BIO-PROBE[®] JC virus probe

The JC Virus BIO-PROBE[®] labeled probe is prepared by nick translation of a clone of the entire JC virus genome. JC virus, a member of the papovavirus family, usually causes subclinical infections. However, in immunocompromised individuals it can cause a fatal demyelinating brain disease, progressive multifocal leukoencephalopathy. JC virus is also neurooncogenic in primates. Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

Citations: 10

View Online »

Ordering Information

Order Online »

ENZ-40847 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 JC Virus BIO-PROBE[®] labeled probe can be used in a variety of hybridization techniques, including Southern blots, Northern blots or dot blots, at a concentration of approximately 50ng/ml. The probe can also be used for *in situ* hybridization applications at concentrations of approximately 0.5 to 5µg/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[®] PLUS detection systems for IHC).

Concentration

20µg/ml

Formulation

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

Technical Info / Product Notes

BIO-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



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