

ZytoLight® SPEC ETV6 Dual Color Break Apart Probe

For the detection of translocations involving
the ETV6 gene at 12p13.2 by fluorescence *in*

situ hybridization (FISH).

The ZytoLight® SPEC ETV6 Dual Color Break Apart Probe is designed for the detection of translocations involving the chromosomal region 12p13.2 harboring the ETV6 gene. The balanced chromosomal translocation t(12;21)(p13.2;q22.1), which leads to ETV6-RUNX1 fusion, represents the most frequent genetic rearrangement (19-27%) in initial childhood B-cell precursor (BCP) acute lymphoblastic leukemia (ALL) and has been associated with good prognosis.

Ordering Information [Order Online »](#)

ZTV-Z-2176-50	50µl
ZTV-Z-2176-200	200µl

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 ETS variant gene 6, TEL

Application FISH

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), 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