

EDA-A1 monoclonal antibody (Renzo-2)

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

Citations: 3

[View Online »](#)

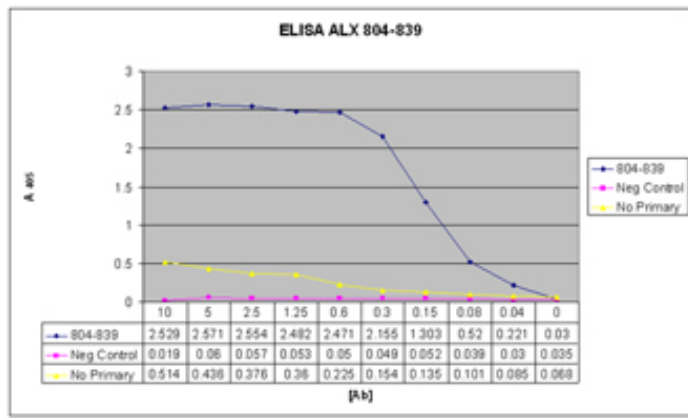
Ordering Information

[Order Online »](#)

ALX-804-839-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



ELISA analysis of the mAb concentration vs Absorbance (A405) of the human EDA-A1 using the mAb (Prod. No. ALX-804-839). Blue, experimental data, substrate EDA-A1(Prod. No. ALX-522-038); Pink, negative control, substrate water; Yellow, EDA-A1 (Prod. No. ALX-522-036) no primary antibody.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ectodysplasin-A1
Application	ELISA, Flow Cytometry, IP
Clone	Renzo-2
Crossreactivity	EDA-A2
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human EDA-A1 (ectodysplasin-A1) extracellular domain (aa 245-391) (Prod. No. ALX-522-038).
Isotype	IgG1
Positive Control	Prod. No. ALX-522-038
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse
UniProt ID	Q92838
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)