

VisionArray® HPV PreCise Master Mix

For the amplification of HPV specific
sequences

The VisionArray HPV PreCise Master Mix is a ready to use PCR Master Mix for amplifying the genomes of HPV types including but not limited to those detected by the VisionArray HPV Chips. The VisionArray PreCise Master Mix contains the VisionArray HPV Primer Mix 2.0, dNTP/dUTP Solution, the VisionArray PreCise Taq DNA Polymerase, the PCR-Buffer, MgCl₂, and the VisionArray Uracil-DNA Glycosylase. During PCR, specific sections of the L1 region of the human papillomavirus (HPV) will be biotinylated. The hybridization and detection of the amplified sequences is subsequently performed using the VisionArray Detection Kit and a VisionArray HPV Chip.

Ordering Information

[Order Online »](#)

ZTV-ES-0007-50	50 tests
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product Notes For the Manual please [click here](#).



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