

# 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 <sup>®</sup> 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 <i>in situ</i> 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 <i>in situ</i> hybridization and detection procedures (compatible with Enzo's POLYVIEW <sup>®</sup> PLUS and SAVIEW <sup>®</sup> PLUS detection systems for IHC).
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.



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)