

ZytoLight® SPEC CDK4/CEN 12 Dual Color Probe

For the detection of the human VHL gene and alpha-satellites of chromosome 12 by

Fluorescence *in situ* hybridization (FISH). The ZytoLight® SPEC CDK4/CEN 12 Dual Color Probe is designed for the detection of CDK4 gene amplifications. Due to amplification of the respective chromosomal region, CDK4 is overexpressed in many human tumors such as soft tissue sarcomas, osteosarcomas, and gliomas. In glioblastomas, the lack of amplification of several genes like CDK4 was recognized to be associated with a longer survival time.

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| ZTV-Z-2103-50 | 50µl |
| ZTV-Z-2103-200 | 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 Cyclin-Dependent Kinase 4/Centromere 12

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