

pDonor-D09 Datasheet

Clone Information

Catalog No.: pDonor-D09

Description: Empty donor vector

Whole Plasmid Size: 7580 bp

Vector: pDonor-D09

Antibiotic: Ampicillin

Cloning Site at 5': XbaI, AvrII, Sall, SwaI

Stable Selection Marker: Puromycin

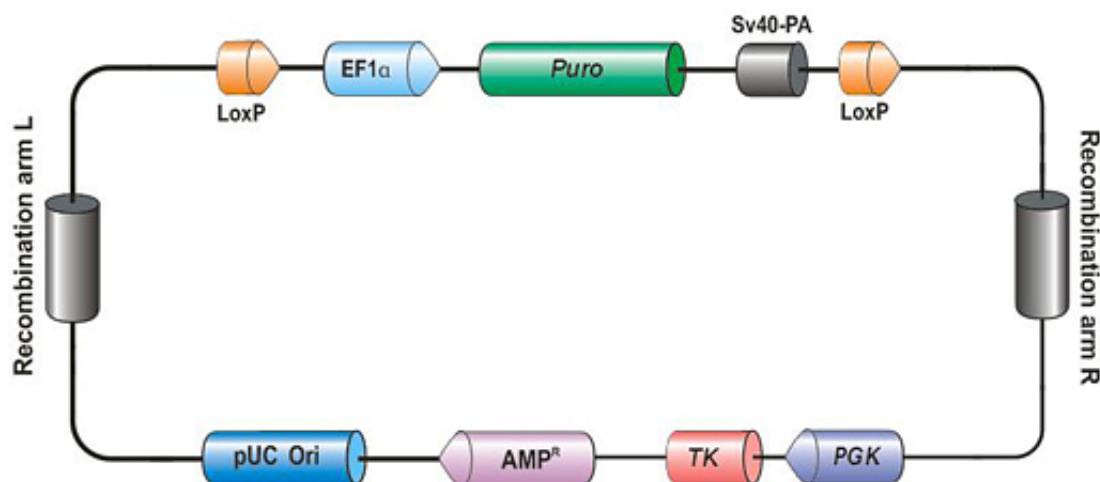
Cloning Site at 3': PacI, EcoRI, Acc65I, KpnI, NheI

Suggested Sequencing Primers:

Forward: 5'-ACAGGAAACAGCTATGACCATG-3'

Reverse: 5'-GTGGATGTGGAATGTGTGCGAG-3'

Vector Information for pDonor-D09



Restriction Enzyme Information for pDonor-D09

Table 1. Restriction Enzymes That Do Not Cleave pDonor-D09

AflII	BaeI	BaeI	BbsI	BbvCI	BpII	BsaBI	BstBI
BstZ17I	Clal	FseI	NsiI	PmeI	PmlI	Ppu10I	SfiI
SgrAI	SrfI						

Table 2. Restriction Enzymes That Cleave pDonor-D09 Once

Acc65I	7547	AfeI	4321	AgeI	2154	AhdI	6221	AhlI	2288	AjuI	4835	AleI	7517	Aor51HI	4321
ArsI	477	Ascl	64	AsiGI	2154	AsiSI	100	Asp718I	7547	AspA2I	2009	AspEI	6221	AsuNHI	7558
AvrII	2009	BamHI	1933	BarI	4540	BclI	36	BcuI	2288	BlnI	2009	BlpI	2316	BmeRI	6221
BmtI	7562	BoxI	3690	Bpu1102I	2316	BsaAI	2879	BshTI	2154	Bsp1720I	2316	Bsp68I	2931	BspOI	7562
BspQI	7226	BstBAI	2879	BstPAI	3690	BstSNI	2879	BtuMI	2931	CciNI	80	CellI	2316	Cfr9I	3745
CpoI	763	CspAI	2154	CspCI	419	CspI	763	DrI	6221	Eam1105I	6221	Eco105I	2879	Eco47III	4321
EcoRI	7539	FauNDI	5047	FbaI	36	HpaI	1356	KpnI	7551	Ksp22I	36	KspAI	1356	LguI	7226
MfeI	172	MluI	2665	MunI	172	NdeI	5047	NheI	7558	NotI	80	NruI	2931	OliI	7517
PacI	7533	PaeR7I	55	PalAI	64	PciI	7109	PciSI	7226	PinAI	2154	Ppu21I	2879	PscI	7109
PshAI	3690	RgaI	100	RruI	2931	Rsr2I	763	RsrII	763	Sall	2018	SapI	7226	SbfI	50
SdaI	50	SfaAI	100	Sfr274I	55	SgfI	100	SgsI	64	Slal	55	Smal	3747	Smil	2030
SnaBI	2879	SpeI	2288	Sse8387I	50	StrI	55	Swal	2030	TspMI	3745	XbaI	1998	XhoI	55
XmaI	3745	XmaJI	2009												

