

Restriction Enzymes

Single buffer system: One for All - All for One

Achieve 100 % enzyme activity and complete digestion in **JBS Universal Buffer (UB)**. Save time with **JBSpeed restriction enzymes**.



Restriction Enzyme	Cat.-No.	Units		Reaction Conditions			Heat inactivation	Recognition Sequence (5'→3')	Methylation sensitivity	Compatible cohesive ends
		S-pack	L-pack	Buffer concentration	Reaction time	Reaction temperature				
AccI	EN-E2004	500	5x 500	1x	60 min	37 °C	20 min @ 80 °C	GTJMK AC	CpG	CviAI
AluI	EN-101	600	5x 600	1x	60 min	37 °C	20 min @ 65 °C	AGJCT	-	[blunt]
ApaI	EN-E2015	5000	5x 5000	1x	60 min	25 °C	20 min @ 65 °C	G GGCJC	CpG dcm	BanII, BspI286I
ApaLI	EN-172	2000	5x 2000	0.5x	60 min	37 °C	-	GJTGCA C	CpG	SfiI
AsuII	EN-102	3500	5x 3500	1x	60 min	37 °C	20 min @ 65 °C	TTJCG AA	CpG	AccI, AclI, AclI, BsaHI, BstBI, ClaI, MspI, NarI, TaqI
AvaI	EN-E2006	2000	5x 2000	1x	60 min	37 °C	20 min @ 65 °C	CJYCGR G	CpG	-
AvaII	EN-E2007	1000	5x 1000	1x	60 min	37 °C	20 min @ 65 °C	GJGWC C	CpG dcm	PpuMI
BamHI	EN-103	7500	5x 7500	1x	60 min	37 °C	20 min @ 80 °C	GJGATC C	-	BclI, DpnII, BglII, BstYI
BanII	EN-E2060	1000	5x 1000	1x	60 min	37 °C	20 min @ 65 °C	G RGCYC	-	ApaI
BclI	EN-104	2500	5x 2500	1x	60 min	50 °C	-	TJGATC A	dam	BamHI, BstYI, BglII, MboI
BglI	EN-105	2000	5x 2000	1x	60 min	37 °C	20 min @ 65 °C	GCCN NNNJNGGC	CpG	-
BglII	EN-106	1300	5x 1300	1x	60 min	37 °C	-	AJGATC T	-	BamHI, BstYI, BclI, DpnII
BseAI	EN-107	650	5x 650	1x	60 min	55 °C	-	TJCCGG A	CpG dcm dam	AgeI, Aval, AccIII, XmaI
BseBI	EN-108	4500	5x 4500	1x	60 min	60 °C	-	CCJW GG	-	-
BseCI	EN-109	3500	5x 3500	1x	60 min	55 °C	-	ATJCG AT	CpG dam	AccI, AclI, AclI, BsaHI, BstBI, ClaI, MspI, NarI, TaqI
BshFI	EN-110	7000	5x 7000	1x	60 min	37 °C	20 min @ 80 °C	GGJCC	-	[blunt]
BsiSI	EN-111	2200	5x 2000	1x	60 min	55 °C	-	CJCG G	-	AccI, AclI, AclI, BsaHI, BstBI, ClaI, MspI, NarI, TaqI
BssAI	EN-112	250	5x 250	1x	60 min	65 °C	-	RJCCGG Y	-	-
BssHII	EN-E2115	200	5x 200	1x	60 min	50 °C	20 min @ 80 °C	GJCGC C	CpG	MluI, AscI
BstEII	EN-144	1750	5x 1750	1x	60 min	60 °C	-	GJGTNAC C	-	-
BstXI	EN-E2118	1000	5x 1000	1x	60 min	50 °C	20 min @ 65 °C	CCAN NNNNJTGG	dcm	-
CspAI	EN-113	150	5x 150	1x	60 min	37 °C	20 min @ 65 °C	AJCCGG T	-	AgeI, Aval, AccIII, XmaI
DpnI	EN-160	200	5x 200	1x	60 min	37 °C	20 min @ 80 °C	GAJTC	CpG	BamHI, BstYI, BclI, BglII
DraI	EN-E2145	2000	5x 2000	1x	60 min	37 °C	20 min @ 65 °C	TTTJAAA	- EcoKI	[blunt]
EcoRI	EN-114	15000	5x 15000	1x	60 min	37 °C	20 min @ 65 °C	GJAATC C	CpG	ApoI, MfeI
EcoRV	EN-115	3000	5x 3000	1x	60 min	37 °C	20 min @ 80 °C	GATJATC	CpG	[blunt]
FokI	EN-E2710	500	5x 500	1x	60 min	37 °C	20 min @ 65 °C	GGATGNNNNNNNNNNN	CpG dcm	-
HincII	EN-E2200	1000	5x 1000	1x	60 min	37 °C	20 min @ 65 °C	GTJIRAC	CpG EcoKI	-
HindIII	EN-116	7500	5x 7500	1x	60 min	37 °C	20 min @ 80 °C	AJAGCT T	-	-
HinfI	EN-117	2500	5x 2500	1x	60 min	37 °C	20 min @ 80 °C	GJANT C	CpG	-
HpaI	EN-118	750	5x 750	1x	60 min	37 °C	-	GTTJAAAC	CpG	[blunt]
