

Furin convertase (human) monoclonal antibody (MON-152)

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

Citations: 27

[View Online »](#)

Ordering Information

[Order Online »](#)

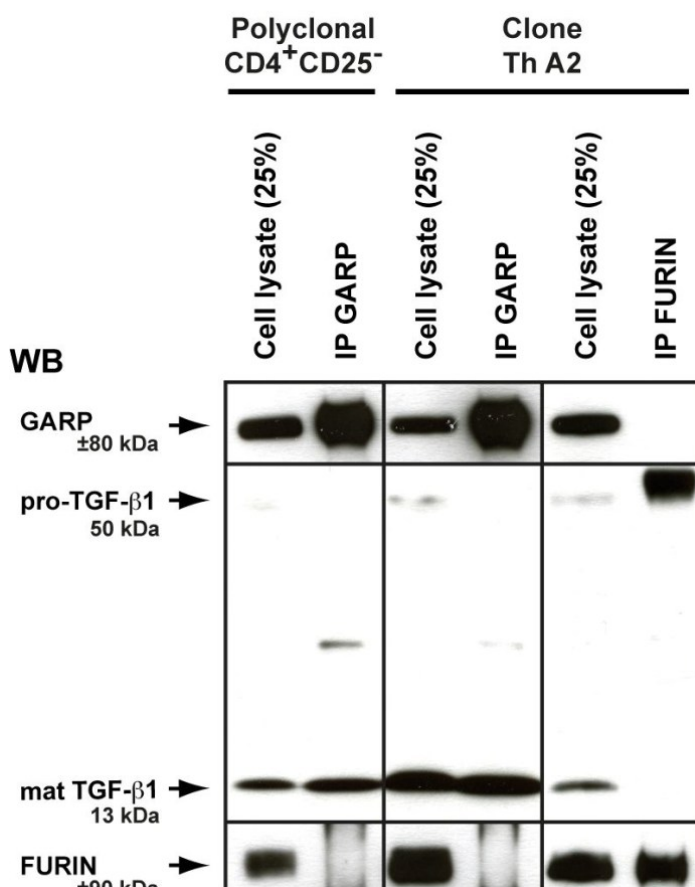
ALX-803-017-R100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

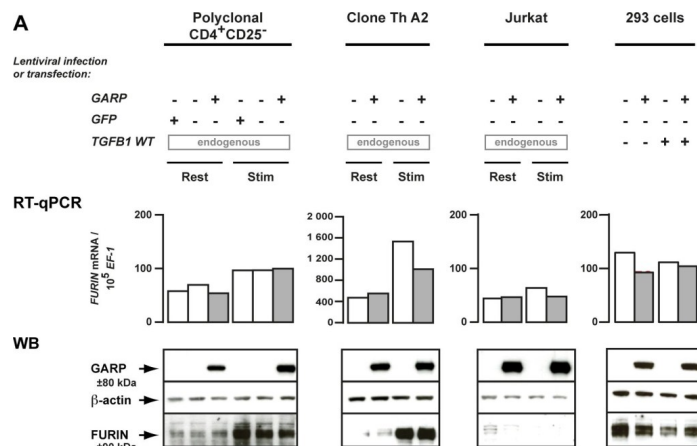
C

GARP-transduced cells:



GARP does not increase FURIN expression or activity, and does not co-immunoprecipitate with FURIN.A. Expression of FURIN mRNA and protein were analyzed by RT-qPCR and WB in the human cells described in Figure 2. B. FURIN activity was measured 24 hours after transfection of 293 cells. Lysates of transfected cells were incubated with a FURIN fluorogenic substrate directly (top panel), or after capture on plastic-coated anti-FURIN antibody (bottom panel), to measure FURIN-like or FURIN specific activity, respectively. Graphs show mean fluorescence intensity at the indicated time (min) after addition of the substrate. The FURIN inhibitor Dec-RVKR-CMK was added to one condition to verify the specificity of the assay. C. Lysates of cells described in Figure 2 were immunoprecipitated with anti-GARP (IP GARP) or anti-FURIN (IP FURIN) antibodies. Immunoprecipitation products or total cell lysates (25% of input used for IPs) were analyzed by WB with anti-GARP, anti-TGF-β or anti-FURIN antibodies, as indicated.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: GARP is regulated by miRNAs and controls latent TGF-β1 production by human regulatory T cells. *PLoS One* (2013)



GARP does not increase FURIN expression or activity, and does not co-immunoprecipitate with FURIN.A. Expression of FURIN mRNA and protein were analyzed by RT-qPCR and WB in the human cells described in Figure 2. B. FURIN activity was measured 24 hours after transfection of 293 cells. Lysates of transfected cells were incubated with a FURIN fluorogenic substrate directly (top panel), or after capture on plastic-coated anti-FURIN antibody (bottom panel), to measure FURIN-like or FURIN specific activity, respectively. Graphs show mean fluorescence intensity at the indicated time (min) after addition of the substrate. The FURIN inhibitor Dec-RVKR-CMK was added to one condition to verify the specificity of the assay. C. Lysates of cells described in Figure 2 were immunoprecipitated with anti-GARP (IP GARP) or anti-FURIN (IP FURIN) antibodies. Immunoprecipitation products or total cell lysates (25% of input used for IPs) were analyzed by WB with anti-GARP, anti-TGF-β or anti-FURIN antibodies, as indicated.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: GARP is regulated by miRNAs and controls latent TGF-β1 production by human regulatory T cells. *PLoS One* (2013)

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-152
Crossreactivity	Does not cross-react with mouse furin convertase.
Formulation	Liquid. Ascites. Contains no preservatives.
Host	Mouse
Immunogen	Recombinant human furin.
Isotype	IgG1
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 as obtained after injection of MON-152 producing hybridoma cells and isolated using a physiological salt solution (0.9% NaCl).
Species Reactivity	Human
Specificity	Recognizes an epitope in the cysteine-rich region of human furin convertase.
UniProt ID	P09958

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-
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)