## BIO-PROBE<sup>®</sup> BK virus probe

The BK Virus BIO-PROBE<sup>®</sup> labeled probe is prepared by nick translation of a clone of the entire BK virus genome. BK virus, a simian polyoma virus, is related to SV40 and also bears some similarity to the human virus JC virus. BK virus is highly tumorigenic in hamsters. Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

Citations: 10

View Online »

**Ordering Information** 

Order Online »

ENZ-40848

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 BK Virus 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 2.0µg/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<sup>®</sup> PLUS and SAVIEW<sup>®</sup>

PLUS detection systems for inc).

Concentration 20µg/ml

Formulation Liquid. In 10mM TRIS HCI, 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



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com