Table 3. Restriction Enzymes That Cleave pDonor-D09 Twice

AanI	1381	AanI	1994	Acc16I	4975	Acc16I	5998	AccI	2019	AccI	2875	AccIII	760	AccIII	3263
Adel	1218	Adel	4706	Aor13HI	760	Aor13HI	3263	Asp700I	4430	Asp700I	5621	AspI	689	AspI	4050
AssI	4718	AssI	5740	AvII	4975	AvII	5998	Axyl	382	Axyl	1782	BmcAI	4718	BmcAI	5740
BsaXI	1025	BsaXI	7241	Bse21I	382	Bse21I	1782	BseAI	760	BseAI	3263	BsiWI	703	BsiWI	2715
BsmI	1442	BsmI	2472	Bsp13I	760	Bsp13I	3263	Bsp1407I	1739	Bsp1407I	3550	Bsp19I	4058	Bsp19I	4144
BspEI	760	BspEI	3263	BsrGI	1739	BsrGI	3550	BstAUI	1739	BstAUI	3550	BstEII	781	BstEII	4338
BstPI	781	BstPI	4338	Bsu36I	382	Bsu36I	1782	BtgZI	117	BtgZI	3119	Csil	1214	Csil	1626
DrallI	1218	DrallI	4706	Ecl136II	3085	Ecl136II	7509	Eco53kl	3085	Eco53kl	7509	Eco81I	382	Eco81I	1782
Eco91I	781	Eco91I	4338	EcoICRI	3085	EcoICRI	7509	EcoO65I	781	EcoO65I	4338	Fall	1897	Fall	1929
FbII	2019	FbII	2875	Fspl	4975	Fspl	5998	HindIII	335	HindIII	4830	KfII	4052	KfII	4362
Kpn2I	760	Kpn2I	3263	MabI	1214	MabI	1626	MroI	760	MroI	3263	MroXI	4430	MroXI	5621
Mva1269I	1442	Mva1269I	2472	NcoI	4058	NcoI	4144	Nsbl	4975	Nsbl	5998	PaeI	2084	PaeI	3029
PctI	1442	PctI	2472	Pdml	4430	Pdml	5621	Pfl23II	703	Pfl23II	2715	PflFI	689	PflFI	4050
Psil	1381	Psil	1994	Psp124BI	3087	Psp124BI	7511	PspEI	781	PspEI	4338	PspLI	703	PspLI	2715
Psyl	689	Psyl	4050	SacI	3087	SacI	7511	Scal	4718	Scal	5740	SexAI	1214	SexAI	1626
SphI	2084	SphI	3029	Sspl	1573	Sspl	5416	SstI	3087	SstI	7511	Tth111I	1689	Tth111I	4050
Xmil	2019	Xmil	2875	Xmnl	4430	Xmnl	5621	Zrml	4718	Zrml	5740				

