

DIGX[®] HPV type 16/18/31/33/51 probe

A ready-to-use digoxigenin-labeled HPV
type 16/18/31/33/51 probe

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

Citations: 1

[View Online »](#)

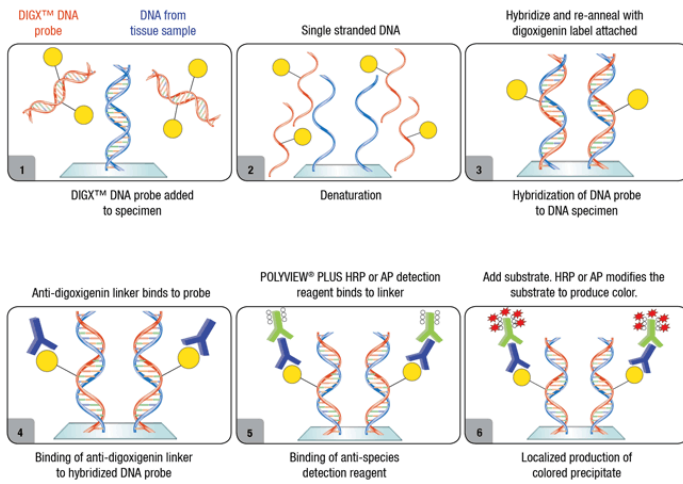
Ordering Information

[Order Online »](#)

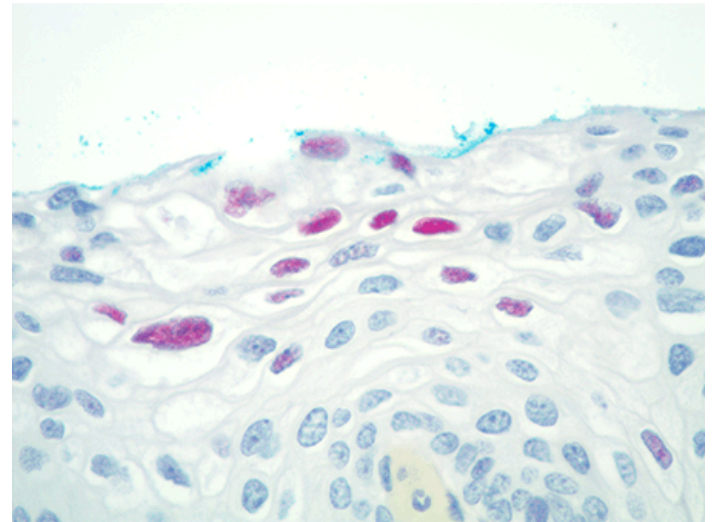
ENZ-GEN114-1000	1ml
ENZ-GEN114-6000	6ml

Manuals, SDS & CofA

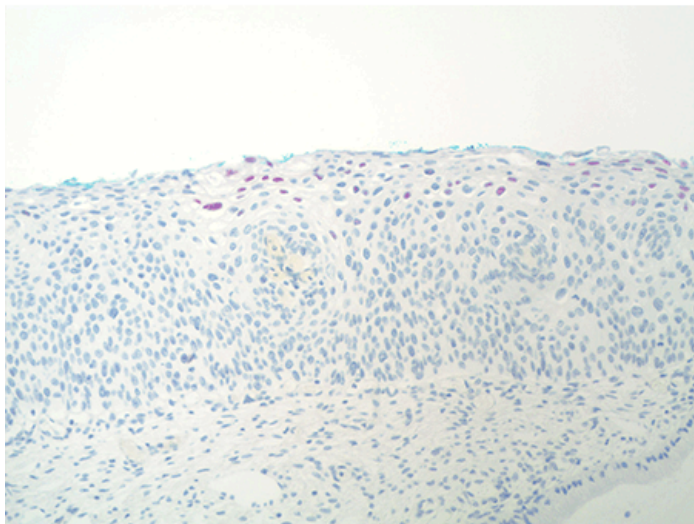
[View Online »](#)



Workflow of hybridization with DIGX[®] HPV Probes and detection with DIGX[®] linkers and POLYVIEW[®] PLUS reagents



DIGX[®] HPV type 16/18/31/33/51 probe was used to assay cervical tissue using a Leica Bond III. Image was acquired with 40x objective.



DIGX[®] HPV type 16/18/31/33/51 probe was used to assay cervical tissue using a Leica Bond III. Image was acquired with 10x 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 [®] HPV type 16/18/31/33/51 probe employs a mixture of the HPV type 16, HPV type 18, HPV type 31, HPV type 33 and HPV type 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 digoxigenin.
-------------------	--

Technical Info / Product Notes	The HPV probe can be optimized for a polymer detection system using diluent buffer, <i>In situ</i> hybridization buffer for HPV probes (Ready-to-Use) (Prod. No. ENZ-33905).
--------------------------------	--

DIGX[®] rabbit anti-digoxigenin linker (Prod. No. ENZ-ABS303) or DIGX[®] mouse anti-digoxigenin linker (Prod. No. ENZ-ABS304) can be used for the detection of the DIGX[®] HPV probes in conjunction with POLYVIEW[®] PLUS reagents and 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.



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)