

# BIO-PROBE<sup>®</sup> Lambda probe

Negative control probe for *in situ* hybridization

The Lambda BIO-PROBE<sup>®</sup> labeled probe is prepared by nick translation of the lambda bacteriophage strain c1857*ind1 Sam7*.

Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

## Ordering Information

[Order Online »](#)

ENZ-40840	2µg
-----------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application Notes** The BIO-PROBE<sup>®</sup> Lambda probe can be used in a variety of hybridization techniques, including Southern blots, Northern blots or dot blots, at concentrations of 50 to 100ng/ml. The probe can also be used as a negative control probe for *in situ* hybridization applications at concentrations of 0.5 to 2.0µg/ml.

The hybridized biotinylated DNA probe can be detected by its interaction with biotin-binding proteins, such as avidin, streptavidin or antibodies coupled to fluorescent dyes or color producing enzymes. Complete kits and protocols are available for *in situ* hybridization and detection procedures (Enzo's SAVIEW<sup>®</sup> PLUS and POLYVIEW<sup>®</sup> PLUS enhanced *in situ* detection systems).

**Concentration** 20µg/ml

**Formulation** Liquid. In 10mM TRIS-HCl, pH 7.5, containing 1mM EDTA.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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