

ZytoLight® SPEC IRF4,DUSP22 Dual Color Break Apart Probe

The ZytoLight® SPEC IRF4,DUSP22 Dual Color Break Apart Probe is designed to detect translocations involving the DUSP22 and IRF4 genes. Rearrangements of the IRF4/DUSP22 chromosomal region have been detected in various B-cell and T-cell lymphomas. Large B-cell lymphoma (LBC) with IRF4 rearrangement is considered a distinct new provisional entity. Most cases have IG/IRF4 fusions and have a favorable prognosis. Moreover, IRF4 translocation has a high specificity for cutaneous cutaneous anaplastic large cell lymphoma (ALCL) supporting the clinical utility of FISH for IRF4 in the differential diagnosis of T-cell lymphoproliferative disorders.

Ordering Information

[Order Online »](#)

ZTV-Z-2210-50	50µl
ZTV-Z-2210-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dual specificity phosphatase 22, JKAP, Interferon regulatory factor 4, MUM1

Application FISH

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), using 10 µl probe solution per test.

Technical Info / Product Notes 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