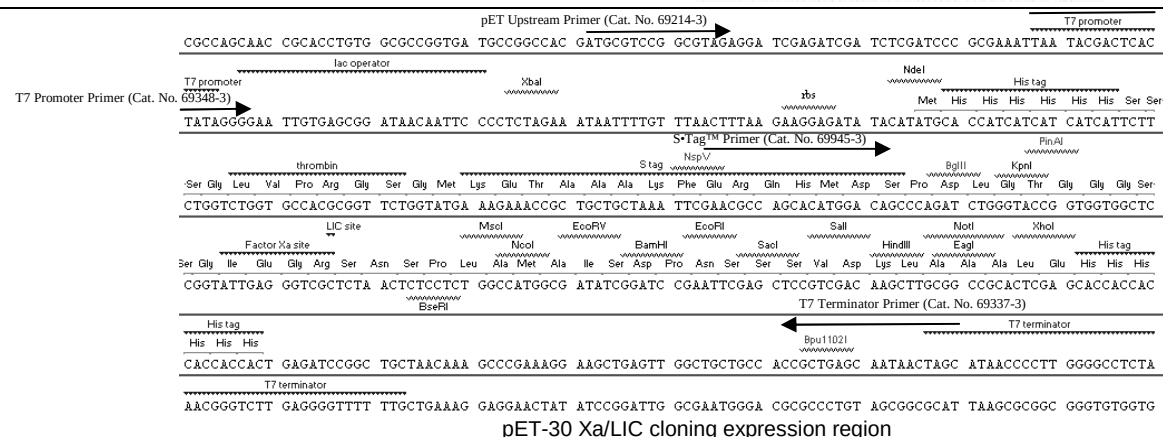
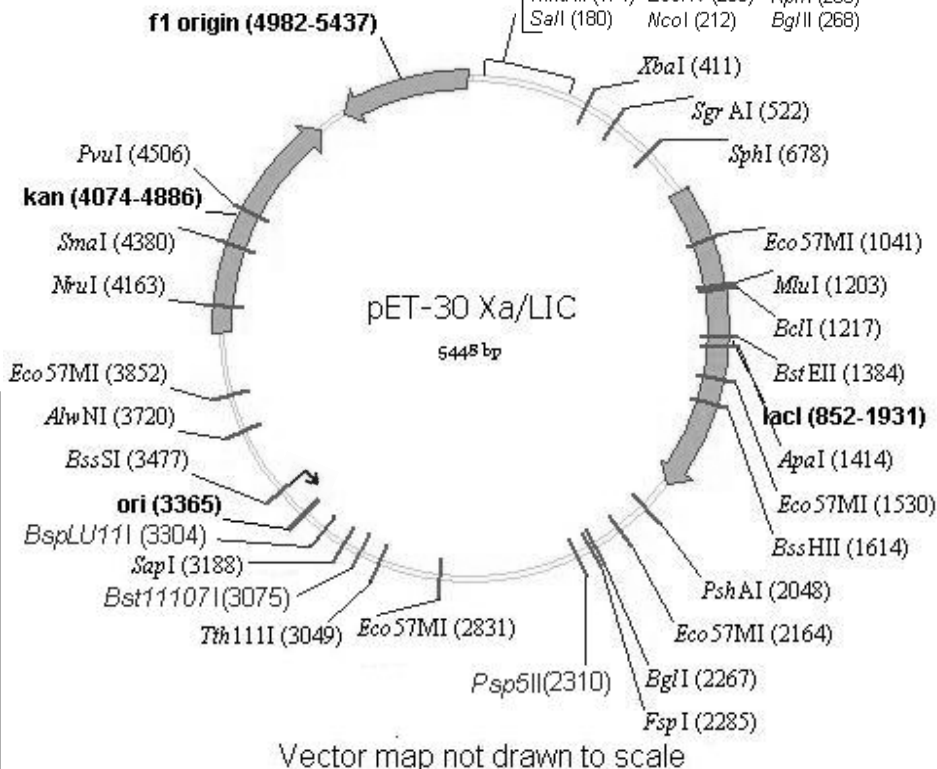


