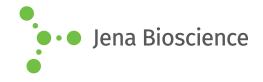
DATA SHEET

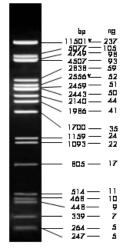




■ \(\lambda\text{DNA/ Pstl Digest}\)

ready-to-use

Cat. No.	Amount
M-105S	1 ml (100 ng/μl)
M-105L	5 x 1 ml (100 ng/μl)



 $5~\mu l$ loaded onto 1.5 % agarose

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

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

Shelf Life: 12 months

Form: liquid, DNA fragment mix in 10 mM Tris-HCl pH 7.5, 10 mM EDTA,

 $8\ \%$ (w/v) Glycerol, Tartrazine and Xylene Cyanol FF

Color: green

Concentration: 100 ng/µl

Description:

The PstI digest of λ DNA yields the following discrete fragments (in base pairs): 11501*, 5077, 4749, 4507, 2838, 2556*, 2459, 2443, 2140, 1986, 1700, 1159 1093, 805, 514, 468, 448, 339, 264, 247, 216, 211, 200, 164, 150, 94, 87, 72, 15.

The cohesive ends (12b cos site of bacteriophage) of fragments 11501 bp and 2556 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.

Recommended Load:

5 μl per lane

Preparation:

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