

ZytoLight® SPEC SMARCB1/22q12 Dual Color Probe

For the detection of the human SMARCB1 gene and chromosome 22 specific

sequences by fluorescence *in situ*
hybridization (FISH)

The ZytoLight® SPEC SMARCB1/22q12 Dual Color Probe is designed for the detection of deletions of the chromosomal region harboring the SMARCB1 gene. Rhabdoid tumors are highly malignant neoplasms that typically arise in infancy and early childhood. They are classified as atypical teratoid/rhabdoid tumors when they occur in the CNS or as malignant rhabdoid tumors when they are found in renal or extra-renal sites.

Ordering Information

[Order Online »](#)

ZTV-Z-2178-50	50µl
---------------	------

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 INI1, SNF5, BAF47

Application FISH

Quantity Sufficient for 5 tests, 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