

# ZytoFast® EBV Probe (Digoxigenin)

For the detection of Epstein-Barr-Virus (EBV)  
EBER RNA by chromogenic *in situ*  
hybridization (CISH)

The ZytoFast EBV-CISH System is designed for the detection of Epstein-Barr virus (EBV) RNA in paraffin-embedded tissue sections or cell samples. EBV is a member of the gamma-herpesvirus group and one of the most common viruses in humans. Transmission of EBV requires close, intimate contact with a person excreting the virus in its saliva. EBV is the etiological agent of infectious mononucleosis and has been implicated in the pathogenesis of an increasing number of human malignancies such as Burkitt's lymphoma, nasopharyngeal carcinoma, and polyclonal lymphomas.

## Ordering Information

[Order Online »](#)

ZTV-T-1114-400	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** Epstein-Barr Virus

**Application** CISH

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

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

**Specificity** Epstein-Barr virus RNA

**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)