## BIO-PROBE<sup>®</sup> Cytomegalovirus probe

The Cytomegalovirus (CMV) BIO-PROBE<sup>®</sup> labeled probe is prepared by nick translation of cloned fragments of the CMV Towne strain. The fragments include a total of 30-31 kb of DNA, approximately 20% of the CMV genome. The probe is specific for cytomegalovirus DNA. It does not hybridize to DNAs of other herpesviruses (i.e., herpes simplex virus or Epstein-Barr virus) but it has been found to hybridize to the DNA of CMV clinical isolates. Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

Citations: 13

View Online »

**Ordering Information** 

Order Online »

**ENZ-40835** 2μg

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

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

after receipt when stored as recommended.

Long Term Storage -20°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application Notes** 

The Cytomegalovirus (CMV) BIO-PROBE<sup>®</sup> labeled 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 for in situ hybridization applications at concentrations of 0.5 to 1.0µg/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

20µg/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.

Last modified: May 29, 2024

