



# pET-17b Restriction Sites

Enzyme	# Sites	Locations	Enzyme	# Sites	Locations	Enzyme	# Sites	Locations		
AatII	1	3233	DpnI	19		Scal	1	2791		
AccI	1	1188	DraI	3	2177 2196 2888	ScrFI	11			
AceIII	4	927 1068 1370 2610	DrdI	2	1111 1526	SfaNI	13			
Acil	44		Dsal	1	390	Sfcl	5	169 332 1683 1874 2552		
AflIII	1	1418	EaeI	4	147 191 387 2699	SpeI	1	199		
AluI	16		EagI	2	147 191	SspI	1	3115		
AlwI	13		Eam1105I	1	2311	StyI	1	59		
Alw21I	6	215 412 1236 1736 2897	EarI	2	1302 3106	TaqI	4	142 361 1518 2962		
		2982	Ecil	3	1492 1638 2466	TaqII	5	1320 2659 2844 2997 3014		
Alw44I	3	1232 1732 2978	Eco47III	1	672	TfiI	4	229 468 972 1393		
AlwNI	1	1834	Eco57I	2	1966 2978	Thal	12			
ApoI	1	174	EcoO109I	4	55 382 424 3287	TseI	20			
AvaI	2	141 368	EcoRI	1	174	Tsp45I	5	857 1070 1165 2567 2778		
Avall	5	382 424 703 2449 2671	EcoRII	4	384 1444 1565 1578	Tsp509I	7	174 298 348 2178 2484		
BamHI	1	205	EcoRV	1	166			2739 3304		
BanI	2	217 2259	FauI	6	348 433 714 900 1121	Tth111I	1	1163		
BanII	1	215			1131	Tth111III	4	879 2008 2015 2047		
BbsI	2	536 3289	FokI	8	637 699 777 963 1104	UbaII	11			
BbvI	20				2277 2458 2745	VspI	2	347 2483		
BccI	4	162 2348 2472 2759	FspI	2	399 2533	XbaI	1	306		
Bce83I	5	23 1509 1807 2048 2916	GdiII	3	147 191 2699	XhoI	1	141		
BceII	2	179 1920	HaeI	4	389 1433 1444 1896	XmnI	2	976 2910		
BcgI	6	224 258 995 1029 2816	HaeII	4	591 674 1296 1666					
		2850	HaeIII	13						
Bfal	8	72 200 262 307 432	HgaI	5	954 1111 1529 2107 2837	Enzymes that do not cut pET-17b:				
		1913 2166 2501	HgiEI	1	2004	AflII	AgeI	ApaI	ApaBI	AscI
BglI	1	2431	HhaI	17		AvrII	BaeI	BclI	BmgI	BsaXI
BglII	1	364	Hin4I	3	210 2310 2384	BseRI	BsmI	BspMI	BsrGI	BssHII
BpmI	2	945 2381	HincII	1	2852	BstEII	Bsu36I	Clal	DrallI	DrdII
Bpu10I	1	524	HindIII	1	223	EcoNI	FseI	Hpal	MluI	MunI
Bpu1102I	1	82	Hinfl	8	229 339 468 972 1318	NarI	NcoI	NgoAIV	NruI	NsiI
Bsal	2	331 2372			1393 1789 2306	NspV	PacI	PfiMI	PmeI	PmlI
BsaAI	1	1170	HphI	8	463 1038 1047 2154 2381	PshAI	RleAI	RsrII	SacII	Sall
BsaBI	3	164 363 615			2797 3003 3038	SexAI	SfiI	SgfI	SgrAI	SmaI
BsaHI	2	2848 3230	KpnI	1	221	SnaBI	SphI	SrfI	Sse8387I	StuI
BsaJI	3	59 390 1578	Maell	8	489 513 743 1169 2121	SunI	Swal	XcmI		
BsaWI	5	4 607 1624 1771 2602			2537 2910 3230					
Bsbl	2	1134 2854	MaeIII	14						
BscGI	9	46 394 779 1112 1745	MbolI	8	536 1289 2080 2151 2906					
		2091 2312 2336 2858			2984 3093 3289					
Bsgl	1	578	MmeI	2	1633 1817					
Bsil	3	1591 2975 3282	MnlI	19						
BsiEI	6	150 194 1334 1758 2681	MscI	1	389					
		2830	MseI	16						
BsII	10	52 391 754 845 887	MslI	6	405 600 991 2563 2722					
		1067 1440 1458 1624 1903			3081					
BsmAI	4	331 1059 2372 3148	MspI	16						
BsmBI	1	1059	MspA1I	6	86 1009 1128 1760 2005					
BsmFI	1	689			2946					
BsoFI	31		MwoI	15						
Bsp24I	8	231 263 1911 1943 2089	NciI	7	428 756 1062 1097 1798					
		2121 3215 3247			2494 2845					
Bsp1286I	6	215 412 1236 1736 2897	NdeI	1	268					
		2982	NheI	1	261					
BspEI	2	4 607	NlaIII	16						
BspGI	1	944	NlaIV	13						
BspLU11I	1	1418	Nott	1	147					
BsrI	16		Nspl	3	763 1055 1422					
BsrBI	3	146 1351 3152	Pfi1108I	1	2329					
BsrDI	2	2372 2546	PleI	4	347 1312 1797 2300					
BsrFI	1	2391	Psp5II	2	382 424					
Bst1107I	1	1189	Psp1406I	3	743 2537 2910					
BstXI	2	160 186	PstI	2	173 2556					
BstYI	10	134 205 364 610 2059	PvuI	1	2681					
		2070 2156 2168 2936 2953	PvuII	1	1009					
Cac8I	12		RcaI	3	2138 3146 3251					
CjeI	14		RsaI	3	219 1224 2791					
CjePI	14		SacI	1	215					
CviJI	52		SapI	1	1302					
CviRI	14		Sau96I	10	55 382 424 703 890					
Ddel	10	82 103 524 686 1226			2353 2432 2449 2671 3287					
		1693 2102 2268 2808 3234	Sau3AI	19						