

# DIGX<sup>®</sup> HPV type 6/11 probe

A ready-to-use digoxigenin-labeled HPV type 6/11 probe

The DIGX<sup>®</sup> HPV type 6/11 probe is a mixture of digoxigenin-labeled HPV type 6 and HPV type 11 specific probes in buffered formamide and hybridization enhancers.

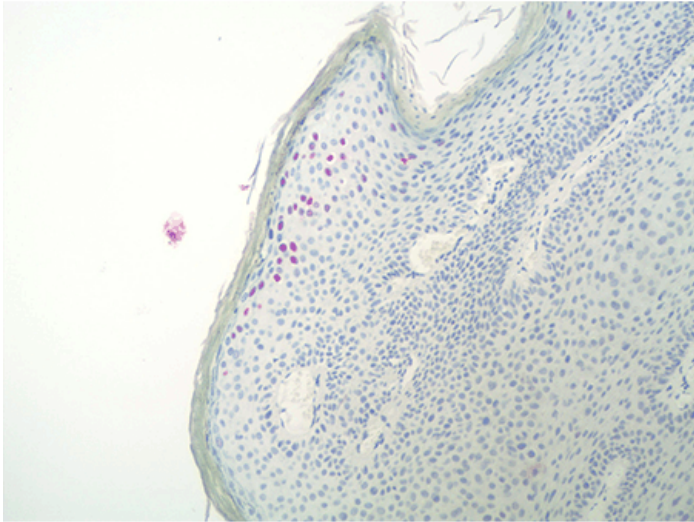
## Ordering Information

[Order Online »](#)

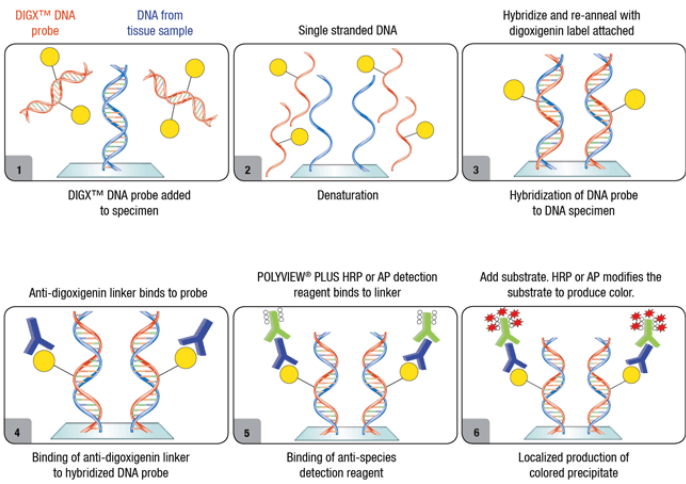
ENZ-GEN112-1000	1ml
ENZ-GEN112-6000	6ml

## Manuals, SDS & CofA

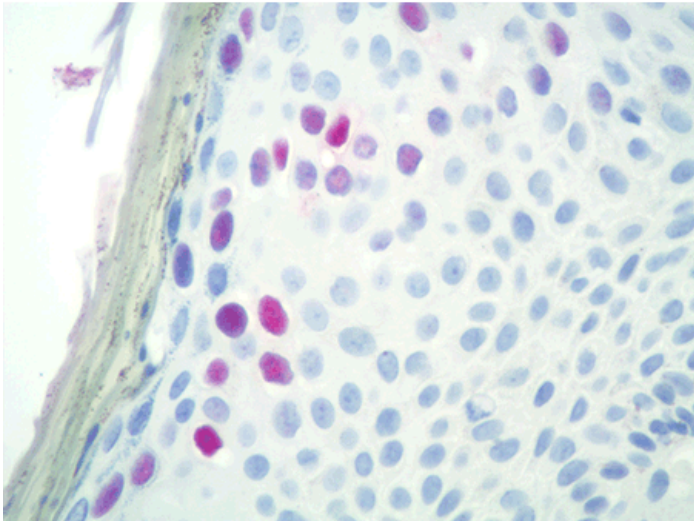
[View Online »](#)



DIGX<sup>®</sup> HPV type 6/11 probe was used to assay cervical tissue using a Leica Bond III. Image was acquired with 10x objective.



Workflow of hybridization with DIGX<sup>®</sup> HPV Probes and detection with DIGX<sup>®</sup> linkers and POLYVIEW<sup>®</sup> PLUS reagents



DIGX<sup>®</sup> HPV type 6/11 probe was used to assay cervical tissue using a Leica Bond III. Image was acquired with 40x objective.

## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended. This product is ready-to-use and additional dilution is not recommended.
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 DIGX<sup>®</sup> HPV type 6/11 probe employs a mixture of the HPV type 6 and HPV type 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 digoxigenin.

### Technical Info / Product Notes

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

DIGX<sup>®</sup> rabbit anti-digoxigenin linker (Prod. No. ENZ-ABS303) or DIGX<sup>®</sup> mouse anti-digoxigenin linker (Prod. No. ENZ-ABS304) can be used for the detection of the DIGX<sup>™</sup> HPV probes in conjunction with POLYVIEW<sup>®</sup> PLUS reagents or kits.

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-  
fr@enzolifesciences.com](mailto:info-fr@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)