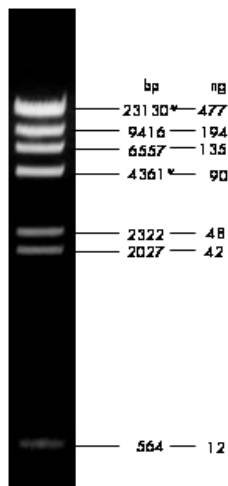




λDNA/ HindIII Digest

ready-to-use

Cat. No.	Amount
M-101S	100 µg
M-101L	5 x 100 µg



1 µg loaded onto 1.0 % agarose

For *in vitro* use only!

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 HindIII digest of λ DNA yields the following 8 discrete fragments (in base pairs): 23130*, 9416, 6557, 4361*, 2322, 2027, 564, 125.

The cohesive ends (12b cos site of bacteriophage lambda) of fragments 23130 bp and 4361 bp may be separated by heating to 65 °C for 5 min.

Note:

Dilute in TE or other buffer of minimal ionic strength. DNA may denature if diluted in dH₂O and subsequently heated.

Recommended Load:

5 µl per lane

Preparation:

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