

Lambda DNA/*EcoR* I/ *Hind* III Digest ready-to-use

Cat. No.	Amount
M-103S	100 µg
M-103L	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

0.2 µg/µl

Description

The *EcoR* I and *Hind* III digest of λ DNA yields the following 13 discrete fragments (in base pairs): 21226*, 5148, 4973, 4268, 3530*, 2027, 1904, 1584, 1375, 947, 831, 564, 125.

Preparation

Lambda DNA was completely digested by *EcoR* I and *Hind* III, 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 21226 bp and 3530 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.

