

UNC5A (human) monoclonal antibody (Dettwy-1)

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

Ordering Information

[Order Online »](#)

ALX-804-864-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling For long term storage, aliquot and freeze at -20°C. Avoid freeze/thaw cycles.

Short Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Netrin receptor UNC5A precursor, UNC5H1, UNC-5 homolog A, UNC-5 homolog 1

Application ELISA, Flow Cytometry

Clone Dettwy-1

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Mouse

Immunogen Extracellular domain of human UNC5A (aa 26-306) fused to the Fc portion of human IgG1.

Isotype IgG1

Purity Detail Protein G-affinity purified.

Source Purified from concentrated hybridoma tissue cell culture supernatant.

Species Reactivity Human

UniProt ID Q6ZN44

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

.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: May 29, 2024
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com