

Na⁺/K⁺ ATPase α 3 subunit monoclonal antibody (XVIF9-G10)

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

Citations: 4

[View Online »](#)

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 $\alpha 3$
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.
Isotype	IgG1
Purity Detail	Diluted ascites.
Recommendation Dilutions/Conditions	Immunohistochemistry (5 μ g/ml)Western Blot (1 μ g/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 ⁺ /K ⁺ ATPase.
UniProt ID	F1P767

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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