

# ZytoFast® HPV type 31/33 Probe (Digoxigenin)

## For the detection of human Human Papilloma Virus (HPV) type 31/33 DNA by chromogenic

**In situ hybridization (CISH)**  
The ZytoFast HPV-CISH System is designed for the detection and discrimination of human papilloma virus (HPV) DNA in paraffin-embedded tissue sections or cell samples. Until now, approximately 10–15 HPV types are associated with lesions that can progress to cancer. The infection with the HPV hr-types can lead to development of cancer of the cervix, vulva, vagina, anus, or penis.

### Ordering Information

[Order Online »](#)

ZTV-T-1057-400	400µl
----------------	-------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Protect from light.

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Human Papilloma Virus Type 31/33

**Application** CISH

**Formulation** Solution of salt (containing synthetic polynucleotides) in water

**Quantity** Sufficient for 40 tests, using 10 µl probe solution per test.

**Specificity** Human papilloma virus types 31/33 DNA

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).  
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)