# **BIO-PROBE**<sup>®</sup> Lambda probe

Negative control probe for in situ hybridization

The Lambda BIO-PROBE® labeled probe is prepared by nick translation of the lambda bacteriophage strain c1857ind1 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.

Last modified: May 29, 2024

