

ZytoFast® HPV High-Risk (HR) Types Probe (Digoxigenin)

For the detection of human Human Papilloma Virus (HPV) type

The ZytoFast HPV-CISH System is designed for the detection and discrimination of human papilloma virus (HPV) DNA. Most of the approx. 30 identified genital HPV types, predominantly types 6 and 11, are called (low-risk) types, and may cause mild Pap test abnormalities or genital warts. Until now, approximately 10–15 HPV types are associated with lesions that can progress to cancer. These cancer-associated HPV types are designated as high-risk HPV types. 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-1140-400	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 16/18/31/33/35/39/45/51/52/56/58/59/66/68/82

Application CISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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

Specificity HPV type 16/18/31/33/35/39/45/51/52/56/58/59/66/68/82

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

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