

BIO-PROBE[®] Blur 8 probe

Positive control probe for human DNA
sequences in both *in situ* and membrane
hybridization applications

The Blur 8 BIO-PROBE[®] labeled probe is prepared by nick translation of a clone of the human *alu* repeat sequence. Fragment size range: 100-1000 base pairs (as estimated by agarose gel electrophoresis).

Citations: 5

[View Online »](#)

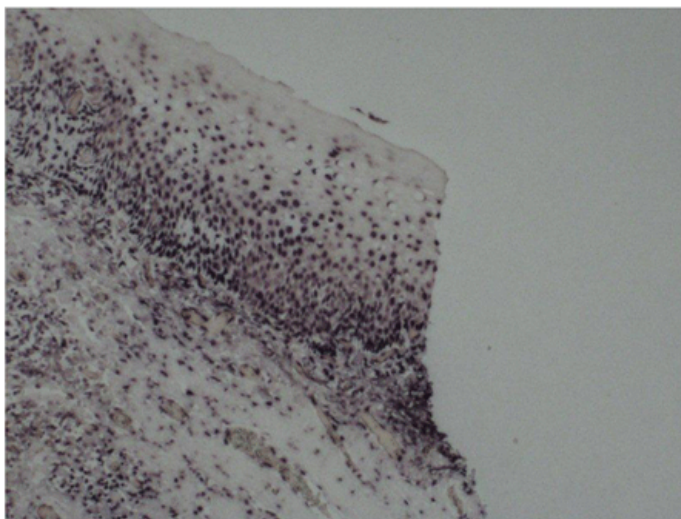
Ordering Information

[Order Online »](#)

ENZ-40849	2μg
-----------	-----

Manuals, SDS & CofA

[View Online »](#)



Blur 8 BIOPROBE® staining in tonsil tissue. Blur 8 BIOPROBE® was hybridized to tonsil tissue. SIMPLY SENSITIVE® AP-NBT/BCIP in situ detection system (ENZ-32870) was used for visualization.

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 Blur 8 BIO-PROBE[®] labeled probe can serve as a positive control probe for human DNA sequences in both *in situ* and membrane hybridization applications. It can be used at a concentration of approximately 50ng/ml for Southern blots, Northern blots or dot blots. The probe can also be used for *in situ* hybridization applications at concentrations of approximately 0.5 to 5.0µg/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.

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.

Last modified: May 29, 2024



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)