

ZytoLight® SPEC CCND1 Break Apart/2q11/CEN 6 Quadruple Color

Probe

The ZytoLight® SPEC CCND1 Break Apart/2q11/CEN 6 Quadruple Color Probe is designed for an accurate identification of renal cell carcinoma (RCC) subtypes by the simultaneous detection of rearrangements affecting the CCND1 gene at 11q10.3 and enumeration of chromosomes 2 and 6 in tumor cells. Clear cell RCC, papillary RCC, chromophobe RCC, and renal oncocytomas are the most frequent renal cell tumor subtypes.

For the detection of translocations involving the CCND1 gene at 11q10.3 as well as for the detection of human chromosomes 2 and 6 by fluorescence *in situ* hybridization (FISH).

Ordering Information

[Order Online »](#)

| | |
|----------------|-------|
| ZTV-Z-2118-200 | 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 Cyclin D1, PRAD1, Chromosome 2, Centromere 6

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