

PATHO-GENE[®] HPV

type 6/11 probe

A mixture of biotin-labeled HPV 6 and HPV 11-specific probes for *in situ* hybridization

The PATHO-GENE[®] HPV type 6/11 probe is a mixture of biotin-labeled HPV 6 and HPV 11-specific probes in buffered formamide and hybridization enhancers.

Citations: 40

[View Online »](#)

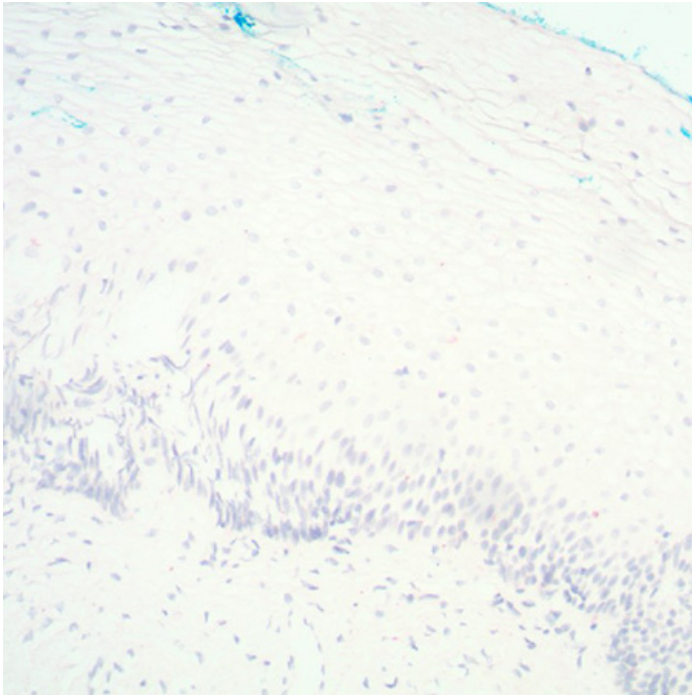
Ordering Information

[Order Online »](#)

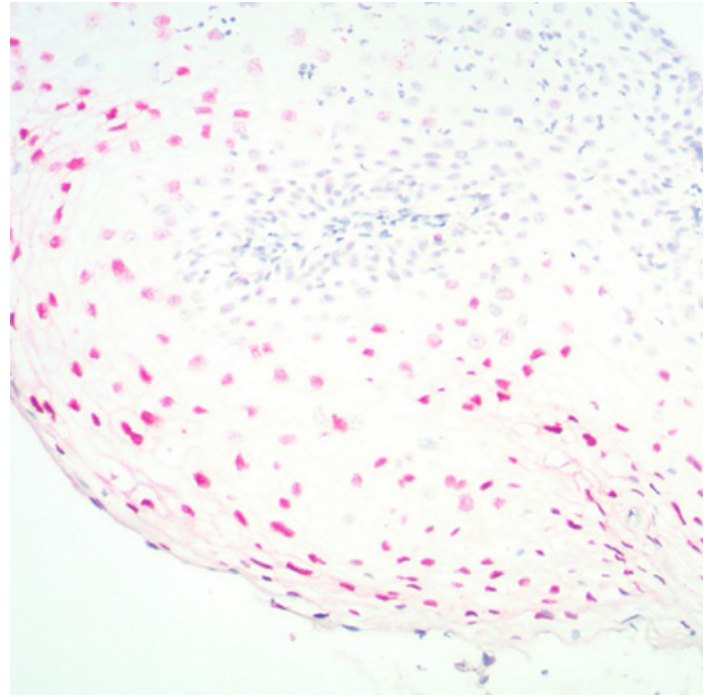
ENZ-32885	1ml
ENZ-32885-6000	6ml

Manuals, SDS & CofA

[View Online »](#)



Negative detection of HPV 6/11 with ENZ-32885 PATHO-GENE® HPV 6/11 probes in cervical FFPE tissue.



Positive detection of HPV 6/11 (red) with ENZ-32885 PATHO-GENE® HPV 6/11 probes in vulva 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.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	The ENZO PATHO-GENE [®] HPV type 6/11 probe employs a mixture of the HPV 6 and HPV 11-specific probes that detect HPV types that have been associated with benign lesions presenting as condylomas and mild dysplasia. 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	<p>HPV probe can be optimized for polymer detection system using diluent buffer, <i>In situ</i> hybridization buffer for HPV probes (Ready-to Use) (Prod. No. ENZ-33905).</p> <p>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.</p>
--------------------------------	--



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)