

ZytoLight® SPEC USP6 Dual Color Break Apart Probe

For the detection of translocations involving the USP6 gene at 17p13.2 by fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC USP6 Dual Color Break Apart Probe is designed to detect translocations involving the USP6 gene. Translocations affecting USP6 have been initially found in primary aneurysmal bone cysts (ABC), a benign, but locally aggressive bone lesion. More recently, nodular fasciitis (NF), another mesenchymal lesion, has been tested positive for USP6 rearrangements. For both lesions it is assumed that the detection of USP6 rearrangements by FISH represents a valuable diagnostic tool.

Ordering Information [Order Online »](#)

ZTV-Z-2151-50	50µl
ZTV-Z-2151-200	200µl

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ubiquitin-specific peptidase 6, TRE2, TRE17

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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