

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

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

Citations: 14

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 3 months when stored at +4°C.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	ICC, IP, WB
<b>Clone</b>	MON-148
<b>Formulation</b>	Liquid. Ascites fluid. Contains no preservatives.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human furin.
<b>Isotype</b>	IgG2a
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Source</b>	Ascites fluid from Balb/c mice obtained after ingestion of MON-148 producing hybridoma cells and isolated using a physiological salt solution (0.9% NaCl).
<b>Species Reactivity</b>	Human, Mouse
<b>Specificity</b>	Recognizes epitope region 2 (R2) in the catalytic domain of furin convertase.

UniProt ID

P09958

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)