

BIO-PROBE® BK virus probe

The BK Virus BIO-PROBE® 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 [®] 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 <i>in situ</i> 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 <i>in situ</i> 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 HCl, 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.



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)