



Enzyme	Cleavage Site 5 → 3	JBS Enzyme	Enzyme	Cleavage Site 5 → 3	JBS Enzyme	Enzyme	Cleavage Site 5 → 3	JBS Enzyme
Aat I	AGG↓CCT	SseB I	BstMA I	CTGCA↓G	Pst I	Nco I	C↓CATGG	Nco I
Acc I	GT↓MKAC	Acc I	BstMB I	↓GATC	{Dpn I}, Mbo I, Sau3A I	Nde I	CA↓TATG	Nde I
Acc III	T↓CCGGA	BseA I	BstN I	CC↓WGG	BseB I	Nde II	↓GATC	{Dpn I}, Mbo I, Sau3A I
Acc113 I	AGT↓ACT	Sca I	BstO I	CC↓WGG	BseB I	NgoM IV	G↓CCGGC	{Nae I}
Acc65 I	G↓GTACC	{Kpn I}	BstP I	G↓GTNACC	BstE II	Nhe I	G↓CTAGC	Nhe I
Afa I	GT↓AC	Rsa I	BstSN I	TAC↓GTA	SnaB I	Not I	GC↓GGCCGC	Not I
Afl I	G↓GWCC	Sin I	BstX I	CCANNNN↓NTGG	BstX I	Nru I	TCG↓CGA	Nru I
Afl II	C↓TTAAG	MspC I	Bsu15 I	AT↓CGAT	BseC I	Nsp III	C↓YCGRG	Ava I
Age I	A↓CCGGT	CspA I	BsuR I	GG↓CC	BshF I	Nsp V	TT↓CGAA	Asu II
Aha III	TTT↓AAA	Dra I	BsuTU I	AT↓CGAT	BseC I	Pae I	CGATG↓C	Sph I
Ahl I	A↓CTAGT	Spe I	BtsC I	GGATG (2/0)	{Fok I}	PaeR7 I	C↓TCGAG	Sla I
Ajn I	↓CCWGG	{BseB I}	CciN I	GC↓GGCCGC	Not I	Pal I	GG↓CC	BshF I
Alu I	AG↓CT	Alu I	Cfr10 I	R↓CCGGY	BssA I	Pau I	G↓CCGCGC	BssH II
Alw44 I	G↓TGCAC	ApaI I	Cfr13 I	G↓GNCC	PspP I	Pce I	AGG↓CCT	SseB I
Ama87 I	C↓YCGRG	Ava I	Cfr42 I	CCGC↓GG	SgrB I	Pdi I	GCC↓GGC	Nae I
Apa I	GGGCC↓C	Apa I	Cfr9 I	C↓CCGGG	{Sma I}	PfiF I	GACN↓NNGTC	Tth111 I
ApaL I	G↓TGCAC	ApaL I	Cla I	AT↓CGAT	BseC I	Pho I	GG↓CC	BshF I
AsiA I	A↓CCGGT	CspA I	Cpo I	CG↓GWCCG	Rsr II	PinA I	A↓CCGGT	CspA I
Asp I	GACN↓NNGTC	Tth111 I	Csp I	CG↓GWCCG	Rsr II	Ple19 I	CGAT↓CG	Pvu I
Asp718 I	G↓GTACC	{Kpn I}	Csp45 I	TT↓CGAA	Asu II	Psp124B I	GAGCT↓C	Sst I
AspS9 I	G↓GNCC	PspP I	Csp6 I	G↓TAC	{Rsa I}	Psp6 I	↓CCWGG	{BseB I}
Asu II	TT↓CGAA	Asu II	CspA I	A↓CCGGT	CspA I	PspA I	C↓CCGGG	{Sma I}
AsuNH I	G↓CTAGC	Nhe I	Din I	GGC↓GCC	{Nar I}	PspE I	G↓GTNACC	BstE II
Ava I	C↓YCGRG	Ava I	Dpn I	GA↓TC	Dpn I, Mbo I, Sau3A I	PspG I	↓CCWGG	{BseB I}
Ava II	G↓GWCC	Sin I	Dpn II	↓GATC	{Dpn I}, Mbo I, Sau3A I	PspOM I	G↓GGCCC	{Apa I}
BamH I	G↓GATCC	BamH I	Dra I	TTT↓AAA	Dra I	PspP I	G↓GNCC	PspP I
Ban II	GRG↓C	Ban II	Ecl136 II	GAG↓CTC	{Sst I}	Pst I	CTGCA↓G	Pst I