HpaII	EN-E2260	2000	5x 2000	1x	60 min	37 °C	20 min @ 65 °C	CJCG G	CpG	AccI, AclI, ClaI, BstBI, TaqI, AclI, BsaHI, HinPII, NarI
KpnI	EN-119	3500	5x 3500	1x	60 min	37 °C	-	G GTACJC	-	-
MboI	EN-120	300	5x 300	1x	60 min	37 °C	20 min @ 65 °C	JGATC	CpG dam	BamHI, BstYI, BclI, BglII
MbolI	EN-E2284	250	5x 250	1x	60 min	37 °C	20 min @ 65 °C	GAAGANNNNNN NJ	CpG dam	-
MluI	EN-E2287	1000	5x 1000	1x	60 min	37 °C	20 min @ 65 °C	AJCGC T	CpG EcoKI	AscI, BssHII
MnlI	EN-E2289	250	5x 250	1x	60 min	37 °C	20 min @ 65 °C	CCTNNNNNN NJ	CpG	-
MspCI	EN-121	1300	5x 1300	2x	60 min	37 °C	20 min @ 65 °C	CJTAA G	-	AflII, SmlI
NaeI	EN-122	300	5x 300	1x	60 min	37 °C	20 min @ 65 °C	GCCJGGC	CpG	[blunt]
NarI	EN-E2291	200	5x 200	0.5x	60 min	37 °C	20 min @ 65 °C	GGJCG CC	CpG	AccI, AclI, ClaI, BstBI, TaqI, AclI, HinPII, BsaHI, HpaII
NcoI	EN-123	600	5x 600	1x	60 min	37 °C	20 min @ 65 °C	CJCATG G	-	BspHI, FstI, PciI
NdeI	EN-E2293	2000	5x 2000	1x	60 min	37 °C	20 min @ 65 °C	CAJTA TG	-	AseI, BfaI, Csp6I, MseI
NheI	EN-146	550	5x 550	1x	60 min	37 °C	20 min @ 65 °C	GJCTAG C	CpG	AvrII, SpeI, StyI, XbaI
NotI	EN-124	300	5x 300	1x	60 min	37 °C	20 min @ 65 °C	GCJGGC GC	CpG	EagI, EaeI
NruI	EN-125	700	5x 700	1x	60 min	37 °C	20 min @ 65 °C	TCGJCGA	CpG dam	[blunt]
PspPI	EN-126	900	5x 900	1x	60 min	25 °C	15 min @ 55 °C	GJGNC C	dcm	-
PstI	EN-127	8000	5x 8000	1x	60 min	37 °C	15 min @ 55 °C	C TGCAJG	-	NsiI, SbfI
PvuI	EN-E2320	200	5x 200	1x	60 min	37 °C	20 min @ 80 °C	CG ATJCG	CpG dam	AsiSI, PacI, BsiEI
PvuII	EN-128	4500	5x 4500	1x	60 min	37 °C	-	CAGJCTG	-	[blunt]
RsaI	EN-129	1000	5x 1000	1x	60 min	37 °C	20 min @ 65 °C	GTJAC	CpG	[blunt]
SalI	EN-130	2000	5x 2000	2x	60 min	37 °C	20 min @ 65 °C	GJTCGA C	CpG	PspXI, XhoI
Sau3AI	EN-150	500	5x 500	1x	60 min	37 °C	20 min @ 65 °C	JGATC	CpG	BamHI, BclI, BglII, XhoII
Scal	EN-131	1200	5x 1200	1x	60 min	37 °C	20 min @ 80 °C	AGTJACT	-	[blunt]
SfiI	EN-132	400	5x 400	1x	60 min	50 °C	-	GCCN NNNJNGGCC	CpG dcm	-
SgrBI	EN-133	1600	5x 1600	1x	60 min	37 °C	20 min @ 65 °C	CC GCJGG	CpG	BsiEI, SacII
Slal	EN-134	5000	5x 5000	1x	60 min	37 °C	20 min @ 65 °C	CJTCGA G	CpG	AvaI, SalI, SmlI, XhoI
SmaI	EN-135	1100	5x 1100	1x	60 min	25 °C	20 min @ 65 °C	CCCJGGG	CpG	[blunt]
SnaBI	EN-136	350	5x 350	1x	60 min	37 °C	20 min @ 80 °C	TACJGTA	CpG	[blunt]
SpeI	EN-E2398	500	5x 500	1x	60 min	37 °C	20 min @ 65 °C	AJCTAG T	- EcoKI	AvrII, NheI, StyI, XbaI
SphI	EN-137	250	5x 250	2x	60 min	37 °C	20 min @ 65 °C	G CATGJC	-	NlaIII, NspI
SseBI	EN-138	1500	5x 1500	1x	60 min	37 °C	20 min @ 65 °C	AGGJCTT	dcm	[blunt]
SspI	EN-139	600	5x 600	1x	60 min	37 °C	20 min @ 65 °C	AATJATT	-	[blunt]
SstI	EN-140	1600	5x 1600	1x	60 min	37 °C	20 min @ 65 °C	G AGCTJC	-	BanII, SacI
StyI	EN-141	6000	5x 6000	2x	60 min	37 °C	20 min @ 65 °C	CJCWGG G	-	AvrII
TaqI	EN-142	3500	5x 3500	1x	60 min	65 °C	20 min @ 80 °C	TJCG A	dam	AccI, ClaI, BstBI, AclI, AclI, BsaHI, HinPII, HpaII, NarI
Tth111I	EN-E2420	1000	5x 1000	1x	60 min	65 °C	-	GACN JN NGTC	-	-
XbaI	EN-143	3500	5x 3500	1x	60 min	37 °C	20 min @ 65 °C	TJCTAG A	dam	AvrII, NheI, SpeI, StyI

