

# ZytoLight® SPEC FGFR2 Dual Color Break Apart Probe

For the detection of translocations involving  
the FGFR2 gene at 10q26.13 by fluorescence

*In situ hybridization (FISH).*  
The ZytoLight® SPEC FGFR2 Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal region 10q26.13 harboring the FGFR2 gene. Translocations and inversions affecting FGFR2 have been detected in several solid tumors, including e.g. breast cancer, lung cancer, and the intrahepatic subtype of cholangiocarcinoma.

## Ordering Information

[Order Online »](#)

ZTV-Z-2169-50	50µl
ZTV-Z-2169-200	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 2, FGF Receptor 2, BEK

**Application** FISH

**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](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)