

# ZytoLight® SPEC CRLF2 Dual Color Break Apart Probe

For the detection of translocations involving  
the CRLF2 gene at Xp22.33 and Yp11.32 by

**Fluorescence *in situ* hybridization (FISH).**  
The ZytoLight® SPEC CRLF2 Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal regions Xp22.33 and Yp11.32 harboring the CRLF2 gene. Approximately 7% of patients with B-cell precursor ALL and more than 50% of B-ALL in children with Down syndrome harbor alterations involving the CRLF2 gene.

## Ordering Information

[Order Online »](#)

ZTV-Z-2201-50	50µl
---------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cytokine receptor-like factor 2, CRL2, TSLPR

**Application** FISH

**Quantity** Sufficient for 5 tests, using 10 µl probe solution per test.

**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)