

ZytoLight® SPEC IKZF1/CEN 7 Dual Color Probe

For the qualitative detection of human IKZF1 gene deletions and alpha satellites of

Chromosome 7 by Fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC IKZF1/CEN 7 Dual Color Probe is designed for the detection of deletions affecting the IKZF1 gene. The IKZF1 gene is located on 7p12.2 and encodes a zinc finger transcription factor, which is required for normal hematopoietic differentiation and proliferation, particularly in lymphoid lineages. Genomic deletions affecting the IKZF1 gene are found in approximately 15% of pediatric and ~40% of adult B-cell precursor acute lymphoblastic leukemia (B-ALL) cases.

Ordering Information

[Order Online »](#)

ZTV-Z-2304-50	50µl
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application FISH

Quantity Sufficient for 5 tests 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