

ZytoLight® SPEC ERG Dual Color Break Apart Probe

For the detection of translocations involving
the ERG gene at 21q22.2 by fluorescence *in*

situ hybridization (FISH).
The ZytoLight® SPEC ERG Dual Color Break Apart Probe is designed to detect aberrations involving the ERG gene at 21q22.2 frequently detected in prostate cancers. ERG rearrangements have been observed in 40-60% of prostate cancers identified via prostate-specific antigen screening.

Ordering Information

[Order Online »](#)

ZTV-Z-2138-200	200µl
----------------	-------

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 v-ETS Erythroblastosis Virus E26 Oncogene Homolog

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 20 tests, 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