

pET-32 Ek/LIC Restriction Sites

Enzyme	# Sites	Locations				
AccI	2	180	3409			
AccIII	8	1305	2033	2364	3148	3289
		3591	4831	5515		
Acil	85					
AfIII	2	1538	3639			
AluI	25					
AlwI	17					
Alw21I	10	159	190	431	1038	1522
		2633	3457	3957	5118	5203
Alw44I	5	427	1518	3453	3953	5199
AlwNI	1	4055				
ApaI	1	1749				
ApaBI	1	1222				
ApoI	5	192	287	1813	5477	5488
AvaI	2	158	228			
Avall	8	606	2090	2466	2554	2645
		2924	4670	4892		
BamHI	1	198				
BanI	11					
BanII	6	190	235	922	936	1749
		5750				
BbsI	4	1684	2023	2397	2757	
BbvI	29					
BccI	14					
Bce83I	7	21	2352	2522	3730	4028
		4269	5137			
BceII	5	1057	1398	2025	4141	5701
BcgI	14					
BclI	1	1552				
Bfal	7	70	747	2653	4134	4387
		4722	5826			
BglI	2	2602	4652			
BglIII	1	258				
BmgI	1	1747				
BpmI	5	1376	1865	2499	3166	4602
Bpu10I	1	2745				
Bpu1102I	1	80				
BsaI	1	4593				
BsaAI	2	3391	5675			
BsaBI	3	811	821	2836		
BsaHI	6	861	882	996	1495	2178
		5069				
BsaJI	10	57	211	228	486	513
		975	981	2173	2611	3799
BsaWI	7	2	1857	2360	2828	3845
		3992	4823			
BsaXI	2	2197	5623			
Bsbl	2	3355	5582			
BscGI	14					
BseRI	1	234				
BsgI	3	1389	1589	2799		
Bsil	2	3812	5196			
BsiEI	6	169	2323	3555	3979	4902
		5051				
BsII	25					
BsmAI	7	1235	1640	1766	2153	3280
		4593	5369			
BsmBI	2	2153	3280			
BsmFI	4	999	2540	2910	5890	
BsoFI	50					
Bsp24I	10	828	860	1379	1411	1681
		1713	4132	4164	4310	4342
Bsp1286I	15					
BspEI	2	2	2828			
BspGI	1	3165				
BspLU11I	1	3639				
BsrI	25					
BsrBI	4	767	3572	5373	5819	
BsrDI	4	1585	1951	4593	4767	
BsrFI	8	380	848	857	1224	2436
		2596	4612	5776		
BssHII	1	1949				

Enzyme	# Sites	Locations				
Bst1107I	1	3410				
BstEII	1	1719				
BstXI	3	1340	1469	1592		
BstYI	12					
CacBI	39					
CjeI	26					
CjePI	18					
Clal	2	525	815			
CviJI	90					
CvIRI	28					
DdeI	11					
DpnI	32					
DraI	3	4398	4417	5109		
DraIII	1	5675				
DrdI	3	3332	3747	5630		
DrdII	3	374	1261	5680		
Dsal	4	211	486	975	2611	
EaeI	7	166	366	382	846	978
		2212	4920			
EagI	1	166				
Eam1105I	1	4532				
EarI	3	1156	3523	5327		
EcII	4	1315	3713	3859	4687	
Eco47III	3	943	2444	2893		
Eco57I	2	4187	5199			
EcoNI	1	1073				
EcoO109I	3	53	971	2645		
EcoRI	1	192				
EcoRII	9	384	512	1261	1576	2116
		2173	3665	3786	3799	
EcoRV	1	206				
FauI	17					
FokI	13					
FspI	2	2620	4754			
GdIII	6	166	382	846	978	2212
		4920				
HaeI	6	368	1266	2587	3654	3665
		4117				
HaeII	14					
HaeIII	26					
HgaI	13					
HgiEII	3	436	1136	4225		
HhaI	46					
Hin4I	4	203	1437	4531	4605	
HincII	2	181	2044			
HindIII	1	173				
HinfI	16					
HpaI	1	2044				
HphI	18					
KpnI	1	255				
MaeII	15					
MaeIII	18					
MbolI	15					
MluI	1	1538				
MmeI	3	3854	4038	5652		
MnlI	29					
MscI	1	368				
MseI	28					
MsiI	9	1590	1878	1908	2626	2821
		3212	4784	4943	5302	
MspI	33					
MspA1I	10	84	300	1568	2138	2231
		3230	3349	3981	4226	5167
MwoI	41					
NarI	4	861	882	996	2178	
NciI	14					
NcoI	1	211				
NdeI	2	363	708			
NgoAIV	4	848	2436	2596	5776	
NlaIII	26					
NlaIV	27					
NotI	1	166				

Enzyme	# Sites	Locations				
NspI	4	1013	2984	3276	3643	
NspV	1	285				
Pfi1108I	2	2425	4550			
PfiMI	2	277	1120			
PleI	10	483	795	1087	1174	1970
		3533	4018	4521	5610	5618
PshAI	1	2383				
Psp5II	1	2645				
Psp1406I	6	1200	2568	2964	4758	5131
		5460				
PstI	1	4777				
PvuI	1	4902				
PvuII	3	2138	2231	3230		
RcaI	3	936	4359	5367		
RsaI	5	253	659	1685	3445	5012
RsrII	1	606				
SacI	1	190				
Sall	1	179				
SapI	1	3523				
Sau96I	19					
Sau3AI	32					
Scal	1	5012				
ScrFI	23					
SfaNI	21					
Sfci	5	780	3904	4095	4773	5894
SgrAI	1	857				
SmaI	1	230				
SphI	1	1013				
SrfI	1	230				
SspI	2	5336	5467			
StyI	2	57	211			
TaqI	18					
TaqII	10	446	1446	1664	2337	3541
		4880	5065	5218	5235	5579
TfiI	6	583	2217	2519	2689	3193
		3614				
Thal	35					
Tsel	29					
Tsp45I	8	1719	2547	3078	3291	3386
		4788	4999	5848		
Tsp509I	18					
Tth111I	1	3384				
Tth111II	6	1377	2070	3100	4229	4236
		4268				
UbaII	21					
VspI	4	795	2223	2282	4704	
XbaI	1	746				
XcmI	3	1394	1910	1928		
XhoI	1	158				
XmnI	3	405	3197	5131		

Enzymes that do not cut pET-32 Ek/LIC:

AatII	AflII	AgeI	AscI	AvrII
BaeI	BsmI	BspMI	BsrGI	Bsu36I
FseI	MunI	NheI	NruI	NsiI
PacI	PmeI	PmlI	RleAI	SacII
SexAI	SfiI	SgfI	SnaBI	SpeI
Sse8387I	StuI	SunI	Swal	