

PATHO-GENE[®] HPV

Type 6 Probe

Analyte Specific Reagent.

The PATHO-GENE[®] HPV Type 6 Probe is a biotin-labeled DNA probe in buffer containing formamide, salts and dextran sulfate.

Ordering Information

[Order Online »](#)

ENZ-GEN150-1000	1ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Do not use after expiration date.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status ASR - Analyte Specific Reagent

Product Details

Application Notes

The ENZO PATHO-GENE® HPV Type 6 Probe is a biotin-labeled DNA probe designed to hybridize to HPV Type 6 DNA sequences. The probe is denatured and hybridized to the tissue section fixed and pretreated on the microscope slide. Specific hybridization between the HPV DNA probe and DNA in the specimen is determined by the detection of biotin.

Technical Info / Product Notes

1. This product is an Analyte Specific Reagent (ASR). Analytical and performance characteristics are not established.
2. The probe has affinity to HPV 6 genotype and may cross-react with known or uncharacterized viral types. It is the responsibility of the user to validate any test for its specific use.
3. Specimens and all materials exposed to them should be considered potentially biohazardous, handled and disposed with proper precautions. Consult local and/or state authorities for the recommended method of disposal. Avoid contacting the skin and mucous membranes with reagents and specimens. Follow standard laboratory precautions to prevent exposure to eyes and skin. If reagents or specimens come in contact with sensitive areas, wash with copious amounts of water.
4. Avoid microbial contamination of reagents as it may cause incorrect results.
5. See also the product Safety Data Sheet (SDS).

This product or the use of this product is covered by one or more Enzo patents, including, but not limited to the following: U.S. Patent No. 4,994,373; Canadian Patent No. 1,309,672; 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)