Proteinase inhibitor 9 (human) monoclonal antibody (PI9-17)

The human intracellular serpin proteinase inhibitor 9 (PI9) is the only human protein able to inhibit the activity of the serine protease granzyme B. Granzyme B is expressed by cytotoxic lymphocytes and induces rapid target cell apoptosis.

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

Ordering Information

Order Online »

ALX-804-457-C125

125µg

Manuals, SDS & CofA

View Online »

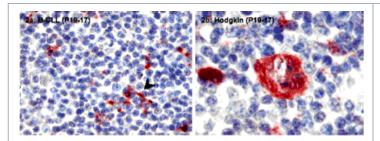


Figure 2: The following tumors were stained for PI9: a) B-CLL and b) classical Hodgkin disease, nodular sclerosing subtype.

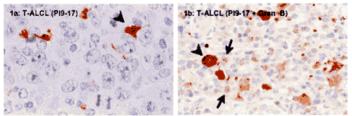


Figure 1: Tissue sections of T-ALCL a) stained with MAb to PI9 (Prod. No. <u>ALX-804-457</u>) and b) double-stained with MAb to PI9 (red) (Prod. No. ALX-804-457) and granzyme B (brown). In panel b the arrows point to tumor-infiltrating CTLs positive for granzyme B and the arrowhead denotes an example of a PI9⁺ neoplastic cell.""""

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PI9, Serpin B9, CAP3, Cytoplasmic antiproteinase 3

Application IHC (PS), WB

Application Notes Detects bands of ~42kDa (native PI9) and ~67kDa (PI9

complexed to granzyme B) by Western blot.

Not suitable for Flow Cytometry.

Clone PI9-17

Crossreactivity Does not cross-react with PI6, PI8 or PAI-2.

Formulation Liquid. In serum free culture supernatant containing 0.7%

BSA and 0.1% sodium azide.

Host Mouse

Immunogen Recombinant human full-length PI9 (proteinase inhibitor

9).

lsotype lgG1

Positive Control YT-Indy cells (a natural killer leukemia cell line) can be

used as positive control. In other cell lines a non-specific

protein band of ~63kDa can be detected.

Recommendation Dilutions/Conditions

Immunohistochemistry (formaldehyde-fixed paraffinembedded human tissues after pretreatment with 10mM citric acid pH 6.0 for 20min; use ~1:100. For fixed cytospin preparations fix cells for 10min in 10% formalin and then treat them as paraffin sections)Western Blot (standard procedures)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human

UniProt ID

P50453

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024

