

Toll-like receptor 3 monoclonal antibody (40C1285.6)

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

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-362-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

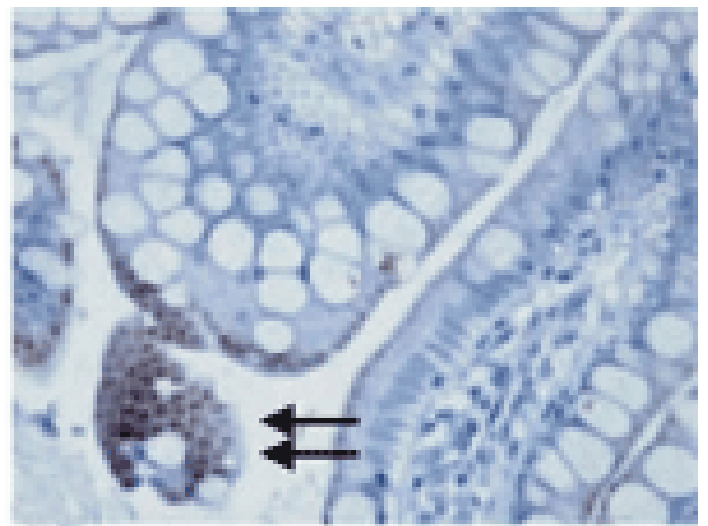
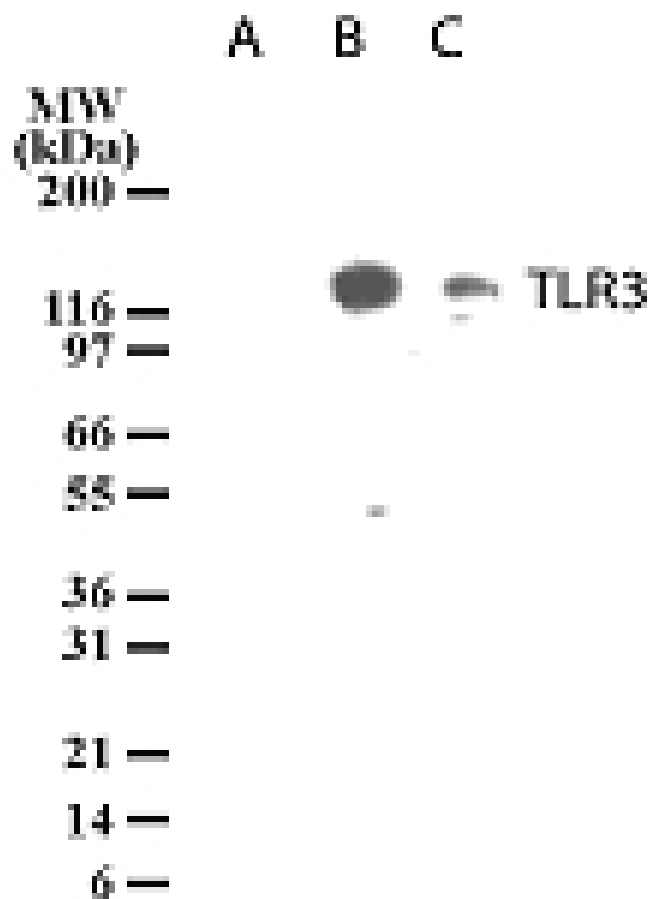


Figure: Western blot analysis of TLR3 in lysates from untransfected 293 cells (lane A), 293 cells transfected with human TLR3 cDNA (lane B), and 20µg/lane human intestine tissue lysate (lane C).

Handling & Storage

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

Alternative Name	TLR3, CD283
Application	Flow Cytometry, ICC, IHC (PS), IP, WB
Clone	40C1285.6
Formulation	Liquid. In PBS containing 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 55-70 (VLNLTHNQLRRLPAAN) of human TLR3 (Toll-like receptor 3) cytoplasmic portion.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-4µg/106 cells for intracellular staining)Immunoprecipitation (2µg/106 transfected cells)Western Blot (1-3µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Dog, Human, Mouse
UniProt ID	O15455

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)