

PATHO-GENE[®] HPV

type 16/18/31/33/51

probe

HPV probe that can be optimized for use with
polymer detection system or used as

optimized RTU with Enzo's Streptavidin
detection methods.

The PATHO-GENE[®] HPV type 16/18/31/33/51 probe is a mixture of biotin-labeled HPV-16, HPV- 18, HPV-31, HPV-33 and HPV-51 specific probes in buffered formamide and hybridization enhancers.

Citations: 5

[View Online »](#)

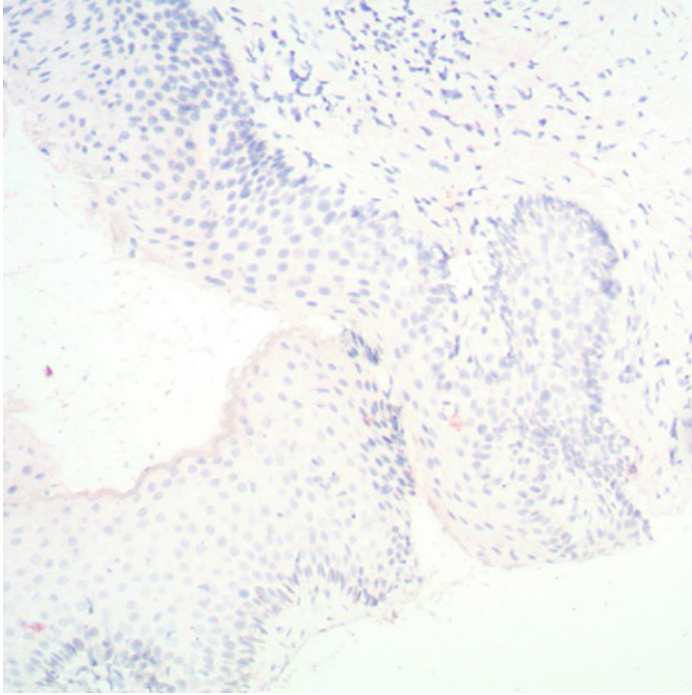
Ordering Information

[Order Online »](#)

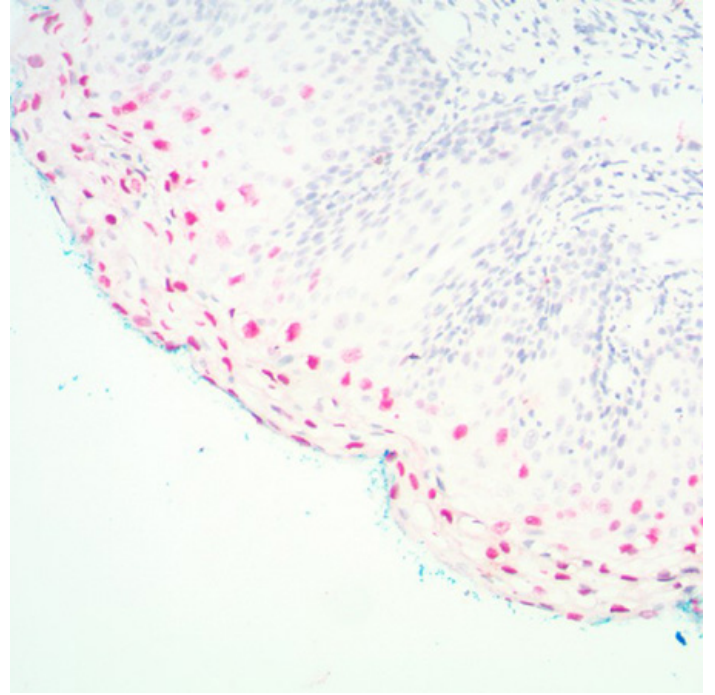
ENZ-32882-1000	1ml
ENZ-32882-6000	6ml

Manuals, SDS & CofA

[View Online »](#)



Negative detection of HPV 16/18/31/33/51 with ENZ-32882 PATHO-GENE® HPV 16/18/31/33/51 probes in vulva FFPE tissue.



Positive detection of HPV 16/18/31/33/51 with ENZ-32882 PATHO-GENE® HPV 16/18/31/33/51 probes in HPV infected cervical FFPE tissue.

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored as recommended. Following dilution in appropriate buffer system, product is stable for at least 2 days.
Handling	Allow contents to warm up to room temperature prior to use.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes

The PATHO-GENE[®] HPV type 16/18/31/33/51 probe employs a mixture of the HPV-16, HPV -18, HPV-31, HPV-33 and HPV-51 specific probes that detect HPV types that have been associated with either condyloma or cervical intraepithelial neoplasia (CIN) and carcinoma in situ (CIS). The probe mixture 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

HPV probe can be optimized for polymer detection system using diluent buffer, *In situ* hybridization buffer for HPV probes (Ready-to-Use) (Prod. No. ENZ-33905).

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.

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)