

Enzyme	#	Locations					Enzyme	#	Locations					Enzyme	#	Locations				
AarI	1	2680					BsgI	6	2870	4587	8049	8994	11668	NaeI	8	622	2398	5492	5616	5623
AatII	1	203							11868							6432	10348	12405		
Acc65I	1	7934					BsiEI	19						NarI	4	11075	12257	12371	12392	
AccI	5	3278	4463	4522	9621	10390	BsiHKAII	10	328	993	4983	7325	8817	NcoI	2	3537	9914			
AcII	3	3753	8326	12053					10399	10407	10817	11737	12221	NdeI	2	2993	7732			
AfeI	2	7293	12312				BsiWI	1	220					NgoMIV	8	620	2396	5490	5614	5621
AflII	3	1037	9335	9575			BsmAI	12								6430	10346	12403		
AflIII	5	1871	7700	7874	10209	11713	BsmBI	5	2318	3614	4167	6669	11098	NheI	1	97				
AgeI	4	1810	6554	7784	8677		BsmFI	26						NotI	1	239				
AhdI	2	2845	10213				BsmI	4	1833	2813	3444	3851		NruI	4	924	1444	4264	10401	
AleI	2	3119	9270				Bsp1286I	20						NsiI	1	77				
Alol	2	4791	8428				BspCNI	25						NspI	5	75	1104	4451	10213	12246
AlwNI	8	2764	2938	4351	5068	7836	BspEI	5	264	477	1451	3842	8494	PacI	1	289				
		7856	8878	10740			BspHI	5	4814	7508	8162	9358	12315	PciI	1	10209				
ApaI	2	9699	11510				BspMI	6	2321	2680	3130	5193	6741	PfIMI	8	138	1656	2764	3613	4180
ApalI	1	11733							10508							4953	8026	12138		
AscI	1	208					BsrBI	6	14	581	2547	3073	4280	PfoI	3	2901	9478	12145		
Asel	7	4769	7726	8218	10489	10971			4902					PmlI	1	216				
		11030	12458				BsrDI	11						PpuMI	5	2404	2761	2935	4348	10585
AsiSI	1	189					BsrFI	19						PshAI	9	4499	9744	9787	9830	9873
AvaI	7	180	274	321	890	5012	BsrGI	1	6507							9916	9959	10045	10131	
		6218	8844				BssHII	5	208	1406	8191	8735	11302	PsiI	2	5184	5849			
AvrII	1	310					BssSI	3	7884	9534	10097			PspOMI	2	9695	11506			
BaeI	3	1007	8222	10488			BstAPI	4	739	4565	10372	12037		PstI	3	2691	6752	8293		
BamHI	1	1975					BstBI	1	229					PvuI	2	189	8600			
BanI	14						BstEII	3	1813	10331	11531			PvuII	5	3942	6702	8708	11024	11117
BanII	9	181	654	9699	10399	10407	BstXI	4	7812	11667	11790	11919		SacI	3	10399	10407	10817		
		10817	11510	12323	12337		BstYI	15						SacII	1	5629				
BbeI	4	11078	12260	12374	12395		BstZ17I	1	4464					Sall	2	3277	10389			
BbsI	5	7777	7966	8843	11228	11567	Bsu36I	1	294					SapI	3	1799	7330	8540		
BceAI	30						BtgI	19						SbfI	1	6752				
BcgI	6	867	1234	2701	4464	10364	BtlI	4	1774	8286	10986	11354		Scal	1	3425				
		11393					Clal	2	194	12438				SexAI	2	3207	10327			
BciVI	3	1683	5614	11261			DraI	7	303	915	3583	3922	5197	SfiI	8	502	2687	2822	6280	6748
BclI	2	8525	11699						7360	10705						8289	10224	12471		
BfrBI	1	75					DraIII	2	4565	10550				SfiI	1	2435				
BglI	3	2435	5898	10347			DrdI	5	1770	4341	5981	6876	8172	SfoI	4	11076	12258	12372	12393	
BglII	2	5873	7940				EaeI	21						SgrAI	6	2396	2645	5614	7784	8941
BlpI	2	400	7206				EagI	9	239	725	2289	5050	5095			12394				
Bme1580I	7	2865	3375	5123	5399	9699			5492	5619	5623	5684		SmaI	2	182	6220			
		11510	11737				EarI	8	1799	3153	6376	6416	7330	SmlI	13					
BmgBI	6	2617	2629	2847	5486	7764			8540	8950	12096			SnaBI	1	8358				
		10524					Ecil	10	1078	1206	2128	3080	4416	SpeI	1	9449				
Bmrl	6	9745	9874	10915	11555	11792			5255	5552	6142	6724	11928	SphI	1	75				
		12189					Eco57MI	15						SrfI	1	182				
BmtI	1	101					EcoICRI	3	10397	10405	10815			Sse8387I	1	6752				
BpII	2	8575	8857				EcoNI	2	6196	12181				SspI	7	1372	3530	5189	6332	9166
Bpml	7	3722	7191	7604	8570	9501	EcoO109I	8	427	2404	2761	2935	3229			9193	9685			
		11392	11881						4348	10585	12281			StuI	1	5901				
Bpu10I	3	4066	6654	7849			EcoRI	2	2051	3838				StyI	7	310	422	3537	4632	5813
BpuEI	8	464	1225	3059	5315	5361	EcoRV	4	1122	6855	7084	11267				6460	9914			
		5533	8304	10905			FseI	1	5625					Swal	1	303				
BsaAI	6	216	4149	7875	8358	8463	FspAI	1	6479					TaqII	15					
		9200					FspI	2	6479	10305				TatI	3	1155	3423	6507		
BsaBI	6	977	2351	10481	10565	12434	HaeII	14						TspGWI	13					
		12444					HincII	9	2620	3279	5522	5557	6252	Tth111I	1	7998				
BsaHI	8	200	1929	2649	11075	11758			8933	10356	10391	11211		XbaI	2	31	5919			
		12257	12371	12392			HindIII	1	232					XcmI	4	5424	11328	11346	11862	
BsaI	1	9537					HpaI	2	10356	11211				XhoI	2	321	5012			
BsaWI	15						KasI	4	11074	12256	12370	12391		XmaI	2	180	6218			
BsaXI	7	2029	4770	7577	8295	8980	KpnI	1	7938					XmnI	4	6385	6455	9166	10650	
		9474	11045				MfeI	4	810	1234	1744	7714		ZraI	1	201				
BseRI	1	8427					MluI	2	1871	11713										
BseYI	7	1589	3251	5694	8617	9374	MscI	3	3575	5476	8145									
		11179	11314				MslI	16												

Enzymes that do not cut pETcoco™-1:
BbvCI Fall PmeI PpiI PsrI
RsrII SanDI

© 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.merck4biosciences.com
bioscienceshelp@
emdchemicals.com