

Galectin-3 monoclonal antibody (A3A12)

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

Citations: 2

[View Online »](#)

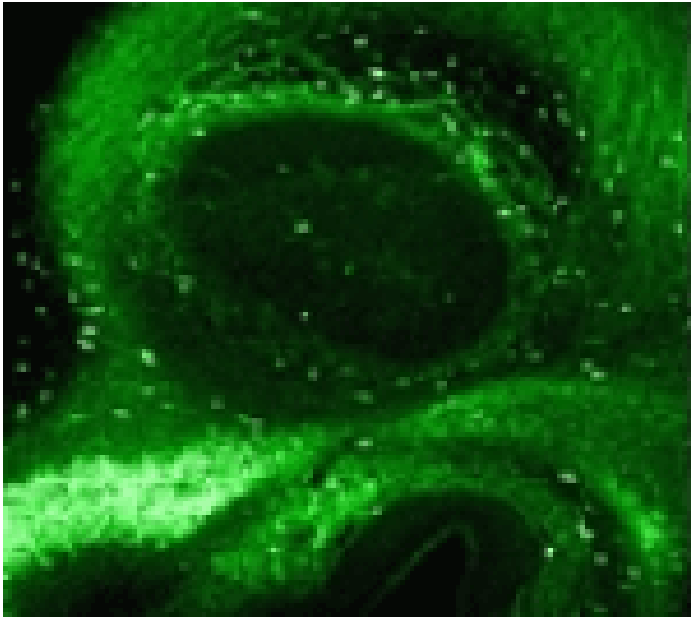
Ordering Information

[Order Online »](#)

ALX-804-284-C100	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

Application ELISA, Flow Cytometry, IHC (PS), WB

Application Notes Detects a band of ~30kDa by Western blot.

Clone A3A12

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Recombinant human galectin-3.

Isotype IgG1

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 6µg/ml)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P47211

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



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-
fr@enzolifesciences.com](mailto:info-fr@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)