

ZytoLight® SPEC ERG/ TMPRSS2 TriCheck™ Probe

For the detection of **ERG-TMPRSS2 rearrangements by fluorescence *in situ* hybridization (FISH).**

The ZytoLight® SPEC ERG/TMPRSS2 TriCheck™ Probe is designed to detect deletions between the ERG and the TMPRSS2 gene at 21q22 resulting in the TMPRSS2-ERG fusion frequently observed in prostate cancer. Furthermore, the triple color approach allows the detection of other translocations affecting either of these genes.

Ordering Information

[Order Online »](#)

ZTV-Z-2135-200	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/Transmembrane Protease, Serine 2 TriCheck™

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