

In situ hybridization buffer for HPV probes (Ready-to-Use)

Optimized buffer to use with Enzo's HPV collection of probes

In situ hybridization buffer for HPV probes (Ready-to Use) provides a high quality hybridization medium for *In situ* hybridization analyses using either DNA or oligonucleotide probes labeled with biotin-, or fluorescein-modified nucleotides.

Ordering Information

[Order Online »](#)

ENZ-33905-0050	50ml
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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

Technical Info / Product Notes	Diluent buffer used for optimizing Enzo's HPV line of probes for polymer detection application. These include HPV 16/18/31/33/51 probe (Prod. No. ENZ-32882), HPV Screening probe (Prod. No. ENZ-32884), HPV 6/11 probe (Prod. No. ENZ-32885), HPV 16/18 probe (Prod. No. ENZ-32886) and HPV 31/33/51 probe (Prod. No. ENZ-32887).
--------------------------------	--



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com