

Performance Chart for Restriction Enzymes



Enzyme	Supplied NEBuffer	% Activity in NEBuffers			Incub. Temp. (°C)	Inact. Temp. (°C)	Dil.	Unit Substrate	Methylation Sensitivity	Notes
		r1.1	r2.1	r3.1						
BamHI	rCutsmart	50	50	100	37	80	A	λ DNA	OC	
BclII	rCutsmart	100	100	100	37	80	A	λ DNA	OC	
BglII	rCutsmart	100	100	100	37	80	A	λ DNA	OC	
BspI	rCutsmart	100	100	100	37	80	A	λ DNA	OC	
BstI	rCutsmart	100	100	100	37	80	A	λ DNA	OC	

Chart Legend and Notes:

- Recombinant
- Time-Saver qualified (cleaves substrate in 5 – 15 min under recommended conditions)
- Engineered for maximum performance
- dam methylation sensitivity
- dcm methylation sensitivity
- CpG methylation sensitivity (applies to eukaryotic genomic DNA only)
- Indicates that the restriction enzyme requires two or more sites for cleavage
- Supplied with its own unique reaction buffer that is different from the four standard NEBuffers. Compatibility with the four standard NEBuffers is indicated by the chart.
- Supplied with a separate vial of S-adenosylmethionine (SAM) or Adenosine 5'-Triphosphate (ATP). To obtain 100% activity, SAM or ATP should be added to the 1X reaction mix as specified on the product data card.

The following notes appear with any enzymes having ligation efficiencies lower than 100% as assessed by ligation and recutting:

- Ligation is < 10%
- Ligation is 25% – 75%
- Recutting after ligation is < 5%
- Recutting after ligation is 50% – 75%
- Ligation and recutting after ligation is not applicable since the enzyme is either a nicking enzyme, is affected by methylation, or if the enzyme cleaves outside its recognition sequence.

The following notes appear with any enzymes when star activity is a concern:

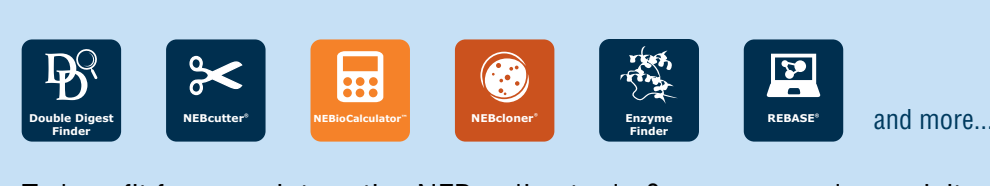
- Star activity may result from extended digestion, high enzyme concentration or a glycerol concentration of > 5%.
- Star activity may result from extended digestion.
- Star activity may result from a glycerol concentration of > 5%.
- May exhibit star activity in this buffer.

*For added flexibility, NEB offers an isochizomer or HF enzyme, supplied with rCutsmart Buffer.

To address the need for BSA-free reagents, NEB has switched our BSA-containing reaction buffers to Recombinant Albumin (rAlbumin)-containing buffers. We are also in the process of transitioning our enzyme formulations to contain rAlbumin. NEB has rigorously tested these changes and has not seen a difference in performance with these changes. Learn more at www.neb.com/BSA-free.

NEBuffer Compositions (1X)

- NEBuffer r1.1: 10 mM Bis Tris Propane-HCl, 10 mM MgCl₂, 100 μg/ml Recombinant Albumin (pH 7.0 @ 25°C)
- NEBuffer r2.1: 50 mM NaCl, 10 mM Tris-HCl, 10 mM MgCl₂, 100 μg/ml Recombinant Albumin (pH 7.9 @ 25°C)
- NEBuffer r3.1: 100 mM NaCl, 50 mM Tris-HCl, 10 mM MgCl₂, 100 μg/ml Recombinant Albumin (pH 7.9 @ 25°C)
- rCutsmart Buffer: 50 mM Potassium acetate, 20 mM Tris-acetate, 10 mM Magnesium acetate, 100 μg/ml Recombinant Albumin (pH 7.9 @ 25°C)



To benefit from our interactive NEB online tools & resources, please visit www.neb.com/nebtools

Activity of DNA Modifying Enzymes in rCutsmart Buffer

A selection of DNA modifying enzymes were assayed in rCutsmart Buffer and functional activity was compared to the activity in their supplied buffers. Reactions were set up according to the recommended reaction conditions, with rCutsmart Buffer (plus required supplement) replacing the supplied buffer.

Tech Tip: When supplements are required, one can simply add the supplied buffer of the respective modifying enzyme at 1X concentration to the rCutsmart Buffer to achieve appropriate activity for most applications – no change of buffers needed.

Enzyme	Activity in rCutsmart	Required Supplements	Enzyme	Activity in rCutsmart	Required Supplements
Phosphatases:			Nucleases, other:		
Antarctic Phosphatase	+++	Requires Zn ²⁺	DNAse I (RNase-free)	+++	Requires Ca ²⁺
Shrimp Alkaline Phosphatase (SAP)	+++		DNAse I-XT	+++	Requires Ca ²⁺
Quick CIP	+++		Exonuclease III (Nth)	+++	Requires Ca ²⁺
Ligases:			Exonuclease VII	+++	
T4 DNA Ligase	+++	Requires ATP	Exonuclease V (Rec BCD)	+++	Requires ATP
Hi-T4 [®] DNA Ligase	+++	Requires ATP	Exonuclease III	+++	
SaI [®] -T4 [®] DNA Ligase	+++	Requires ATP	Exonuclease I	+++	
E. coli DNA Ligase	+++	Requires NAD	PPase	+++	
T3 DNA Ligase	+++	Requires ATP + PEG	Lambda Exonuclease	+++	
T7 DNA Ligase	+++	Requires ATP + PEG	McrBC	+++	
Polymerases:			Micrococcal Nuclease	+++	Requires Ca ²⁺
T4 DNA Polymerase	+++	Requires ATP	RecI	+++	
DNA Polymerase I, Large (Klenow) Fragment	+++	A λ DNA	T5 Exonuclease	+++	
DNA Polymerase I	+++		T7 Exonuclease	+++	
DNA Polymerase Klenow Exo-	+++		Thermobifida EosI	+++	
Bst DNA Polymerase	+++		Thermobifida USER II	+++	
phi29 DNA Polymerase	+++	Requires DTT	Thermobifida USER III	+++	
T7 DNA Polymerase (unmodified)	+++	Requires DTT	Thermobifida OGG	+++	
Transferrase/Kinases:			USER [®] Enzyme	+++	
T4 Polynucleotide Kinase	+++	Requires ATP + DTT			
T4 PNK (3' phosphatase minus)	+++	Requires ATP + DTT			
CpG Methyltransferase (M. SssI)	+++				
GpC Methyltransferase (M. CviPI)	+++	Requires DTT			
T4 Phage β-glucosyltransferase (14-50T)	+++				

+++ full functional activity
 ++ 50-100% functional activity
 + 0-50% functional activity