

# BIO-PROBE<sup>®</sup> Epstein-Barr virus (EBV) probe

Biotinylated Epstein-Barr virus (EBV) DNA probe for *in situ* hybridization

The BIO-PROBE<sup>®</sup> Epstein-Barr virus (EBV) labeled probe is prepared by nick translation of the 3Kb Bam HI "W" region (the internal repeat) of the EBV genome. The Bam HI clone is contained in pUC 18 vector. The probe does not cross hybridize with the other herpes virus DNAs, herpes simplex virus DNA or cytomegalovirus DNA.

## Ordering Information

[Order Online »](#)

ENZ-40836	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> Epstein-Barr virus (EBV) probe can be used in a variety of hybridization techniques, including Southern blots, Northern blots or dot blots, at concentrations of 50 to 100 ng/mL. The probe can also be used for *in situ* hybridization applications at concentrations of 0.5 to 2.0 µg/mL. Biotinylated probes have been shown to hybridize to homologous DNA at the same rate and to the same extent as non-biotinylated probes. 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.

**Concentration** 20 µg/mL

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

**Technical Info / Product Notes** BIO-PROBE<sup>®</sup> is a trademark of Enzo Life Sciences, Inc. Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.



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)