

ZytoLight® SPEC SPI1 Dual Color Break Apart Probe

For the qualitative detection of translocations involving the human SPI1

The ZytoLight® SPEC SPI1 Dual Color Break Apart Probe is designed to detect rearrangements affecting the chromosomal region 11p11.2 harboring the SPI1 gene. SPI1 is a member of the ETS family of transcription factors and is essential for the normal development of hematopoietic stem cells. SPI1 rearrangement were detected in some pediatric T-ALL cases.

Ordering Information

[Order Online »](#)

ZTV-Z-2291-50	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