Table 4. Restriction Enzymes That Cleave pDonor-D09 Three Times

AasI	2871	AasI	5138	AasI	7007	Acc36I	39	Acc36I	2551	Acc36I	4385	AccB7I	1873	AccB7I	2816
AccB7I	4706	AccI	294	AccI	5619	AccI	5992	AcsI	1266	AcsI	1412	AcsI	7539	Alw44I	5052
Alw44I	5549	Alw44I	6795	ApaLI	5052	ApaLI	5549	ApalI	6795	Apol	1266	Apol	1412	Apol	7539
Asel	6046	Asel	7281	Asel	7340	Ball	1022	Ball	3291	Ball	3585	BasI	1873	BasI	2816
BasI	4706	BciVI	1316	BciVI	5384	BciVI	6911	BfuAI	39	BfuAI	2551	BfuAI	4385	BfuI	1316
BfuI	5384	BfuI	6911	BgIII	2582	BgIII	4656	BgIII	4664	Bpu10I	476	Bpu10I	2420	Bpu10I	2432
BpvUI	100	BpvUI	4956	BpvUI	5852	Bse3DI	2327	Bse3DI	5987	Bse3DI	6161	BseMI	2327	BseMI	5987
BseMI	6161	BsePI	64	BsePI	1102	BsePI	3732	BseYI	2392	BseYI	3078	BseYI	6805	BsgI	297
BsgI	2275	BsgI	2318	BspMI	39	BspMI	2551	BspMI	4385	BsrDI	2327	BsrDI	5987	BsrDI	6161
BssHII	64	BssHII	1102	BssHII	3732	BstXI	1734	BstXI	4422	BstXI	7519	Bsul	1316	Bsul	5384
Bsul	6911	Bvel	39	Bvel	2551	Bvel	4385	Cfr42I	861	Cfr42I	4360	Cfr42I	7518	DrdI	2871
DrdI	5138	DrdI	7007	DseDI	2871	DseDI	5138	DseDI	7007	Eco147I	978	Eco147I	1787	Eco147I	2337
Eco32I	2870	Eco32I	2974	Eco32I	4592	EcoRV	2870	EcoRV	2974	EcoRV	4592	Gsal	2396	Gsal	3082
Gsal	6809	KroI	531	KroI	2734	KroI	3639	Kspl	861	Kspl	4360	Kspl	7518	MisI	1022
MisI	3291	MisI	3585	MluNI	1022	MluNI	3291	MluNI	3585	Mmel	4567	Mmel	6716	Mmel	6900
MroNI	531	MroNI	2734	MroNI	3639	MscI	1022	MscI	3291	MscI	3585	Msp20I	1022	Msp20I	3291
Msp20I	3585	Mvrl	100	Mvrl	4956	Mvrl	5852	NaeI	533	NaeI	2736	NaeI	3641	NgoMIV	531
NgoMIV	2734	NgoMIV	3639	Paul	64	Paul	1102	Paul	3732	Pcel	978	Pcel	1787	Pcel	2337
Pdil	533	Pdil	2736	Pdil	3641	PflMI	1873	PflMI	2816	PflMI	4706	Ple19I	100	Ple19I	4956
Ple19I	5852	PshBI	6046	PshBI	7281	PshBI	7340	Psp1406I	294	Psp1406I	5619	Psp1406I	5992	Ptel	64
Ptel	1102	Ptel	3732	PvuI	100	PvuI	4956	PvuI	5852	PvuII	331	PvuII	4925	PvuII	7289
SacII	861	SacII	4360	SacII	7518	Sfr303I	861	Sfr303I	4360	Sfr303I	7518	SgrBI	861	SgrBI	4360
SgrBI	7518	SseBI	978	SseBI	1787	SseBI	2337	StuI	978	StuI	1787	StuI	2337	Van91I	1873

Van91I	2816	Van91I	4706	Vnel	5052	Vnel	5549	Vnel	6795	Vspl	6046	Vspl	7281	Vspl	7340
Xapl	1266	Xapl	1412	Xapl	7539	Xcml	2840	Xcml	4033	Xcml	4830				
