

ZytoLight® SPEC KRAS/CEN 12 Dual Color Probe

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

Fluorescence *in situ* hybridization (FISH). The ZytoLight® SPEC KRAS/CEN 12 Dual Color Probe is designed for the detection of KRAS gene amplifications found e.g. in lung cancer. Mutations of KRAS are frequently found in epithelial malignancies and lead to activation of the downstream mitogen-activated protein kinase resulting in unchecked cellular proliferation and tumor progression.

Ordering Information

[Order Online »](#)

ZTV-Z-2115-200	200µl
----------------	-------

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 v-Ki-ras2 Kirsten rat sarcoma viral oncogene homolog, Centromere 12

Application FISH

Quantity Sufficient for 20 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