

Toll-like receptor 5 (human) polyclonal antibody

Note: Human Toll-like Receptor 5 (TLR5) is also described as Toll/Interleukin-1 Receptor-like Gene 3 (TIL3).

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---|
| Handling | For maximum product recovery after thawing, centrifuge the vial before opening the cap. 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 | TLR5, Toll/Interleukin-1 receptor-like gene 3, TIL3 |
| Application | ELISA, Flow Cytometry, IHC, WB |
| Formulation | Affinity purified antibody in 10 mM KHPO ₄ , 140 mM NaCl, BSA @ 1mg/ml and 0.1% sodium azide. |
| Host | Goat |
| Immunogen | Synthetic peptide corresponding to aa 151-181 (D151LSKNQIRSLYLHPSFGKLNLSLKSIDFSSNQ181) of human TLR5 (Toll-like receptor 5). |
| Purity Detail | Epitope-affinity purified IgG. |
| Quality Control | Immunocytochemistry on human peripheral blood leukocytes. |
| Recommendation Dilutions/Conditions | ELISA (1:37,500)Western Blotting (1:50)IHC (1:250)Flow Cytometry (1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity | Human |
| UniProt ID | O60602 |
| 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)