PMCA ATPase monoclonal antibody (5F10)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

Ordering Information

Order Online »

ALX-804-049-R100

100µl

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 NotesDetects 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



info-