

ZytoLight® SPEC ABL1 Dual Color Break Apart Probe

For the detection of translocations involving
the ABL1 gene at 9q34.12 by fluorescence *in*

situ hybridization (FISH).

The ZytoLight® SPEC ABL1 Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal region 9q34.12 harboring the ABL1 gene. The translocation t(9;22)(q34.1;q11.2) results in BCR/ABL1 fusion and is observed in approx. 90% of patients with chronic myeloid leukemia and in approx. 25% of adults with acute lymphoblastic leukemia.

Ordering Information

[Order Online »](#)

ZTV-Z-2199-50	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 ABL proto-oncogene 1

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

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