

Data Sheet

Lambda DNA/*Bst*E II Digest ready-to-use

Cat. No.	Amount
M-106S	100 µg
M-106L	500 µg

For *in vitro* use only
 Expiry: 18 months
 Storage: -20°C

Form

Liquid. Supplied in Jena Bioscience 6x Green Gel Loading Buffer (PCR-257-gr).

Concentration

0.4 µg/µl

Description

The *Bst*E II digest of λ DNA yields the following discrete fragments (in base pairs): 8453*, 7242, 6369, 5687*, 4822, 4324, 3675, 2323, 1929, 1371, 1264, 702, 224, 117.

Preparation

Lambda DNA was completely digested by *Bst*E I, phenol extracted, ethanol precipitated, dissolved in 10 mM Tris-HCl (pH 8.0) and 1 mM EDTA.

Note

Dilute in TE or other buffer of minimal ionic strength. DNA may denature if diluted in dH₂O and subsequently heated.

*The cohesive ends (12b *cos* site of bacteriophage) of fragments 8453 bp and 5687 bp may anneal to and form the additional band. These fragments may be separated by heating to 65°C for 5 min and then cooling on ice for 3 min.