Ban III	AT↓CGAT	BseC I	Eco105 I	TAC↓GTA	SnaB I	Psy I	GACN↓NNGTC	Tth111 I
Bbe I	GGCG↓C	{Nar I}	Eco130 I	C↓CWWGG	Sty I	Pvu I	CGAT↓CG	Pvu I
Bbu I	GCATG↓C	Sph I	Eco147 I	AGG↓CCT	SseB I	Pvu II	CAG↓CTG	Pvu II
Bcl I	T↓GATCA	Bcl I	Eco24 I	GRG↓C	Ban II	Rsa I	GT↓AC	Rsa I
Bco I	C↓YCGRG	Ava I	Eco32 I	GAT↓ATC	EcoRV	Rsr II	CG↓GWCCG	Rsr II
Bcu I	A↓CTAGT	Spe I	Eco47 I	G↓GWCC	Sin I	Rsr2 I	CG↓GWCCG	Rsr II
Bfr I	C↓TTAAG	MspC I	Eco88 I	C↓YCGRG	Ava I	Sac I	GAGCT↓C	Sst I
BfuC I	↓GATC	{Dpn I}, Mbo I, Sau3A I	Eco91 I	G↓GTNACC	BstE II	Sac II	CCGC↓GG	SgrB I
Bgl I	GCCNNNN↓NGGC	Bgl I	EcoICR I	GAG↓CTC	{Sst I}	Sal I	GTCGAC	Sal I
Bgl II	A↓GATCT	Bgl II	EcoO65 I	G↓GTNACC	BstE II	Sau3A I	↓GATC	{Dpn I}, Mbo I, Sau3A I
Bme18 I	G↓GWCC	Sin I	EcoR I	G↓AATTC	EcoR I	Sau96 I	G↓GNCC	PspP I
BmeT110 I	CY↓CGRG	{Ava I}	EcoR II	↓CCWGG	{BseB I}	Sca I	AGT↓ACT	Sca I
Bmt I	GCTAG↓C	{Nhe I}	EcoR I	G↓AATTC	EcoR I	Sfi I	GGCCNNNN↓NGGCC	Sfi I
Bpu14 I	TT↓CGAA	Asu II	EcoR V	GAT↓ATC	EcoRV	Sfo I	GGC↓GCC	{Nar I}
BpvU I	CGAT↓CG	Pvu I	EcoT14 I	C↓CWWGG	Sty I	Sfr274 I	C↓TCGAG	Sla I
Bsa29 I	AT↓CGAT	BseC I	EcoT38 I	GRG↓C	Ban II	Sfr303 I	CCGC↓GG	SgrB I
Bse118 I	R↓CCGGY	BssA I	Ege I	GGC↓GCC	{Nar I}	Sfu I	TT↓CGAA	Asu II
BseA I	T↓CCGGA	BseA I	Ehe I	GGC↓GCC	{Nar I}	SgrB I	CCGC↓GG	SgrB I
BseB I	CC↓WGG	BseB I	Erh I	C↓CWWGG	Sty I	Sin I	G↓GWCC	Sin I
BseC I	AT↓CGAT	BseC I	FauND I	CA↓TATG	Nde I	Sla I	C↓TCGAG	Sla I
BseG I	GGATG (2/0)	{Fok I}	Fba I	T↓GATCA	Bcl I	Sma I	CCC↓GGG	Sma I
BseP I	G↓CCGCGC	BssH II	Fbl I	GT↓MKAC	Acc I	SnaB I	TAC↓GTA	SnaB I
BshF I	GG↓CC	BshF I	Fok I	GGATG (9/13)	Fok I	SpaH I	GCATG↓C	Sph I
BshT I	A↓CCGGT	CspA I	FriO I	GRG↓C	Ban II	Spe I	A↓CTAGT	Spe I
BsiHKC I	C↓YCGRG	Ava I	Fun II	G↓AATTC	EcoR I	Sph I	GCATG↓C	Sph I
BsiS I	C↓CGG	BsiS I	Hae III	GG↓CC	BshF I	SseB I	AGG↓CCT	SseB I
BsoB I	C↓YCGRG	Ava I	Hap II	C↓CGG	BsiS I	Ssp I	AAT↓ATT	Ssp I
Bsp106 I	AT↓CGAT	BseC I	Hinc II	GTY↓RAC	Hinc II	Sst I	GAGCT↓C	Sst I
Bsp119 I	TT↓CGAA	Asu II	Hind II	GTY↓RAC	Hinc II	Stu I	AGG↓CCT	SseB I
