

TN-AAVS1-R Datasheet

Clone Information

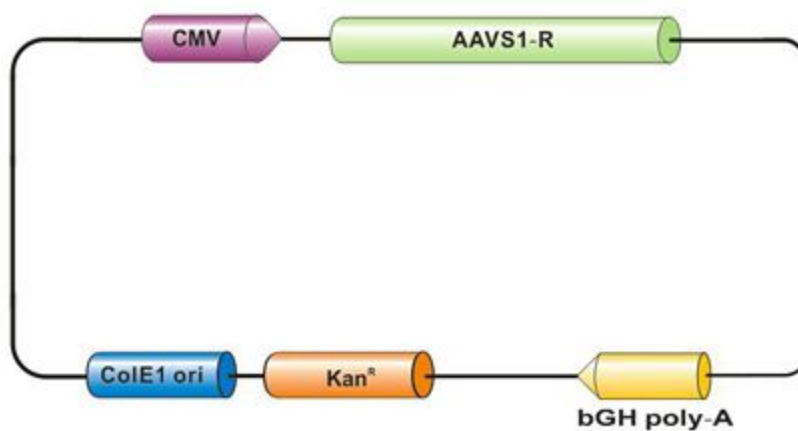
Catalog No.: TN-AAVS1-R

Description: Right TALEN clone for human AAVS1 safe harbor site

Target site: TTTCTGTCACCAATCCT

Resistance: Kanamycin

Whole plasmid size: 5973 bp



Restriction Enzyme Information for TN-AAVS1-R

Table 2. Restriction Enzymes That Cleave TN-AAVS1-R Once

PciI	64	MluI	29	PmlI	5185	RsrII	4844	NmeAIII	4552	PfiFI	4446	Tth111I	4446	EagI	4234
BbsI	3992	XbaI	3766	PaeR7I	3760	PspXI	3760	TliI	3760	XhoI	3760	BglII	3748	HindIII	3416
BseRI	3196	BamHI	3154	NdeI	285	SnaBI	391	Ec053kI	617	SacI	619	BsaI	677	NheI	696

BmtI 700 AfIII 709 BspDI 791 ClaI 791 Acc65I 855 KpnI 859 Bsu36I 917 FspAI 980
StuI 1126 BpuI 1150
