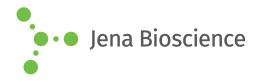
DATA SHEET





Dpnl

Neoschizomers: BfuCl, Bsp143l, BstENII, BstKTl,BstMBl, DpnII, Kzo9l, Ndell JBSpeed Restriction Enzyme

Cat. No.	Amount						
EN-160S	200 Units						
EN-160L	5 x 200 Units						
5'		G	А	↓	т	С	3'
3'		С	т	↑	A	G	5'

Unit Definition: One unit is the amount of enzyme required to completely digest 1 μ g of pBR322 (22 sites) in 1 hour in a total reaction volume of 50 μ l. Enzyme activity was determined in the recommended reaction buffer.

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Form: liquid (Supplied in 10 mM Tris-HCl pH 7.4, 400 mM KCl, 0.1 mM EDTA, 1 mM DTT, 200 $\mu g/ml$ BSA and 50 % [v/v] glycerol)

Concentration: 10 units/µl

Source: Diplococcus pneumoniae G41, recombinant, E. coli

Supplied with: 10x Universal Buffer (UB)

Recommended 50 µl assay

5 μl	10x Universal Buffer (UB)		
1 µg	pure DNA ¹ or PCR product ²		
10 units	enzyme		
fill up to 50 µl	PCR grade water		

¹ Supercoiled or high molecular weight DNA (e.g. plant genomic DNA) may require longer incubation time or higher amount of enzyme.
 ² Some enzymes may require additional DNA bases flanking the

² Some enzymes may require additional DNA bases flanking the restriction site for complete digestion.

Protocol:

- The enzyme should not exceed 10 % of total reaction volume.
- Add enzyme as last component. Mix components well before adding enzyme. After enzyme addition, mix gently by pipetting. Do not vortex.
- Incubate 5 to 10 min. at 37 °C.
- Stop reaction by alternatively:
 Addition of 2.1 μl EDTA pH 8.0 [0.5 M], final 20 mM
 Heat Inactivation (20 min. at 80 °C)
 - Spin Column DNA Purification (e.g. PCR Purification Kit, Cat.-No. PP-201S/L)

- Gel Electrophoresis and Single Band Excision (e.g. Agarose Gel Extraction Kit, Cat.-No. PP-202S/L)

- Phenol-Chloroform Extraction or Ethanol Precipitation.

Double Digestion - Buffer Compatibility:

B1 - 75-100 % Relative Activity

- B2 75-100 % Relative Activity
- B3 50-75 % Relative Activity
- B4 10 % Relative Activity
- B5 75-100 % Relative Activity
- 1x UB 100 % Relative Activity (recommended)

Please note that the optimum digestion condition for this enzyme is 1x UB. Within the Universal Buffer (UB) system, the most majority of our enzymes display 100% Relative Activity in 1x UB and only few either in 0.5x or 2x UB. If optimum condition for second enzyme is different than the recommended for the first enzyme, we suggest carrying out first the restriction at the higher recommended concentration of UB and dilute the reaction volume to the adequate UB concentration for further proceeding with the second restriction.



Jena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com