

ZytoLight® SPEC CIC Dual Color Break Apart Probe

For the qualitative detection of translocations involving the human CIC gene

The ZytoLight® SPEC CIC Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 19q13.2 harboring the CIC gene. Rearrangements involving the CIC gene are frequently found in EWSR1-negative small blue round cell tumors (SBRCT) which arise in soft tissues of children and young adults and have been described as aggressive tumors with an inferior overall survival compared to Ewing sarcoma (EWS).

Ordering Information

[Order Online »](#)

| | |
|---------------|------|
| ZTV-Z-2285-50 | 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