

ZytoFast® human Ig- kappa/Ig-lambda CISH Kit (Biotin/Digoxigenin)

The ZytoFast Ig kappa/Ig lambda CISH System is designed for the detection of Ig-kappa (κ) and/or Ig-lambda (λ) light chain mRNA in paraffin-embedded tissue sections or cell samples. Determination of kappa-to-lambda ratio is useful to distinguish between neoplastic and reactive lymphoid proliferations. Polyclonal expression of κ or λ light chains is considered to reflect a reactive hyperplasia in contrast to the monoclonal expression in malignant lymphoma, the most common hematologic malignancy encountered in the Western world.

Ordering Information

[Order Online »](#)

| | |
|---------------|----------|
| ZTV-T-1005-40 | 40 tests |
|---------------|----------|

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

Contents Ig-kappa/Ig-lambda Probe (Digoxigenin/Biotin labeled); RNA (+)-Control Probe (Biotin labeled); RNA (-)-Control Probe (Biotin labeled); Pepsin Solution; 20x Wash Buffer TBS; HRP-anti-Digoxigenin; AP-Streptavidin; AEC Solution; NBT/BCIP, good for 40 tests.

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