

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](#).

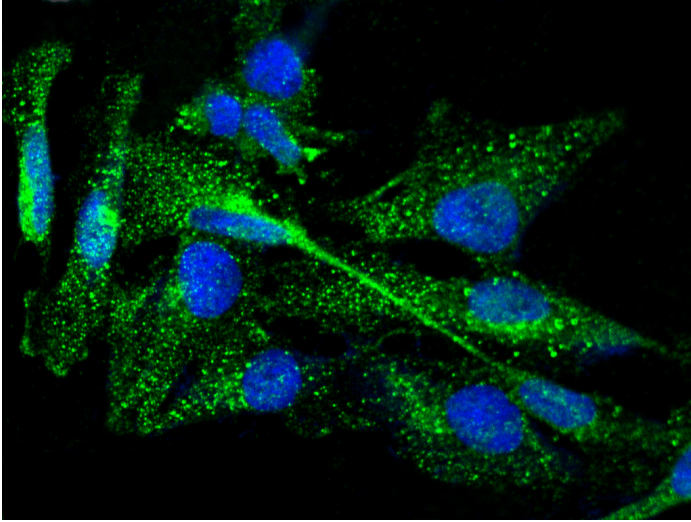
Ordering Information

[Order Online »](#)

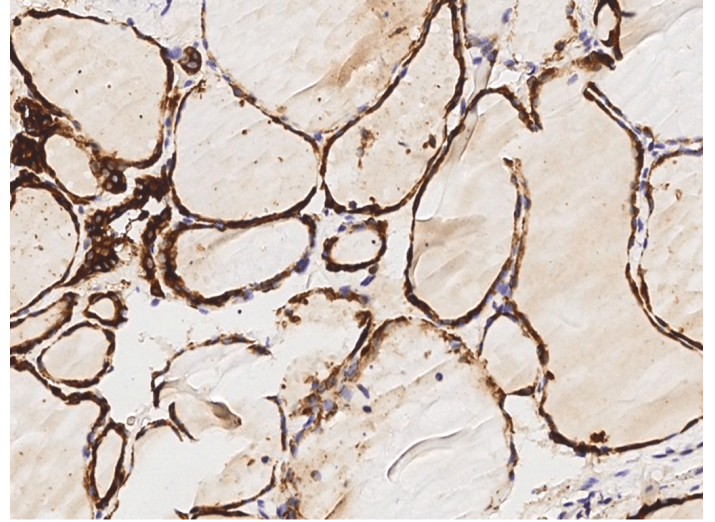
ENZ-ABS1040-0100	100µl
------------------	-------

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 .



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)