Bsp120 I	G↓GGCCC	{Apa I}	Hind III	A↓AGCTT	Hind III	Sty I	C↓CWWGG	Sty I
Bsp13 I	T↓CCGGA	BseA I	Hinf I	G↓ANTC	Hinf I	Taq I	T↓CGA	Taq I
Bsp143 I	↓GATC	{Dpn I}, Mbo I, Sau3A I	Hpa I	GTT↓AAC	Hpa I	Tel I	GACN↓NNGTC	Tth111 I
Bsp19 I	C↓CTAGG	Nco I	Hpa II	C↓CGG	BsiS I	Tli I	C↓TCGAG	Sla I
Bsp68 I	TCG↓CGA	Nru I	Kas I	G↓CGCC	{Nar I}	Tth111 I	GACN↓NNGTC	Tth111 I
BspANI	GG↓CC	BshF I	Kpn I	GGTAC↓C	Kpn I	Vha464 I	C↓TTAAG	MspC I
BspC I	CGAT↓CG	Pvu I	Kpn2 I	T↓CCGGA	BseA I	VpaK11B I	G↓GWCC	Sin I
BspD I	AT↓CGAT	BseC I	Ksp I	CCGC↓GG	SgrB I	Xba I	T↓CTAGA	Xba I
BspE I	T↓CCGGA	BseA I	Ksp22 I	T↓GATCA	Bcl I	Xho I	C↓TCGAG	Sla I
BspT I	C↓TTAAG	MspC I	KspA I	GTT↓AAC	Hpa I	Xma I	C↓CCGGG	{Sma I}
BspT104 I	TT↓CGAA	Asu II	Kzo9 I	↓GATC	{Dpn I}, Mbo I, Sau3A I	XmaC I	C↓CCGGG	{Sma I}
BspX I	AT↓CGAT	BseC I	Mbo I	↓GATC	{Dpn I}, Mbo I, Sau3A I	Xmi I	GT↓MKAC	Acc I
BsrF I	R↓CCGGY	BssA I	Mbo II	GAAGA (8/7)	Mbo II	Zho I	AT↓CGAT	BseC I
BssA I	R↓CCGGY	BssA I	Mlu I	A↓CGCGT	Mlu I	Zrm I	AGT↓ACT	Sca I
BssH I	C↓TCGAG	Sla I	Mly113 I	GG↓CGCC	Nar I			
BssH II	G↓CCGCGC	BssH II	Mnl I	CCCT (7/6)	Mnl I			
BssT1 I	C↓CWWGG	Sty I	Mro I	T↓CCGGA	BseA I			
Bst2U I	CC↓WGG	BseB I	MroN I	G↓CCGGC	{Nae I}			
Bst98 I	C↓TTAAG	MspC I	Msp I	C↓CGG	BsiS I			
BstB I	TT↓CGAA	Asu II	MspC I	C↓TTAAG	MspC I			
BstE II	G↓GTNACC	BstE II	Mva I	CC↓WGG	BseB I			
BstEN II	↓GATC	{Dpn I}, Mbo I, Sau3A I	Mvr I	CGAT↓CG	Pvu I			
BstF5 I	GGATG (2/0)	{Fok I}	Nae I	GCC↓GGC	Nae I			
BstKI I	GAT↓C	{Dpn I}, {Mbo I}, {Sau3A I}	Nar I	GG↓CGCC	Nar I			

Enzyme available from Jena Bioscience Isoschizomer available from Jena Bioscience Neoschizomer available from Jena Bioscience

- Single Letter Code: R = A or G, Y = C or T, M = A or C, K = G or T, S = C or G, W = A or T, H = A or C or T, B = C or G or T, V = A or C or G, D = A or G or T, N = A or C or G or T
- Isoschizomers have same recognition sequence and cutting pattern.
- Neoschizomers (same recognition sequence but different cutting pattern) are indicated with brackets {enzyme}.