

Mineralocorticoid (aldosterone) receptor monoclonal antibody (H10E4C9F)

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

Citations: 8

[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 MR

Application FUNC, ICC, IHC (PS), WB

Application Notes Detects a band of ~116kDa by Western blot.
Functional Application: Blocks the binding of aldosterone to the MR.
Not suitable for immunoprecipitation.

Clone H10E4C9F

Formulation Liquid. Purified ascites in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Aldosterone.

Isotype IgG1

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 1:100; staining is blocked by pre-incubating the sample with aldosterone)Western Blot (1:1,000; ECL: Enhanced detection systems like enhanced chemiluminescence are recommended).Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Produced using the anti-idiotypic method.

Species Reactivity

Chicken, Human, Rabbit, Rat

UniProt ID

P08235 (human), P22199 (rat)

Worry-free GuaranteeThis 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)