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.

lgG2a 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



info-

eu@enzolifesciences.com