

pET-11a Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	1	5604	BsrFI	8	160 495 504 871 2083	NspI	4	660 3134 3426 3793		
AccI	1	3559				Pfi1108I	2	2072 4700		
AceIII	7	952 1680 2011 3298 3439	BssHII	1	1596	PfiMI	3	767 2635 2684		
		3741 4981	Bst1107I	1	3560	PleI	7	446 734 821 1617 3683		
AcI	89		BstEII	1	1366			4168 4671		
AflIII	2	1185 3789	BstXI	3	987 1116 1239	PshAI	1	2030		
AluI	24		BstYI	11		Psp5II	2	2753 2795		
AlwI	16		Cac8I	42		Psp1406I	5	847 2215 3114 4908 5281		
Alw21I	8	685 1169 2492 2783 3607	CjeI	28		PstI	1	4927		
		4107 5268 5353	CjePI	26		PvuI	1	5052		
Alw44I	4	1165 3603 4103 5349	Clal	1	24	PvuII	3	1785 1878 3380		
AlwNI	1	4205	CviJI	95		RcaI	4	583 4509 5517 5622		
ApaI	1	1396	CviRI	26		RsaI	4	165 1332 3595 5162		
ApaBI	2	869 2366	Ddel	11		SapI	1	3673		
ApoI	2	1460 5675	DpnI	29		Sau96I	21			
AvaI	1	2739	DraI	3	4548 4567 5259	Sau3AI	29			
Avall	9	1737 2113 2201 2450 2753	DrdI	2	3482 3897	Scal	1	5162		
		2795 3074 4820 5042	DrdII	1	908	ScrFI	24			
BamHI	1	319	Dsal	2	622 2761	SfaNI	24			
BanI	12		EaeI	6	493 625 1859 2253 2758	Sfcl	5	138 431 4054 4245 4923		
BanII	3	569 583 1396			5070	SgrAI	1	504		
BbsI	5	1331 1670 2044 2907 5660	EagI	1	2253	SphI	1	660		
BbvI	27		Eam1105I	1	4682	SspI	1	5486		
BccI	15		EarI	3	803 3673 5477	StyI	2	244 2683		
Bce83I	7	208 1999 2169 3880 4178	Ecil	5	962 2709 3863 4009 4837	TaqI	12			
		4419 5287	Eco47III	3	590 2091 3043	TaqII	8	1093 1311 1984 3691 5030		
Bcefl	5	704 1045 1672 2481 4291	Eco57I	2	4337 5349			5215 5368 5385		
Bcgl	10	315 349 1477 1511 2011	EcoNI	1	720	TfiI	7	1864 2166 2320 2618 2839		
		2045 3366 3400 5187 5221	EcoO109I	5	240 618 2753 2795 5658			3343 3764		
BclI	1	1199	EcoRI	1	5675	Thal	40			
Bfal	7	257 353 398 2803 4284	EcoRII	10	129 908 1223 1763 1820	Tsel	27			
		4537 4872			2372 2755 3815 3936 3949	Tsp45I	9	124 1366 2194 2461 3228		
BglI	3	2249 2483 4802	EcoRV	2	187 1635			3441 3536 4938 5149		
BglII	1	463	FauI	18		Tsp509I	16			
BmgI	1	1394	FokI	14		Tth111I	1	3534		
Bpml	6	1023 1512 2146 2700 3316	FspI	3	2672 2770 4904	Tth111II	7	1024 1717 3250 4379 4386		
		4752	GdIII	5	493 625 1859 2253 5070			4418 5674		
Bpu10I	1	2895	HaeI	8	913 2234 2306 2363 2760	UbaII	23			
Bpu1102I	1	267			3804 3815 4267	VspI	4	446 1870 1929 4854		
BsaI	1	4743	HaeII	13		XbaI	1	397		
BsaAI	1	3541	HaeIII	27		XcmI	3	1041 1557 1575		
BsaBI	3	462 468 2986	Hgal	15		XmnI	2	3347 5281		
BsaHI	8	508 529 643 1142 1825	HgiEII	2	783 4375					
		2520 5219 5601	Hhal	44		Enzymes that do not cut pET-11a:				
BsaJI	10	115 129 244 622 628	Hin4I	5	16 1084 2455 4681 4755	AflII	Agel	AscI	AvrII	BaeI
		1820 2481 2683 2761 3949	HincII	2	1691 5223	BseRI	BsrGI	Bsu36I	DraIII	FseI
BsaWI	7	189 1504 2007 2978 3995	HindIII	1	29	KpnI	MunI	NcoI	NotI	NsiI
		4142 4973	Hinfl	14		NspV	Pacl	PmeI	PmlI	RleAI
BsaXI	1	1844	HpaI	1	1691	RsrII	Sacl	SacII	Sall	SexAI
Bsbl	2	3505 5225	HphI	17		Sfil	Sgfl	Smal	SnaBI	SpeI
BscGI	13		MaeII	12		SrfI	Sse8387I	StuI	SunI	Swal
BsgI	3	1036 1236 2949	MaeIII	18		XhoI				
Bsil	3	3962 5346 5653	MbolI	15						
BsiEI	6	1970 2256 3705 4129 5052	MluI	1	1185					
		5201	MmeI	2	4004 4188					
BsII	22		MnlI	33						
BsmI	1	2673	MscI	1	2760					
BsmAI	7	882 1287 1413 1800 3430	Msel	24						
		4743 5519	Msil	10	1237 1525 1555 2345 2776					
BsmBI	2	1800 3430			2971 3362 4934 5093 5452					
BsmFI	4	646 2187 2412 3060	MspI	35						
BsoFI	50		MspAI	10	271 1215 1785 1878 2455					
Bsp24I	14				3380 3499 4131 4376 5317					
Bsp1286I	11		MwoI	45						
BspEI	2	189 2978	NarI	5	508 529 643 1825 2520					
BspGI	3	2373 2450 3315	NciI	14						
BspLU11I	1	3789	NdeI	1	359					
BspMI	1	2368	NgoAIV	4	495 2083 2243 2597					
BsrI	26		NheI	1	352					
BsrBI	3	418 3722 5523	NlaIII	31						
BsrDI	4	1232 1598 4743 4917	NlaIV	28						
			NruI	1	2288					