

ZytoLight® SPEC CD274, PDCD1LG2/CEN 9 Dual Color Probe

For the detection of the human CD274, PDCD1LG2 gene cluster and the

The ZytoLight® SPEC CD274, PDCD1LG2/CEN9 Dual Color Probe is designed for the detection of CD274, PDCD1LG2 gene cluster amplifications observed in various carcinomas, e.g. classical non-Hodgkin lymphoma and mediastinal large B-cell lymphoma. The genes encode ligands for the PD-1 receptor of T-cells. CD274 is expressed by cancer cells of various tumor types, including melanoma, non-small cell lung cancer, breast cancer, and renal cell carcinomas.

Ordering Information

[Order Online »](#)

ZTV-Z-2179-50	50µl
ZTV-Z-2179-200	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 PDCD1LG1, PDL1, PDL2, CD273

Application FISH

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 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