

λ DNA/ BstE II Digest, S pack

The BstEII digest of λ DNA yields the following discrete fragments (in base pairs): 8453*, 7242, 6369, 5687*, 4822, 4324, 3675, 2323, 1929, 1371, 1264, 702, 224, 117. * The cohesive ends (12b cos site of bacteriophage) of fragments 8453 bp and 5687 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 dH₂O and subsequently heated.

Ordering Information

[Order Online »](#)

JBS-M-106S	100 μ g
------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com