

# 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 »](#)

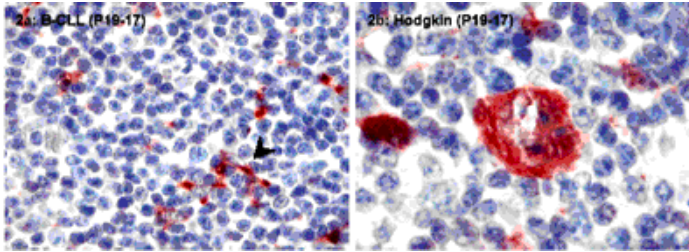
Ordering Information

[Order Online »](#)

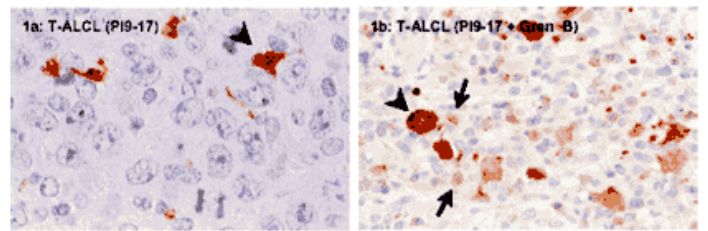
ALX-804-4570	125µg
--------------	-------

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. [ALX-804-457](#)) 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<sup>+</sup> 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).

**Isotype** IgG1

**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 paraffin-embedded 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**

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)