

ZytoDot® 2C SPEC FOXO1 Break Apart Probe

For the detection of translocations involving
the FOXO1 gene at 13q14.11 by chromogenic

In situ hybridization (CISH)
The ZytoDot® 2C SPEC FOXO1 Break Apart Probe is designed for the detection of specific translocations involving the chromosomal region 13q14.11 harboring the FOXO1 gene characteristic for alveolar rhabdomyosarcoma. Among solid tumors of the childhood, rhabdomyosarcoma is the most common soft tissue sarcoma.

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| ZTV-C-3065-100 | 100µl |
|----------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Forkhead Box O1, FKHR

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

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