

ZytoLight® SPEC CREBBP Dual Color Break Apart Probe

For the qualitative detection of
translocations involving the human CREBBP

gene at 16p13.3 by fluorescence in situ hybridization (FISH)
The ZytoLight® SPEC CREBBP Dual Color Break Apart Probe is designed for the detection of translocations involving the chromosomal region 16p13.3 harboring the CREBBP (a.k.a. CBP) gene. Rearrangements of the CREBBP gene have been observed in several hematologic malignancies, including therapy-related and de novo acute myeloid (AML) or lymphoid leukemia (ALL) and myelodysplastic syndrome (MDS).

Ordering Information

[Order Online »](#)

ZTV-Z-2267-50

50µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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