

ZytoLight® SPEC PTEN/CEN 10 Dual Color Probe

For the detection of the human PTEN gene and alpha-satellites of chromosome 10 by

Fluorescence *in situ* hybridization (FISH).

The ZytoLight® SPEC PTEN/CEN 10 Dual Color Probe is designed for the detection of PTEN deletions frequently observed in many tumor types, including renal, melanoma, endometrial, breast, prostate, lung, bladder, and thyroid cancer but also in hematological neoplasms. Deletions affecting the long arm of chromosome 10 have been detected in 30 to 50% of early and advanced stage sporadic melanomas and about 40 to 70% of prostate cancers.

Ordering Information

[Order Online »](#)

ZTV-Z-2078-50	50µl
ZTV-Z-2078-200	200µl

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 Phosphatase And Tensin Homolog, MMAC1, Centromere 10

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

Quantity Sufficient for 5 tests (50 µl size) or 20 tests (200 µl size), 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