T7 promoter	445-463
T7 transcription start	444
His•Tag coding sequence	353-370
S•Tag coding sequence	275-319
Multiple cloning sites (BseR I - <i>Xho</i> I)	158-225
His•Tag coding sequence	140-157
T7 terminator	26-72
<i>lacI</i> coding sequence	852-1931
pBR322 origin	3365
<i>Kan</i> coding sequence	4074-4886
<i>f1</i> origin	4982-5437

All corresponding maps and tables are generated using VectorNTI[®] software. Restriction sites are numbered in accordance with the conventions of the respective software.

TCC
AGGCCATAACTCCCA

<i>Xho</i> I (159)	<i>Sac</i> I (191)	<i>Msc</i> I (217)	<i>Nsp</i> V (295)
<i>Eag</i> I (167)	<i>Eco</i> R I (193)	<i>Bse</i> R I (236)	<i>Nde</i> I (373)
<i>Not</i> I (167)	<i>Bam</i> H I (199)	<i>Pin</i> A I (258)	
<i>Hind</i> III (174)	<i>Eco</i> R V (206)	<i>Kpn</i> I (265)	
<i>Sal</i> I (180)	<i>Nco</i> I (212)	<i>Bal</i> I (268)	



Contact Your Local Distributor
www.emd4biosciences.com
bioscienceshelp@
emdchemicals.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.

