

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

Cat. No.	Amount
M-104S	100 µg
M-104L	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 *Sty* I digest of λ DNA yields the following 11 discrete fragments (in base pairs): 19329\*, 7743, 6223, 4254\*, 3472, 2690, 1882, 1489, 925, 421, 74.

### Preparation

Lambda DNA was completely digested by *Sty* 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 19329 bp and 4254 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.

