

# IP3 receptor type 1 polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

|                      |  |
|----------------------|--|
| Alternative Name     | IP3R1, ITPR1, IP3R-I, Inositol 1,4,5-triphosphate receptor type 1, INSP3R1 |
| Application          | IHC, WB  |
| Application Notes    | Detects a band of ~240kDa by Western blot.                                 |
| Formulation          | Liquid. Neat serum containing 0.05% sodium azide.                          |
| Host                 | Rabbit   |
| Immunogen            | Synthetic peptide corresponding to aa 1829-1848 of human IP3R1.            |
| Species Reactivity   | Dog, Human, Rat  |
| UniProt ID           | Q14643   |
| Worry-free Guarantee | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>       |

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)