

ZytoLight® Glioma 1p/19q Probe Set

For the qualitative detection of deletions involving the human chromosomal region 1p36.31 as well as deletions involving the human chromosomal region 19q13.32-q13.33

The ZytoLight Glioma 1p/19q Probe Set includes FISH probes for the detection of both 1p36 and 19q13 loci and the innovative ZyBlack™ Quenching Solution. Deletions affecting the short arm of chromosome 1 (1p36) and the long arm of chromosome 19 (19q13) are frequently found in human gliomas and are associated with a better survival.

Ordering Information

[Order Online »](#)

ZTV-Z-2272-20

20 tests

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 20 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