## Galectin-3 monoclonal antibody (A3A12)

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

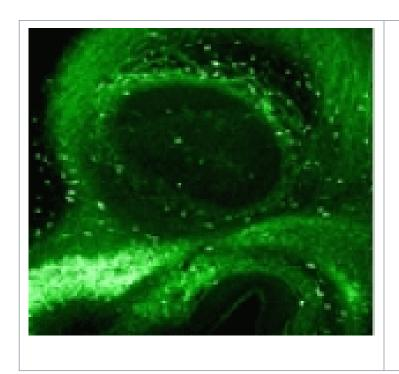
Order Online »

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.

lsotype lgG1

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

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

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