Chart Legend

	JBSpeed enzyme (5-10 min reaction time)
	Cleavage inhibited/impaired by CpG methylation of recognition sequence (eukaryotic genome DNA)
	Cleavage inhibited/impaired by dcm methylation of recognition sequence (bacterial source DNA)
	Cleavage inhibited/impaired by dam methylation of recognition sequence (bacterial source DNA)
	Cleavage inhibited/impaired by EcoKI methylation of recognition sequence (bacterial source DNA)
	Potential inhibition by overlapping CpG methylation site
	Potential inhibition by overlapping dcm methylation site
	Potential inhibition by overlapping dam methylation site
	Potential inhibition by overlapping EcoKI methylation site

Nucleotide 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

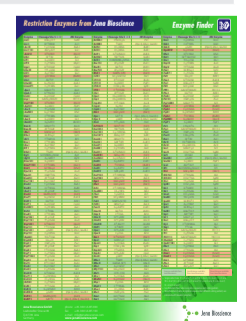
JBS Newsletter:
Learn about new products & special offers!



Find updates of this poster online or order printed copies here.



Check out Neo- & Isoschizomers in the JBS Enzyme Finder!



Jena Bioscience GmbH
Loebstedter Str. 71
07749 Jena
Germany

Phone +49(0)3641-62 85 000
Fax +49(0)3641-62 85 100
info@jenabioscience.com
www.jenabioscience.com



IFTA AG
Certified QMS and EMS according to
DIN EN ISO 9001 and DIN EN ISO 14001
Reg.-No.: ICV03597 034 and ICV03597 534