

PTEN monoclonal antibody (6H2.1)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-2540	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Phosphatase and tensin homolog deleted on chromosome 10

Application ELISA, ICC, IHC, IP, WB

Clone 6H2.1

Formulation Lyophilized. In PBS, pH 7.4.

Host Mouse

Immunogen C-terminal peptide.

Isotype IgG2a kappa

Purity Detail Protein G purified.

Recommendation Dilutions/Conditions If reconstituted with deionized water in 100 µl: WB (1:1'000), IP (1:100), IHC (1:200). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Monkey, Rat

Specificity The antibody recognizes the C-terminal portion of PTEN.

UniProt ID P60484

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: October 22, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)