

ZytoLight® SPEC PDGFRB Dual Color Break Apart Probe

For the detection of translocations involving
the PDGFRB gene at 11q23.3 by

Fluorescence *in situ* hybridization (FISH).
The ZytoLight® SPEC PDGFRB Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 5q32 harboring the PDGFRB gene. Translocations involving the PDGFRB gene are rare genetic disorders and are identified in myelodysplastic/myeloproliferative neoplasms, chronic myeloproliferative disorders, acute myeloid leukemia, and also in atypical chronic myeloid leukemia/chronic myelomonocytic leukemia-like diseases, often with eosinophilia and splenomegaly.

Ordering Information

[Order Online »](#)

ZTV-Z-2197-50	50µl
---------------	------

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 Platelet-derived growth factor receptor-β

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