

DIGX® HPV Type HR-15 Probe

DigX-labeled DNA probe set for ISH detection of 15 high-risk HPV genotypes (types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68, 82)

The DigX HPV HR-15 Probes Set is a ready-to-use in situ hybridization (ISH) probe for chromogenic detection of 15 high-risk human papillomavirus (HPV) genotypes (HPV types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68, and 82) in FFPE tissue sections, fresh tissues, freshly fixed cells, and cell block preparations. The probe is DigX-labeled and compatible with any high-definition chromogen or automated detection system. The signal is brown when HRP/DAB chromogen is used, or red when red AP chromogen is used. All products are Research Use Only (RUO).

Human papillomavirus (HPV) is a group of related DNA viruses. High-risk HPV genotypes are associated with HPV-related conditions of the cervix, oropharynx, and anogenital tract. This probe set simultaneously detects HPV types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68, and 82 in a single hybridization, enabling broad-spectrum identification of active high-risk HPV in tissue specimens. In situ hybridization (ISH) is a molecular technique that uses labeled nucleic acid probes to detect and localize specific DNA sequences directly within tissue sections or cells.

Ordering Information

[Order Online »](#)

ENZ-GEN360-6000	
-----------------	--

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use probe for manual ISH protocols
- Detects 15 high-risk HPV genotypes in a single hybridization: types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68, 82
- Compatible with FFPE tissue sections, fresh tissues, freshly fixed cells, and cell blocks
- Works with any high-definition chromogen or automated detection system
- DigX-labeled — use with anti-DigX antibody-based detection

Handling & Storage

Use/Stability	Stable under recommended storage conditions up to the expiration date indicated on the vial label. Avoid freeze-thaw cycles.
Handling	Protect from light and humidity. Avoid freeze-thaw cycles. Allow all components to reach room temperature before use.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ISH
Recommendation Dilutions/Conditions	Ready to use solution for manual use; Optimization may be required for automated staining systems.
Specificity	HPV high-risk genotypes (15-type panel): types 16, 18, 31, 33, 35, 39, 45, 51, 52, 56, 58, 59, 66, 68, 82
Technical Info / Product Notes	Ready-to-use probe for manual staining. Prior optimization required for automated staining systems. Compatible with Ventana BenchMark and Leica BOND platforms.



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)