

ZytoDot® 2C SPEC MET/CEN 7 Probe

For the detection of the human MET gene and alpha-satellites of chromosome 7 by chromogenic *in situ hybridization (CISH)*.

The ZytoDot® SPEC MET/CEN 7 Probe is designed for the detection of MET gene amplifications found in a variety of human tumors including lung, breast, colorectal, prostate, and gastric carcinomas as well as in gliomas, melanomas and some sarcomas. Several inhibitors of the HGF/MET signaling pathway are being studied and developed as potent therapies to inhibit tumor growth.

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| ZTV-C-3057-400 | 400µl |
|----------------|-------|

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 Met Proto-Oncogene, c-Met, Centromere 7

Application CISH

Quantity Sufficient for 40 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