

ZytoDot® 2C 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 ZytoDot 2C Glioma 1p/19q Probe Set includes CISH probes for the detection of both 1p36 and 19q13 loci. 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. Determination of 1p and 19q status may aid in therapeutic decisions and predict outcome in patients with anaplastic oligodendrogliomas.

Ordering Information

[Order Online »](#)

ZTV-C-3076-10	10 tests
ZTV-C-3076-40	40 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 CISH

Contents ZytoDot 2C SPEC1 p36/1q25 Probe (Prod. No. ZTV-C-3036-100/400)
ZytoDot 2C SPEC 19q13/19p13 Probe (Prod. No. ZTV-C-3037-100/400)

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