

# PTEN monoclonal antibody (6H2.1)

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

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-254-C100	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**      If reconstituted with deionized water in 100 µl: WB (1:1'000), IP (1:100), IHC  
**Dilutions/Conditions**      (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](#).



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)