

ZytoFast® human Ig- kappa/Ig-lambda Permanent CISH Kit

For the simultaneous detection of Ig-kappa (κ) and Ig-lambda (λ) mRNA by chromogenic *In situ* hybridization (CISH)

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 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 malignant lymphoma, the most common hematologic malignancy encountered in the Western world.

Ordering Information

[Order Online »](#)

ZTV-T-1105-40	40 tests
---------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Notes



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)