

ZytoLight® SPEC VHL/CEN 3 Dual Color Probe

For the detection of the human VHL gene
and alpha-satellites of chromosome 3 by

Fluorescence *in situ* Hybridization (FISH).
The ZytoLight® SPEC VHL/CEN 3 Dual Color Probe is designed for the detection of deletions affecting the VHL gene. Loss of heterozygosity at chromosome 3p and inactivation of the VHL gene by deletion or mutation is the most frequent genetic change in sporadic as well as VHL disease-associated conventional renal cell carcinomas (RCC) whereas alterations of this region are rarely seen in papillary and chromophobe RCC.

Ordering Information

[Order Online »](#)

ZTV-Z-2084-200	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 Von Hippel-Lindau Tumor Suppressor/Centromere 4

Application FISH

Formulation Solution of salt (containing synthetic polynucleotides) in water

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