

ZytoLight® SPEC FGFR3 Dual Color Break Apart Probe

For the detection of translocations involving the FGFR3 gene at 4p16.3 by fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC FGFR3 Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal region 4p16.3 harboring the FGFR3 gene. Rearrangements affecting the FGFR3 gene are frequently found in carcinomas of various types including multiple myeloma, bladder cancer, glioblastoma, peripheral T-cell lymphoma, and lung squamous cell carcinoma. Several *in vivo* and *in vitro* studies have demonstrated the therapeutic potential of FGFR inhibitors in cell lines and animal models harboring FGFR3 fusion genes.

Ordering Information

[Order Online »](#)

ZTV-Z-2170-50	50µl
ZTV-Z-2170-200	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 Fibroblast Growth Factor Receptor 3, FGF Receptor 3, JTK4

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