BIO-PROBE® Adenovirus probe

The Adenovirus BIO-PROBE[®] labeled probe is prepared by nick translation of cloned fragments of the adenovirus 5 genome. The fragments constitute a total of 20 kb of DNA, approximately 56% of the genome. The probe is specific for adenovirus and, in addition to hybridization to adenovirus 5 sequences, the probe will cross hybridize with adenovirus types 4, 7, 11, 20, 40 and 41 due to their considerable sequence homology. Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

Citations: 6

View Online »

Ordering Information

Order Online »

ENZ-40834 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 Adenovirus BIO-PROBE[®] 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. Complete kits and protocols are available for *in situ* hybridization and detection procedures (compatible with Enzo's POLYVIEW[®] PLUS and SAVIEW[®]

Concentration 20µg/ml

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

Technical Info / Product NotesBIO-PROBE[®] 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

