

# VisionArray®

## Detection Kit

For qualitative detection of DNA sequences on VisionArray Chips.

The VisionArray Detection Kit is designed for the qualitative detection of specific DNA sequences produced with the VisionArray PreCise Master Mix, and the corresponding VisionArray HPV Chip. The VisionArray Detection Kit contains all reagents that are required for the simple and fast hybridization and detection of PCR products on the VisionArray HPV Chips.

### Ordering Information

[Order Online »](#)

ZTV-VK-0003-50	50 tests
----------------	----------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Contents** Hybridization Solution, 100x Wash Buffer, Detection Solution, Blue Spot Solution, Instructions for use

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



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)