

## Lambda DNA/*EcoR* I Digest ready-to-use

Cat. No.	Amount
M-102S	100 µg
M-102L	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 *EcoR* I digest of λ DNA yields the following 6 discrete fragments (in base pairs): 21226\*, 7421, 5804, 5643, 4878, 3530\*.

### Preparation

Lambda DNA was completely digested by *EcoR* 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 dH<sub>2</sub>O 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.

