

Na⁺/K⁺ ATPase-α monoclonal antibody (M7-PB-E9)

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

Citations: 2

[View Online »](#)

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 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)

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)