## Furin convertase (R2) monoclonal antibody (MON-148)

This antibody is covered by our Worry-Free Guarantee.

Citations: 14

View Online »

**Ordering Information** 

Order Online »

ALX-803-015-R100

100µl

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** Stable for at least 3 months when stored at +4°C.

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Application ICC, IP, WB

Clone MON-148

**Formulation** Liquid. Ascites fluid. Contains no preservatives.

**Host** Mouse

**Immunogen** Recombinant human furin.

lsotype lgG2a

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:200-1:400)Immunoprecipitation

(1:200; 50mM TRIS/HCl, pH 7.4, 150mM NaCl, 1% Triton X-100; 1% sodium deoxycholate, 0.1% SDS)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Source Ascites fluid from Balb/c mice obtained after ingection of

MON-148 producing hybridoma cells and isolated using a

physiological salt solution (0.9% NaCl).

Species Reactivity Human, Mouse

**Specificity** Recognizes epitope region 2 (R2) in the catalytic domain

of furin convertase.

P09958

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com