

PMCA ATPase monoclonal antibody (5F10)

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

Citations: 3

[View Online »](#)

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 Plasma membrane calcium pump ATPase

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

Application Notes Detects a band of ~140kDa by Western blot. Also detects bands at ~95kDa and ~180kDa which may represent aggregation and/or natural proteolytic products of PMCA ATPase.

Clone 5F10

Formulation Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Purified human erythrocyte PMCA ATPase (plasma membrane calcium pump ATPase).

Isotype IgG2a

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 1:500)Immunohistochemistry (paraffin sections, 1:500)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Amphibian, Cat, Chicken, Dog, Human, Monkey, Mouse, Rabbit, Rat, Sheep

Specificity

Recognizes all known (currently four) isoforms of plasma membrane PMCA ATPase. Recognizes an epitope between aa 724-783 of human PMCA ATPase.

UniProt ID

Q00804 (PMCA1), Q01814 (PMCA2)

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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