

PTX3 polyclonal antibody

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-365-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

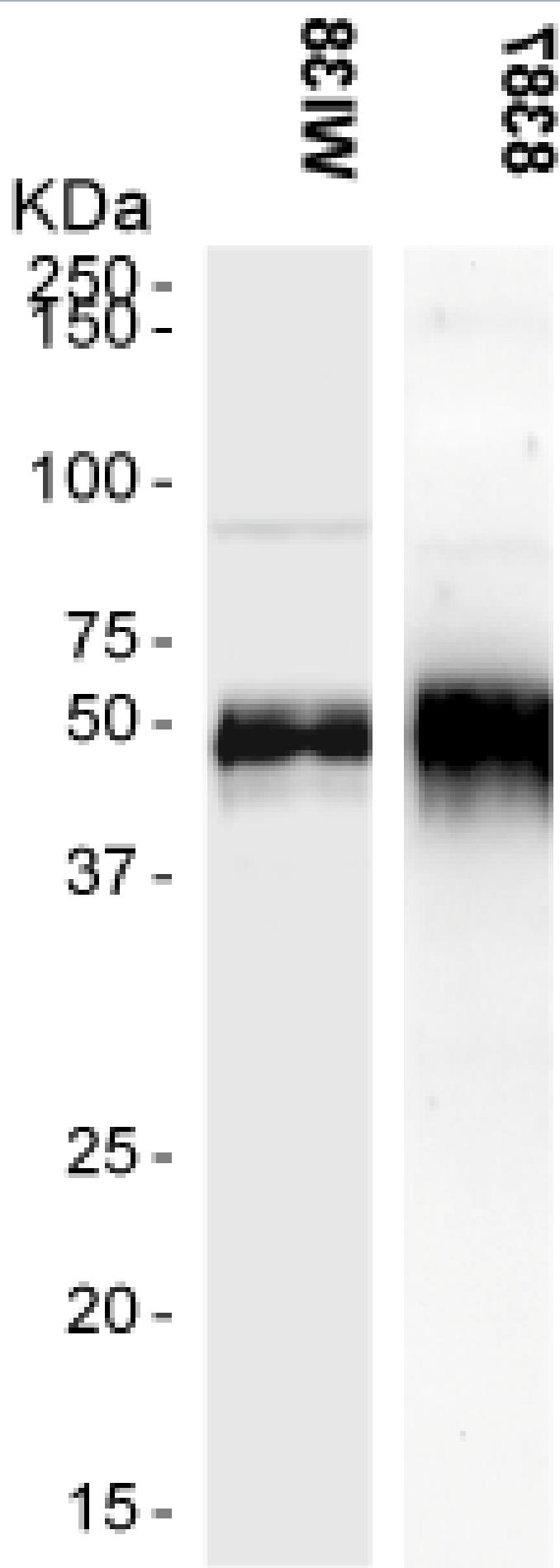


Figure: Lysates of normal lung fibroblasts (WI38) and of human fibrosarcoma (8387) were probed with the PTX3, pAb, (Prod. No. ALX-210-365), over night at +4 °C at a concentration of 500ng/ml in 5% dry milk in PBS+/+ containing 0.05% Tween 20. The signal was revealed with an HRP-conjugated anti-rat IgG secondary antibody in 5% dry milk PBS+/+ containing 0.05%

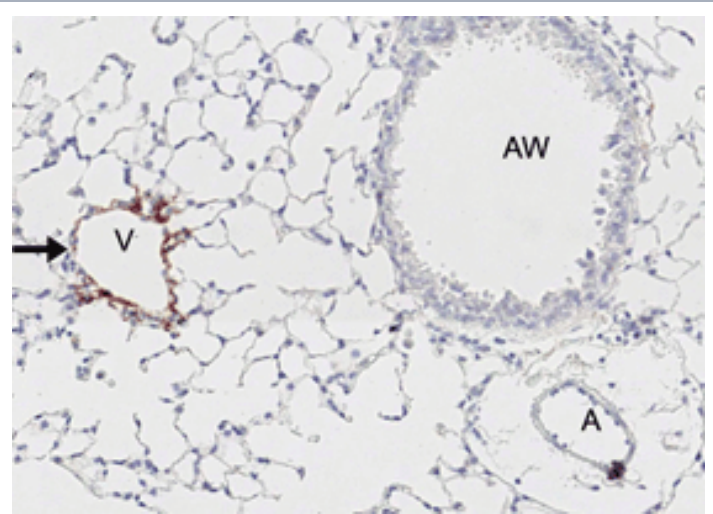


Figure 2: Pulmonary PTX3 expression by immunohistochemistry (IHC) using PTX3, pAb (Prod. No. ALX-210-365).

Method: Photomicrograph of PTX3 stained lung tissue of Cigarette Smoke (CS)-exposed WT mice. Mice were exposed to CS for 4 weeks (subacute).

(V: vein, A: artery, AW: airway. Magnification 200x). In courtesy of Guy Brusselle.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Pentraxin 3, TSG-14, TNF stimulated gene-14

Application ELISA, ICC, IHC (FS), IHC (PS), IP, WB

Application Notes Excellent for use as detection antibody in an ELISA format together with mAb to PTX3 (Prod. No. ALX-804-464 (capture)) and PTX3 standard (Prod. No. ALX-201-124).

Crossreactivity Weakly cross-reacts with bovine PTX3.

Formulation Liquid. In PBS containing 0.1% sodium azide.

Host Rabbit

Immunogen Recombinant human PTX3.

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions ELISA (detection, 1:20,000)Western Blot (500ng/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Chicken, Human, Mouse, Rat

Specificity Recognizes human, mouse and rat PTX3 and chicken homolog PPTX.

UniProt ID P26022

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-be@enzolifesciences.com)

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