

TLR7 monoclonal antibody (4F4)

TLR7 is one of toll-like receptors which are type I transmembrane proteins involved in innate immunity by recognizing microbial conserved structures. These proteins are characterized by multiple copies of leucine rich repeats in the extracellular domain and IL-1 receptor motif in the cytoplasmic domain, and act through adaptor molecules such as MyD88 and TIRAP to activate various kinases and transcription factors.

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

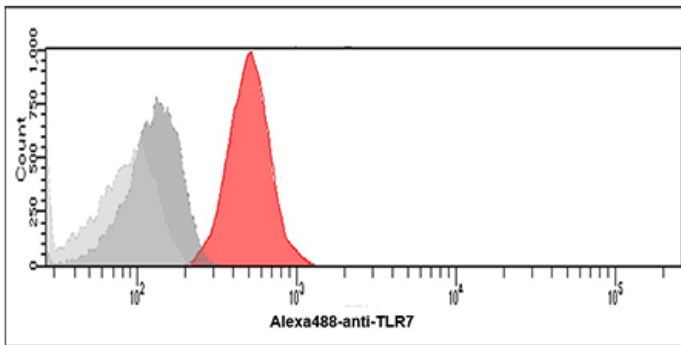
Ordering Information

[Order Online »](#)

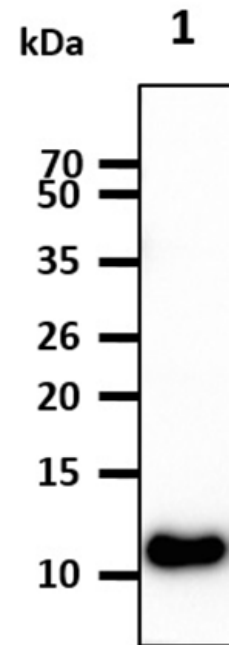
| | |
|-----------------|-------|
| ENZ-ABS742-0100 | 100µl |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of TLR7 in THP-1 cells. The cell was stained at 2-5 μ g for 1 \times 10⁶ cells (red). A Goat anti mouse IgG (Alexa fluor 488) was used as the secondary antibody. Mouse monoclonal IgG was used as the isotype control (dark gray), cells without incubation with primary and secondary antibody was used as the negative control (light gray).



The recombinant protein (50ng) was resolved by SDS-PAGE, transferred to PVDF membrane and probed with anti-human TLR7 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system. Lane 1.: Recombinant human TLR7 protein.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name toll-like receptor 7

Application ELISA, Flow Cytometry, WB

Clone 4F4

Formulation Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.

Host Mouse

Immunogen Recombinant human TLR7 (aa 451-500) purified from E. coli.

Isotype IgG1κ

Purity Detail Protein A affinity purified

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

UniProt ID Q9NYK1

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-
fr@enzolifesciences.com](mailto:info-fr@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)