

Recommended Reaction Conditions for Jena Bioscience Restriction Endonucleases

Restriction enzyme	Tris-HCl (Tris-Acetate) {Bis Tris Propane-HCl}		NaCl (K-Acetate) {KCl}	MgCl ₂ (Mg-Acetate)	DTT	BSA	TX-100	Temp.	Buffer
	mM	pH (25°C)	mM	mM	mM	µg/ml	%	°C	
<i>Alu</i> I	10	7.9	-	10	1	100	-	37	B1
<i>Apa</i> L I	10	7.9	-	10	1	100	-	37	B1
<i>Asu</i> II	10	7.9	50	10	1	100	0.1	37	B2
<i>Bam</i> H I	10	7.9	100	5	1	100	-	37	U
<i>Bcl</i> I	10	7.9	50	10	1	100	-	50	B2
<i>Bgl</i> I	100	7.9	50	5	-	100	0.025	37	U
<i>Bgl</i> II	50	7.9	100	10	1	100	-	37	B3
<i>Bse</i> A I	10	8.0	100	5	1	100	0.02	55	U
<i>Bse</i> B I	10	7.9	50	10	1	100	-	60	B2
<i>Bse</i> C I	50	7.9	100	10	1	100	-	55	B3
<i>Bsh</i> F I	(20)	7.9	(50)	(10)	1	100	-	37	B5
<i>Bsi</i> S I	(33)	7.9	(66)	(10)	0.5	100	0.1	55	U
<i>Bss</i> A I	20	8.5	{100}	3	-	100	0.04	65	U
<i>Bst</i> E II	10	7.4	{100}	5	1	100	0.1	60	U
<i>Csp</i> A I	{10}	7.0	-	10	1	100	-	37	U
<i>Dpn</i> I	(33)	7.9	(66)	(10)	-	100	-	37	U
<i>Eco</i> R I	100	7.9	50	5	-	100	0.025	37	U
<i>Eco</i> R V	10	7.9	50	10	1	100	-	37	B2
<i>Hind</i> III	10	7.9	50	10	1	100	-	37	B2
<i>Hinf</i> I	50	7.9	100	10	1	100	-	37	B3
<i>Hpa</i> I	(20)	7.9	(50)	(10)	1	100	-	37	B5
<i>Kpn</i> I	10	7.0	-	10	1	100	0.01	37	U
<i>Mbo</i> I	10	8.0	{100}	10	1	100	-	37	U
<i>Msp</i> C I	10	7.9	150	10	1	100	-	37	B4
<i>Nae</i> I	10	7.9	-	10	1	100	-	37	B1
<i>Nco</i> I	50	7.9	100	10	1	100	0.02	37	B3
<i>Nhe</i> I	(20)	7.9	(50)	(10)	1	100	-	37	B5
<i>Not</i> I	50	7.9	100	5	1	100	-	37	BU
<i>Nru</i> I	50	8.0	{100}	10	-	100	-	37	U
<i>Psp</i> P I	10	7.9	50	10	1	100	-	25	B2
<i>Pst</i> I	50	7.4	100	10	1	100	-	37	U
<i>Pvu</i> II	10	7.9	50	10	1	100	-	37	B2
<i>Rsa</i> I	10	7.9	50	10	1	100	-	37	B2
<i>Sal</i> I	10	7.9	150	10	1	100	-	37	B4
<i>Sau</i> 3A I	10	7.9	50	10	-	100	-	37	B2
<i>Sca</i> I	10	7.4	100	10	1	100	-	37	B2
<i>Sfi</i> I	10	7.9	50	10	1	100	-	50	B2
<i>Sgr</i> B I	10	7.9	-	10	1	100	0.1	37	B1
<i>Sla</i> I	10	7.9	150	10	1	100	-	37	B4
<i>Sma</i> I	(20)	7.9	(50)	(10)	1	100	-	25	B5
<i>Sna</i> B I	{10}	7.0	-	10	1	100	-	37	U
<i>Sph</i> I	10	7.9	50	10	1	100	-	37	B2
<i>Sse</i> B I	50	7.9	100	10	1	100	-	37	B3
<i>Ssp</i> I	50	7.9	100	10	1	100	-	37	B3
<i>Sst</i> I	10	7.9	-	10	1	100	-	37	B1
<i>Sty</i> I	50	7.9	100	10	1	100	-	37	B3
<i>Taq</i> I	10	8.4	100	10	1	100	-	65	U
<i>Xba</i> I	10	7.9	50	10	1	100	-	37	B2

Requires Triton X-100 for optimal activity. TX-100 is included into the supplied reaction buffer.

Note:

BSA is not included into the supplied reaction buffers and should be added separately from the supplied stock solution (1 mg/ml).

For more information please contact enzymes@jenabioscience.com.

Jena Bioscience GmbH
Löbstedter Str. 80
07749 Jena, Germany
Tel.:+49-3641-6285 000
Fax:+49-3641-6285100
e-mail: info@jenabioscience.com