

TN-AAVS1-L Datasheet

Clone Information

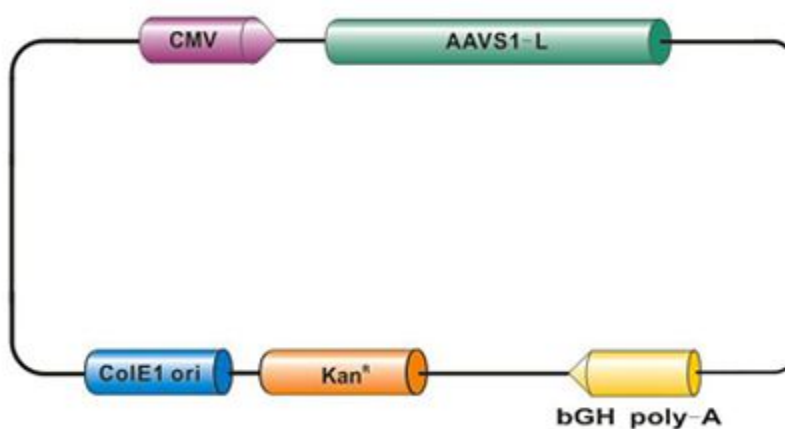
Catalog No.: TN-AAVS1-L

Description: Left TALEN clone for human AAVS1 safe harbor site

Target site: CCCCTCCACCCACAGT

Resistance: Kanamycin

Whole plasmid size: 5973 bp



Restriction Enzyme Information for TN-AAVS1-L

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

PciI	5964	PmlI	5185	RsrII	4844	NmeAIII	4552	PfiFI	4446	EagI	4234	BsaBI	4187	BbsI	3992
XbaI	3766	PaeR7I	3760	BglII	3748	HindIII	3416	BseRI	3196	BamHI	3154	MluI	29	NdeI	285
SnaBI	391	Ec053kI	617	SacI	619	BsaI	677	NheI	696	BmtI	700	AflII	709	BspDI	791

