

# ZytoLight® SPEC

## EWSR1/FLI1 TriCheck

### Probe

#### For the detection of EWSR-FLI1 rearrangements by fluorescence *in situ* hybridization (FISH)

The ZytoLight® SPEC EWSR1/FLI1 TriCheck™ Probe is designed to detect translocations involving the EWSR1 gene and the FLI1 gene. t(11;22)(q24.3;q12.2) is the most frequent translocation in Ewing sarcoma or peripheral primitive neuroectodermal tumors (PNET) juxtaposing the EWSR1 gene next to the FLI1 locus. Detection of the EWSR1 rearrangements by FISH can be used to confirm the diagnosis of Ewing sarcoma/PNET.

#### Ordering Information

[Order Online »](#)

ZTV-Z-2183-50	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** Ewing Sarcoma breakpoint region 1, EWS, Fli-1 proto-oncogene, ETS transcription factor, EWSR2

**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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)