

ZytoLight® SPEC EWSR1 Dual Color Break Apart Probe

For the detection of translocations involving
the EWSR1 gene at 22q12.2 by fluorescence

In situ hybridization (FISH).
The ZytoLight® SPEC EWSR1 Dual Color Break Apart Probe is designed to detect translocations involving the chromosomal region 22q12.2 harboring the EWSR1 gene. Ewing sarcoma is the second most common, highly malignant bone tumor in children and young adults. For prognosis and appropriate treatment it is important to differentiate Ewing sarcoma/PNET from classic neuroblastoma, Wilms tumor, and rhabdomyosarcoma.

Ordering Information

[Order Online »](#)

ZTV-Z-2096-50	50µl
ZTV-Z-2096-200	200µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name EWS, Ewing sarcoma breakpoint region 1

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), 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