BmrI	4403 BmrI	7225 BmuI	302 BmuI	4403 BmuI	7225 BsaXI	3615
BsaXI	4251 BsaXI	5935 Bse3DI	3187 Bse3DI	7031 Bse3DI	7205 BseMI	3187 BseMI	7031 BseMI	7205
BsrDI	3187 BsrDI	7031 BsrDI	7205 Bsul	1011 Bsul	6428 Bsul	7955 DrdI	5913 DrdI	6182
DrdI	8051 Dril	3038 Dril	5492 Dril	7265 DseDI	5913 DseDI	6182 DseDI	8051 Eam1105I	3038
Eam1105I	5492 Eam1105I	7265 MroXI	3924 MroXI	4152 MroXI	6665 PdmI	3924 PdmI	4152 PdmI	6665
PfeI	2401 PfeI	4547 PfeI	8176 PshBI	838 PshBI	7090 PshBI	8229 Psp1406I	6079 Psp1406I	6663
Psp1406I	7036 PsrI	3828 PsrI	4643 PsrI	4710 Tfil	2401 Tfil	4547 Tfil	8176 VneI	6096
VneI	6593 VneI	7839 Vspl	838 Vspl	7090 Vspl	8229 XmnI	3924 XmnI	4152 XmnI	6665
