

# Mineralocorticoid (aldosterone) receptor monoclonal antibody (H10E4C9F)

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-099-R100	100µl
------------------	-------

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 Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

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)