

ZytoLight® SPEC BRAF Dual Color Break Apart Probe

For the detection of translocations involving
the BRAF gene at 7q34 by fluorescence *in*

situ hybridization (FISH)

The ZytoLight® SPEC BRAF Dual Color Break Apart Probe is designed for the detection of rearrangements involving the chromosomal region 7q34 harboring the BRAF gene. Various BRAF translocations were observed in melanocytic nevi, pilocytic astrocytomas, malignant melanoma, prostate and gastric cancer.

Ordering Information

[Order Online »](#)

ZTV-Z-2189-200	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 B-Raf proto-oncogene, BRAF1, NS7

Application FISH

Quantity Sufficient for 20 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