



PRODUCT INFORMATION

# Lambda DNA

Pub. No. MAN0012976  
Rev. Date 06 September 2018 (Rev. B.00)

**#SD0011**

Lot \_\_\_\_ Expiry Date \_\_\_\_\_

Store at -25 °C to -15 °C

Components	#SD0011
Lambda DNA, 0.3 µg/µL	500 µg

**For Research Use Only.** Not for use in diagnostic procedures.

[www.thermofisher.com](http://www.thermofisher.com)

## Description

Linear double-stranded lambda bacteriophage (cI857 Sam7) DNA, 48502 base pairs with a molecular weight of  $31.5 \times 10^6$  Da. Isolated from a heat-inducible lysogenic *E.coli* W3110 strain. GeneBank/EMBL accession numbers NC\_001416, M17233, M24325, V00636, X00906. For map and features see [www.thermoscientific.com/reviewer](http://www.thermoscientific.com/reviewer).

## Applications

- Activity and specificity assays of restriction enzymes.
- Preparation of DNA molecular weight standards.
- Cloning.

## Storage Buffer

10 mM Tris-HCl (pH 7.6) and 1 mM EDTA.

## Important Licensing Information:

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

## PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to [www.thermofisher.com](http://www.thermofisher.com) for Material Safety Data Sheet of the product.

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.