

ZytoLight® SPEC YWHAE Dual Color Break Apart Probe

For the detection of translocations involving
the YWHAE gene at 17p13.3 by fluorescence

In situ hybridization (FISH)
The ZytoLight® SPEC YWHAE Dual Color Break Apart Probe is designed to detect rearrangements involving the chromosomal region 17p13.3 harboring the YWHAE gene. Altered expression of 14-3-3 family proteins is associated with development and progression of cancer. The fusion between YWHAE and one of the FAM22 family members (FAM22A or FAM22B) caused by a t(10;17)(q22;p13) has been identified in the clinically aggressive, high-grade endometrial stromal sarcoma (ESS) as well as in clear cell sarcoma of the kidney (CCSK).

Ordering Information

[Order Online »](#)

ZTV-Z-2175-50	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 14-3-3ε

Application FISH

Quantity Sufficient for 5 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