

# PRODUCT INFORMATION



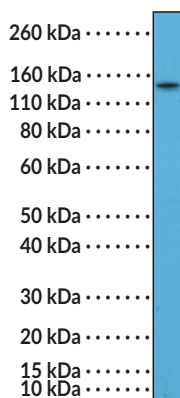
## TPO (C-Term) Rabbit Monoclonal Antibody (Clone RM368)

Item No. 32293

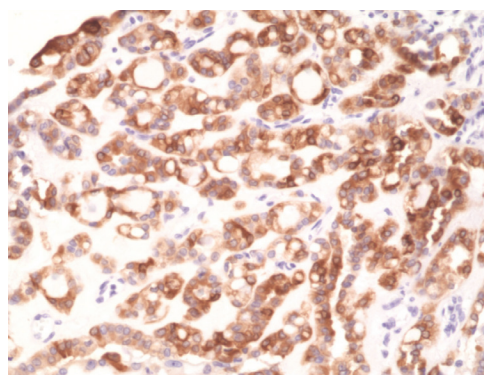
### Overview and Properties

|                            |   |
|----------------------------|---|
| <b>Contents:</b>           | This vial contains 100 µl of protein A-affinity purified monoclonal antibody.   |
| <b>Synonyms:</b>           | Iodide Peroxidase, Thyroid Peroxidase, Thyroperoxidase  |
| <b>Immunogen:</b>          | Peptide from the C-terminal region of human TPO   |
| <b>Cross Reactivity:</b>   | (+) TPO   |
| <b>Species Reactivity:</b> | (+) Human   |
| <b>Form:</b>               | Liquid  |
| <b>Storage:</b>            | -20°C (as supplied)   |
| <b>Stability:</b>          | ≥1 year   |
| <b>Storage Buffer:</b>     | PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide   |
| <b>Clone:</b>              | RM368   |
| <b>Host:</b>               | Rabbit  |
| <b>Isotype:</b>            | IgG   |
| <b>Applications:</b>       | Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:200-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

### Images



WB of TT cell lysate using TPO (C-Term) Rabbit Monoclonal Antibody (Clone RM368) at a dilution of 1:200.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human thyroid tissue using TPO (C-Term) Rabbit Monoclonal Antibody (Clone RM368) at a dilution of 1:1,000.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/20/2024

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Thyroid peroxidase (TPO) is a heme-containing enzyme that is essential for the biosynthesis of thyroid hormones.<sup>1,2</sup> It exists as a homodimer where each monomer is composed of an extracellular N-terminal signal peptide and propeptide, an ectodomain containing MPO-, CCP-, and EGF-like domains, a transmembrane domain, and an intracellular C-terminal domain.<sup>2</sup> TPO is regulated primarily at the transcriptional level by thyroid-stimulating hormone (TSH), which upregulates *TPO* expression. TPO is expressed by thyroid follicular cells and localizes to the apical plasma membrane in its active state.<sup>1,2</sup> It catalyzes the oxidation of iodine, resulting in the formation of monoiodotyrosine (MIT) and diiodotyrosine (DIT) residues on thyroglobulin, a protein that functions as a scaffold for thyroid hormone biosynthesis.<sup>3</sup> TPO subsequently catalyzes the oxidative coupling between MIT and DIT residues on thyroglobulin to produce the thyroid hormones triiodothyronine ( $T_3$ ) and thyroxine ( $T_4$ ).<sup>3</sup> Serum TPO autoantibodies are a hallmark of autoimmune thyroid diseases, including Hashimoto's disease and Graves' disease, and have been found in patients with breast cancer.<sup>2,4</sup> Cayman's TPO (C-Term) Rabbit Monoclonal Antibody (Clone RM368) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Ruf, J. and Carayon, P. Structural and functional aspects of thyroid peroxidase. *Arch. Biochem. Biophys.* **445(2)**, 269-277 (2006).
2. Godlewska, M. and Banga, P.J. Thyroid peroxidase as a dual active site enzyme: Focus on biosynthesis, hormonogenesis and thyroid disorders of autoimmunity and cancer. *Biochimie* **160**, 34-45 (2019).
3. Citterio, C.E., Targovnik, H.M., and Arvan, P. The role of thyroglobulin in thyroid hormonogenesis. *Nat. Rev. Endocrinol.* **15(6)**, 323-338 (2019).
4. Fröhlich, E. and Wahl, R. Thyroid autoimmunity: Role of anti-thyroid antibodies in thyroid and extra-thyroidal diseases. *Front. Immunol.* **8**, 521 (2017).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM