

pET-22b(+) Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AccI	2	180 2985	BsrFI	8	231 424 433 800 2012	NotI	1	166		
AccIII	8	881 1609 1940 2724 2865				NspI	4	589 2560 2852 3219		
		3167 4407 5091	BssHII	1	1525	Pfi1108I	2	2001 4126		
AcI	81		Bst1107I	1	2986	PfiMI	1	696		
AflIII	2	1114 3215	BstEII	1	1295	PleI	9	375 663 750 1546 3109		
AluI	25		BstXI	3	916 1045 1168			3594 4097 5186 5194		
AlwI	16		BstYI	12		PshAI	1	1959		
Alw21I	9	159 190 614 1098 2209	Cac8I	39		Psp5II	1	2221		
		3033 3533 4694 4779	CjeI	24		Psp1406I	6	776 2144 2540 4334 4707		
Alw44I	4	1094 3029 3529 4775	CjePI	18				5036		
AlwNI	1	3631	CviJI	87		PstI	1	4353		
Apal	1	1325	CviRI	24		PvuI	1	4478		
ApaBI	1	798	DdeI	11		PvuII	3	1714 1807 2806		
ApoI	4	192 1389 5053 5064	DpnI	27		RcaI	3	512 3935 4943		
AvaI	1	158	DraI	3	3974 3993 4685	RsaI	3	1261 3021 4588		
Avall	7	1666 2042 2130 2221 2500	DrallI	1	5251	SacI	1	190		
		4246 4468	DrdI	3	2908 3323 5206	Sall	1	179		
BamHI	1	198	DrdII	2	837 5256	SapI	1	3099		
BanI	9	436 457 571 1034 1753	Dsal	3	220 551 2187	Sau96I	18			
		1883 2009 4056 5288	EaeI	6	166 223 422 554 1788	Sau3AI	27			
BanII	5	190 498 512 1325 5326			4496	Scal	1	4588		
BbsI	4	1260 1599 1973 2333	EagI	1	166	ScrFI	19			
BbvI	31		Eam1105I	1	4108	SfaNI	20			
BccI	14		EarI	3	732 3099 4903	Sfcl	5	360 3480 3671 4349 5470		
Bce83I	7	21 1928 2098 3306 3604	Ecil	4	891 3289 3435 4263	SgrAI	1	433		
		3845 4713	Eco47III	3	519 2020 2469	SphI	1	589		
BceII	5	633 974 1601 3717 5277	Eco57I	2	3763 4775	Sspl	2	4912 5043		
Bcgl	10	160 194 1406 1440 1940	EcoNI	1	649	StyI	2	57 220		
		1974 2792 2826 4613 4647	EcoO109I	3	53 547 2221	TaqI	13			
BclI	1	1128	EcoRI	1	192	TaqII	9	1022 1240 1913 3117 4456		
Bfal	7	70 327 2229 3710 3963	EcoRII	7	837 1152 1692 1749 3241			4641 4794 4811 5155		
		4298 5402			3362 3375	TfiI	5	1793 2095 2265 2769 3190		
BglI	3	237 2178 4228	EcoRV	2	217 1564	Thal	34			
BglII	1	392	FauI	17		Tsel	31			
BmgI	1	1323	FokI	10	1160 1169 2434 2496 2574	Tsp45I	8	1295 2123 2654 2867 2962		
Bpml	5	952 1441 2075 2742 4178			2760 2901 4074 4255 4542			4364 4575 5424		
Bpu10I	1	2321	Fspl	2	2196 4330	Tsp509I	18			
Bpu1102I	1	80	GdIII	5	166 422 554 1788 4496	Tth111I	1	2960		
Bsal	1	4169	HaeI	6	225 842 2163 3230 3241	Tth111II	6	953 1646 2676 3805 3812		
BsaAI	2	2967 5251			3693			3844		
BsaBI	3	391 397 2412	HaeII	14		UbaII	20			
BsaHI	6	437 458 572 1071 1754	HaeIII	25		VspI	5	207 375 1799 1858 4280		
		4645	Hgal	12		XbaI	1	326		
BsaJI	7	57 220 551 557 1749	HgiEII	2	712 3801	XcmI	3	970 1486 1504		
		2187 3375	HhaI	45		XhoI	1	158		
BsaWI	7	2 1433 1936 2404 3421	Hin4I	3	1013 4107 4181	XmnI	3	208 2773 4707		
		3568 4399	HinCI	2	181 1620					
BsaXI	2	1773 5199	HindIII	1	173					
Bsbl	2	2931 5158	HinFI	14		Enzymes that do not cut pET-22b(+):				
BscGI	13		HpaI	1	1620	AatII	AflII	AgeI	AscI	AvrII
BseRI	1	260	HphI	16		BaeI	BsmI	BsrGI	Bsu36I	Clal
BsgI	3	965 1165 2375	Maell	15		FseI	KpnI	MunI	NheI	NruI
Bsil	2	3388 4772	MaeIII	18		Nsil	NspV	PacI	PmeI	PmlI
BsiEI	7	169 271 1899 3131 3555	MbolI	14		RleAI	RsrII	SacII	SexAI	SfiI
		4478 4627	MluI	1	1114	Sgfl	Smal	SnaBI	SpeI	SrfI
BsII	20		MmeI	3	3430 3614 5228	Sse8387I	StuI	SunI	Swal	
BsmAI	7	811 1216 1342 1729 2856	MnII	27						
		4169 4945	MscI	1	225					
BsmBI	2	1729 2856	MseI	29						
BsmFI	4	575 2116 2486 5466	MslI	9	1166 1454 1484 2202 2397					
BsoFI	51				2788 4360 4519 4878					
Bsp24I	10	404 436 955 987 1257	MspI	32						
		1289 3708 3740 3886 3918	MspA1I	10	84 267 1144 1714 1807					
Bsp1286I	13				2806 2925 3557 3802 4743					
BspEI	2	2 2404	MwoI	39						
BspGI	1	2741	NarI	4	437 458 572 1754					
BspLU11I	1	3215	NciI	12						
BspMI	1	268	NcoI	1	220					
BsrI	25		NdeI	1	288					
BsrBI	4	347 3148 4949 5395	NgoAIV	5	231 424 2012 2172 5352					
BsrDI	4	1161 1527 4169 4343	NlaIII	25						
			NlaIV	25						