

TRIF polyclonal antibody

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

Citations: 5

[View Online »](#)

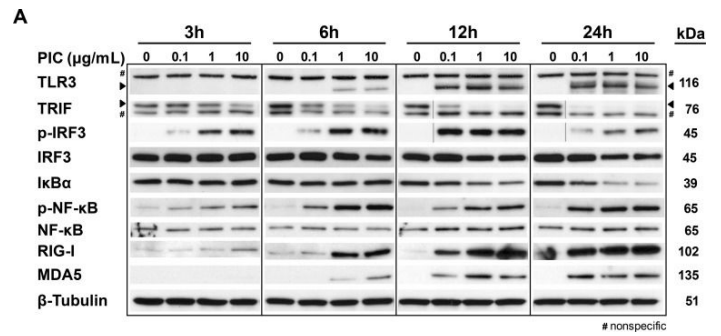
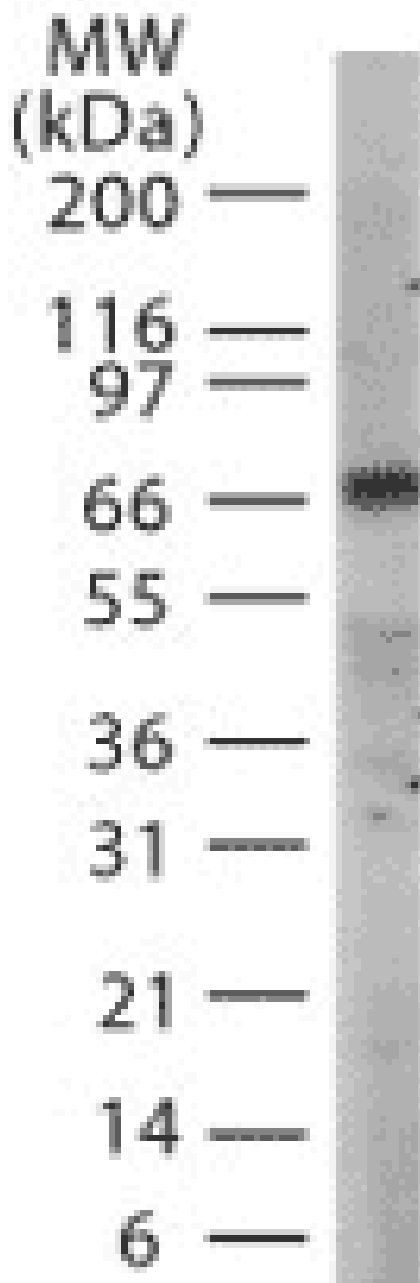
Ordering Information

[Order Online »](#)

ALX-215-0160	200µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)



Poly(I:C) activates the TLR3/TRIF signaling. (A) HLECs were stimulated with Poly(I:C) at 0.1, 1 and 10 $\mu\text{g/mL}$ for 3, 6, 12 and 24 h. Whole cell lysates were analyzed for TLR3, TRIF, phosphorylated-IRF3, total IRF3, I κ B α , phosphorylated-NF- κ B, NF- κ B, RIG-I, MDA5, and β -tubulin (as a loading control) by Western blot. Arrowheads indicate TLR3 (116 kDa, lower band) and TRIF (76 kDa, top band). #, indicates a nonspecific band. Immunoblots are representative of three separate experiments that all showed similar results. (B) Poly(I:C) induces nuclear translocation of NF- κ B. HLEC were stimulated with medium (NT) or 5 $\mu\text{g/mL}$ Poly(I:C) for 24 h. Immunofluorescence staining was performed using antibodies against NF- κ B (red) and nuclei were counterstained with Hoechst 33342 (blue). Images are representative of three separate experiments (400x total magnification).

Image collected and cropped by CiteAb under a CC0-1.0 license from the following publication: Poly(I:C) Induces Human Lung Endothelial Barrier Dysfunction by Disrupting Tight Junction Expression of Claudin-5. *PLoS One* (2016)

Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TICAM-1, Toll-interleukin-1 receptor [TIR] domain-containing adaptor inducing interferon- β [IFN- β], TIR-containing adaptor molecule-1
Application	WB
Formulation	Liquid. Whole antisera with 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 143-158 (A143APGGAIVSELCQALS158) of mouse TRIF [TICAM-1].
Recommendation Dilutions/Conditions	Western Blot (1:500-1:1,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	Q80UF7

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024