## Na<sup>+</sup>/K<sup>+</sup> ATPase-α monoclonal antibody (M7-PB-E9)

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

Order Online »

ALX-804-081-1

400µl

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 Sodium/potassium ATPase subunit α

**Application** ELISA, ICC, IHC, IP, WB

**Application Notes**Detects a band of ~110kDa by Western blot.

Clone M7-PB-E9

Formulation Liquid. In 400µl PBS containing 1mg/ml BSA and 0.05%

sodium azide.

**Host** Mouse

**Immunogen** Purified sheep kidney Na+/K+ ATPase-α.

lgG1

Purity Detail Protein A-affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:100)Western Blot

(1:250)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Bovine, Chicken, Dog, Human, Mouse, Porcine, Rat,

Sheep

**Specificity** Recognizes human, mouse, rat (α3 isoform only), sheep,

bovine, pig, dog and chicken Na<sup>+</sup>/K<sup>+</sup> ATPase-α.

UniProt ID P04074 (α1 subunit)

Last modified: May 29, 2024



info-

eu@enzolifesciences.com