

# Thyroid peroxidase polyclonal antibody

Thyroid peroxidase (TPO) is an enzyme crucial for the synthesis of thyroid hormones. It plays a key role in the iodination of tyrosine residues in thyroglobulin. Anti-TPO antibodies are often used in the diagnosis of autoimmune thyroid diseases, such as Hashimoto's thyroiditis and Graves' disease.

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

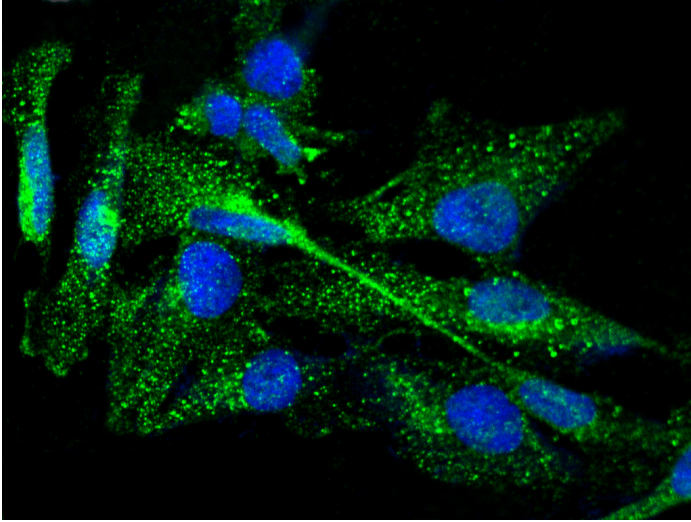
## Ordering Information

[Order Online »](#)

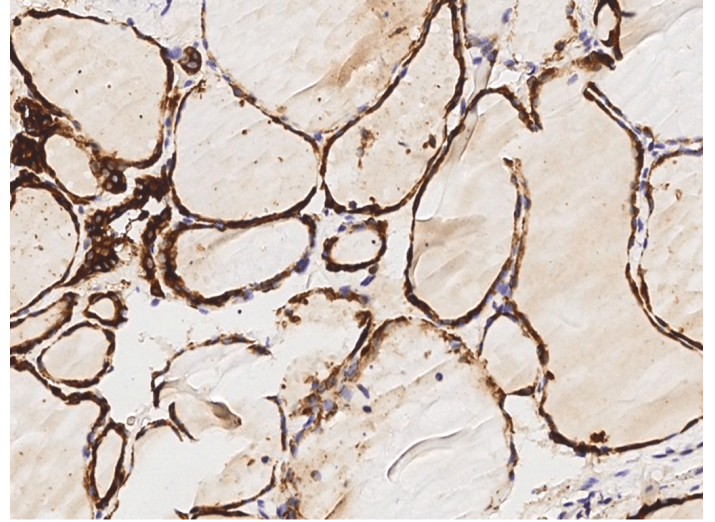
ENZ-ABS1040-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of TPO in U251MG cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human TPO polyclonal antibody (dilution ratio 1:1000) at 4°C overnight. Then cells were stained with the Alexa Fluor 488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to cytoplasm.



Immunochemical staining of human TPO in human thyroid with rabbit polyclonal antibody (1:1000, formalin-fixed paraffin embedded sections).

## 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** TPO

**Application** ELISA, ICC, IF, IHC (PS), WB

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free. Carrier-free.

**Host** Rabbit

**Immunogen** Recombinant human Thyroid peroxidase / TPO.

**Isotype** IgG

**Purity Detail** Protein A and antigen affinity purified.

**Recommendation Dilutions/Conditions** ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:500-1:2'000)Western Blot (1:500-1:2'000)Immunocytochemistry / Immunofluorescence (1:300-1:10'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

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

Last modified: October 7, 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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)