Enzyme	#	Locations						Enzyme	#	Locations						Enzyme	#	Locations					
AccI	2	181	3074					Bst1107I	1	3075						NlaIV	24						
Acil	75							BstEII	1	1384						NotI	1	167					
AflIII	2	1203	3304					BstXI	3	1005	1134	1257				NruI	1	4163					
AluI	22							BstYI	9	133	199	268	767	1979		NsiI	2	4356	4622				
AlwI	13									2496	3945	3956	4755			NspI	4	678	2649	2941	3308		
Alw26I	6	900	1305	1430	1818	2945		Cac8I	40							NspV	1	295					
		4522						ClaI	2	480	4197					PflMI	3	287	785	4769			
AlwNI	1	3720						CviJI	85							PinAI	1	258					
Apal	1	1414						DdeI	11							PleI	9	460	752	839	1635	3198	
ApaLI	3	1183	3118	3618				DpnI	23									3683	4738	5142	5150		
ApoI	7	193	297	1478	4119	4303		DrallI	1	5207						PshAI	1	2048					
		5009	5020					DrdI	3	2997	3412	5162				Psp1406I	4	865	2233	2629	4992		
AvaI	2	159	4378					DsaI	3	212	640	2276				PspSII	1	2310					
AvaII	5	1755	2131	2219	2310	2589		EaeI	5	167	215	511	643	1877		PvuI	1	4506					
BamHI	1	199						EagI	1	167						PvuII	3	1803	1896	2895			
BanI	10	261	337	525	546	660		EarI	3	821	3188	4319				RcaI	3	601	4024	4899			
		1123	1842	1972	2098	5244		Eco47III	3	608	2109	2558				RsaI	4	263	1350	3110	4341		
BanII	6	191	587	601	1414	4161		Eco57I	1	3852						SacI	1	191					
		5282						EcoNI	2	738	4418					Sall	1	180					
BbsI	4	1349	1688	2062	2422			EcoO109I	3	54	636	2310				SapI	1	3188					
BbvI	25							EcoRI	1	193						Sau3AI	23						
BcgI	4	195	1495	2063	2915			EcoRII	9	926	1241	1781	1838	3330		Sau96I	14						
BcgI*	4	161	1529	2029	2881					3451	3464	4394	4751			ScrFI	21						
BclI	1	1217						EcoRV	1	207						SfaNI	23						
BfaI	6	71	412	2318	3799	4106		FauI	17							StcI	4	445	3569	3760	5426		
		5358						Fnu4HI	43							SgfI	1	4506					
BglI	1	2267						FokI	9	1249	1258	2523	2585	2663		SgrAI	1	522					
BglII	1	268								2849	2990	4144	4750			SmaI	1	4380					
BpmI	4	1041	1530	2164	2831			FspI	1	2285						SphI	1	678					
Bpu10I	2	2410	4523					HaeII	14							Sspl	2	4431	4999				
Bpu1102I	1	81						HaeIII	24							StyI	2	58	212				
BsaAI	2	3056	5207					HgaI	11							TaiI	14						
BsaBI	3	476	486	2501				HhaI	46							TaqI	17						
BsaHI	5	526	547	661	1160	1843		HincII	2	182	1709					TfiI	9	1882	2184	2354	2858	3279	
BsaJI	10	58	212	640	646	1838		HindIII	1	174								4417	4473	4645	4736		
		2276	3464	4377	4378	4779		HinfI	18							ThaI	36						
BsaWI	9	3	246	258	1522	2025		HpaI	1	1709						TseI	25						
		2493	3510	3657	4641			HphI	16							Tsp45I	7	1384	2212	2743	2956	3051	
BseRI	1	236						KpnI	1	265								4653	5380				
BsgI	3	1054	1254	2464				MaeIII	16							Tsp509I	21						
BsiEI	5	170	1988	3220	3644	4506		MboII	13							TspRI	13						
BsiHKA I	7	160	191	703	1187	2298		MluI	1	1203						Tth111I	1	3049					
		3122	3622					MnlI	27							VspI	5	460	1888	1947	4705	4894	
BslI	26							MscI	1	217						XbaI	1	411					
BsmI	2	4390	4467					MseI	25							XcmI	3	1059	1575	1593			
BsmBI	3	1818	2945	4522				MSI	6	1255	1543	1573	2291	2486		XhoI	1	159					
BsmFI	4	664	2205	2575	5422					2877						XmnI	2	2862	4895				
Bsp1286I	12							MspI	31														
BspEI	2	3	2493					MspAI	9	85	310	1233	1803	1896									
BspLU11I	1	3304								2895	3014	3646	3891										
BsrI	21							MwoI	39														
BsrBI	4	432	3237	4905	5350			NarI	4	526	547	661	1843										
BsrDI	2	1250	1616					NciI	12														
BsrFI	8	258	513	522	888	2100		NcoI	1	212													
		2261	4460	5308				NdeI	1	373													
BssHII	1	1614						NgoAIV	4	513	2101	2261	5308										
BssSI	1	3477						NlaIII	26														

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

AatII	AflII	AhdI	AscI	AvrII
BsaI	BspMI	BsrGI	Bsu36I	DraI
FseI	MunI	NheI	PacI	PmeI
PmlI	PstI	RsrII	SacI	SanDI
ScaI	SexAI	SfiI	SnaBI	SpeI
SrfI	Sse8387I	StuI	SunI	Swal
UbaEI				

© 2010 EMD Chemicals Inc., an Affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. The Novagen® name and Novagen® logo are registered trademarks of EMD Chemicals Inc. in the United States and in certain other jurisdictions. VectorNTI® is a registered trademark of Life Technologies Corporation in the US and/or certain other countries.

The terms and conditions governing EMD Chemicals' sale of this product are as indicated on our website (www.emdbiosciences.com).

USA and Canada
Tel (800) 628-8470
bioscienceshelp@emdchemicals.com

Germany
Tel 0800 100 3496
techservice@merckbiosciences.de

United Kingdom and Ireland
UK Freephone 0800 622935
Ireland Toll Free 1800 409445
customer.service@merckbiosciences.co.uk

All Other Countries
www.emd4biosciences.com
bioscienceshelp@emdchemicals.com