

Na⁺ /K⁺ ATPase-β1 monoclonal antibody (M17-P5-F11)

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

Ordering Information

[Order Online »](#)

| | |
|-------------------|-------|
| ALX-804-085A-0100 | 100µg |
|-------------------|-------|

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 β 1 |
| Application | ICC, IF, IHC, IHC (PS), IP, WB |
| Clone | M17-P5-F11 |
| Formulation | Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide. |
| Host | Mouse |
| Immunogen | Purified sheep kidney Na ⁺ /K ⁺ ATPase- β . |
| Isotype | IgG2a |
| Purity Detail | Protein A-affinity purified |
| Recommendation Dilutions/Conditions | Immunohistochemistry (1:100)Immunohistochemistry (1:200, paraffin sections)Immunofluorescence (1:100-1:1000)Western Blot (1:1000-1:10,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity | Dog, Human, Mouse, Porcine, Sheep |
| Specificity | Recognizes the β 1 subunit of Na ⁺ /K ⁺ ATPase- β . |
| UniProt ID | P05028 |
| 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)