

# Peroxisomal membrane protein 70 (mouse) polyclonal antibody

Peroxisomes are single membrane organelles which are involved in many important biochemical pathways and are found in almost all eukaryotic cells. Various toxins including phenols, alcohols, and aldehydes are targeted for modification by peroxisomes. Long chain fatty acid metabolism via  $\beta$ -oxidation is also mediated by this organelle. Loss of peroxisomal functions may result in diseases such as Zellweger syndrome, hyperpipecolic acidemia, neonatal adrenoleukodystrophy, and infantile Refsums disease. Peroxisomal membrane protein 70 (PMP70) is an abundant, integral membrane protein of the peroxisome. This protein is induced by treatment with hypolipidemic agents in parallel with peroxisome proliferation and stimulation of the peroxisomal  $\beta$ -oxidation enzymes.

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

## Ordering Information

[Order Online »](#)

ALX-210-205-C100	100 $\mu$ g
------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## 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** ABCD3, PMP70

**Application** Flow Cytometry, WB

**Application Notes** Detects a band of ~70kDa by Western blot.

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 644-659 (CN644YEFKKITED TVEFGS659) of C-terminal rat PMP70 (peroxisomal membrane protein 70). This sequence is completely conserved in mouse PMP70 and differs from human PMP70 by a single aa substitution.

**Purity Detail** Epitope-affinity purified IgG.

**Recommendation Dilutions/Conditions** Western Blot (1:500) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Mouse

**UniProt ID** P16970

**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-  
fr@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)