

ZytoLight® SPEC FGFR1/CEN 8 Dual Color Probe

For the detection of the human FGFR1 gene
and alpha-satellites of chromosome 8 by

Fluorescence *in situ* hybridization (FISH). The ZytoLight® SPEC FGFR1/CEN 8 Dual Color Probe is designed for the detection of FGFR1 gene amplification frequently observed in malignant tumors e.g. breast and prostate cancer and oral squamous cell carcinoma (OSCC).

Ordering Information

[Order Online »](#)

ZTV-Z-2072-50	50µl
ZTV-Z-2072-200	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 FGF Receptor 1, Fibroblast Growth Factor Receptor 1, Centromere 8

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