## Na+/K+ ATPase α3 subunit monoclonal antibody (XVIF9-G10)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

**Ordering Information** 

Order Online »

BML-SA247-0100

100µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** For long-term storage, aliquot the product into individual tubes and freeze at -20°C.

**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 α3

Application IHC, WB

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

Clone XVIF9-G10

**Formulation** Liquid. Ascites fluid containing 0.05% sodium azide.

**Host** Mouse

**Immunogen** Canine cardiac microsomes.

lgG1

Purity Detail Diluted ascites.

Recommendation Dilutions/Conditions Immunohistochemistry (5µg/ml)Western Blot

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

determined individually for each application.

Species Reactivity Dog, Human, Monkey, Rabbit, Rat, Sheep

**Specificity** Recognizes the  $\alpha$ 3 subunit of Na<sup>+</sup>/K<sup>+</sup> ATPase.

UniProt ID F1P767

Last modified: May 29, 2024



info-

eu@enzolifesciences.com