

pBR322 DNA/*Hinf* I Digest ready-to-use

Cat. No.	Amount
M-107S	100 µg
M-107L	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 *Hinf* I digest of pBR322 DNA yields the following 10 discrete fragments (in base pairs): 1632, 517, 504, 396, 344, 298, 221, 220, 154, 75.

Preparation

pBR322 DNA was completely digested by *Hinf* I, phenol/chloroform 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 dH₂O and subsequently heated.

