

ZytoMation®

MET/CEN 7 Dual Color FISH Probe

**For the qualitative detection of human MET
gene amplifications and alpha satellites of**

**Chromosome 7 by fluorescence in situ
hybridization (FISH)**

The ZytoMation MET/CEN 7 Dual Color FISH Probe (PL275) is intended to be used for the qualitative detection of human MET gene amplifications as well as the detection of chromosome 7 alpha satellites in formalin-fixed, paraffin-embedded specimens by fluorescence in situ hybridization (FISH). The probe is intended to be used in combination with the Bond FISH Kit (DS9636) on the automated Bond-MAX or Bond-III system by Leica Biosystems. Interpretation of the results must be made within the context of the patient's clinical history with respect to further clinical and pathologic data of the patient by a qualified pathologist.

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ZTV-Z-2321-5.1ML | 5.1ml |
|------------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

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 FISH

Quantity Sufficient for up to 20 tests using 240 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