

ZytoLight® SPEC JAK2 Dual Color Break Apart Probe

For the qualitative detection of translocations involving the human JAK2

The ZytoLight® SPEC JAK2 Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal region 9p24.1 harboring the JAK2 gene. The JAK family proteins are cytoplasmic tyrosine kinases that are essential in maintaining normal hematopoiesis due to their involvement in the JAK-STAT signaling pathway. Rearrangements which lead to constitutive activation of JAK2 kinase have been described in various hematologic malignancies.

Ordering Information

[Order Online »](#)

ZTV-Z-2294-50	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