

Na⁺/K⁺ 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	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

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 400 μ l PBS containing 1mg/ml BSA and 0.05% sodium azide.
Host	Mouse
Immunogen	Purified sheep kidney Na ⁺ /K ⁺ ATPase- α .
Isotype	IgG1
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 (α 3 isoform only), sheep, bovine, pig, dog and chicken Na ⁺ /K ⁺ ATPase- α .
UniProt ID	P04074 (α 1 subunit)

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](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)