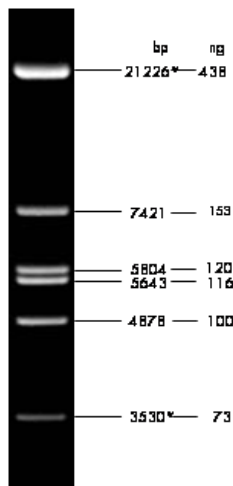




## λDNA/ EcoRI Digest

ready-to-use

Cat. No.	Amount
M-102S	50 µg
M-102L	5 x 50 µg



0.1 µg loaded onto 0.7 % agarose

**For *in vitro* use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at -20 °C

Short term storage (up to 1 month) at 4 °C possible.

**Shelf Life:** 12 months

**Form:** liquid (supplied in Jena Bioscience 6x Green Gel Loading Buffer (#PCR-254-gr))

**Concentration:** 50 ng/µl

### Description:

The EcoRI digest of λ DNA yields the following 6 discrete fragments (in base pairs): 21226\*, 7421, 5804, 5643, 4878, 3530\*.

The cohesive ends (12b cos site of bacteriophage lambda) of fragments 21226 bp and 3530 bp may anneal and form an additional band. These fragments may be separated again by heating to 65 °C for 5 min and then cooling on ice for 3 min.

### 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.

### Recommended Load:

2 - 5 µl per lane

### Preparation:

Lambda DNA (48.502 bp) was completely digested by EcoRI, phenol extracted, ethanol precipitated, dissolved in 10 mM Tris-HCl (pH 8.0) and 1 mM EDTA.