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 β -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 β -oxidation enzymes.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-210-205-C100 100μg

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 NotesDetects 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/ConditionsWestern 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

ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 info-

 INC.
 ENZO LIF

 Phone: 800.942.0430
 (ELS) AG

 info Phone: +4

 usa@enzolifesciences.com
 info

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024