

ZytoDot® 2C SPEC USP6 Break Apart Probe

For the qualitative detection of translocations involving the human USP6

The ZytoDot® 2C SPEC USP6 Break Apart Probe is designed to detect translocations involving the chromosomal region 17p13.2 harboring the USP6 gene. Translocations affecting USP6 have been initially found in primary aneurysmal bone cysts, a benign, but locally aggressive bone lesion. USP6 rearrangements are restricted to spindle cells in primary ABC, indistinguishable from surrounding normal spindle cells.

Ordering Information

[Order Online »](#)

ZTV-C-3077-100	100µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application CISH

Quantity Sufficient for 10 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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)