

Thrombin- antithrombin monoclonal antibody (22)

Thrombin (Factor IIa) is an enzyme, a serine protease, which participates in blood clotting. It is formed when prothrombin (Factor II) is proteolytically cleaved to form thrombin, which is glycoprotein formed by two peptides chains of 36 and 259 amino acids. Thrombin converts soluble fibrinogen into insoluble strands of fibrin and is then neutralized by antithrombin in order to form a thrombin-antithrombin (TAT) complex.

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

Ordering Information

[Order Online »](#)

| | |
|-------------------|-------|
| BPD-ABS-014-22-04 | 400µg |
|-------------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

| | |
|--------------------|---|
| Alternative Name | TAT |
| Application | ELISA, IHC |
| Clone | 22 |
| Formulation | Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide. |
| GenBank ID | 2147 |
| Host | Mouse |
| Immunogen | Thrombin-antithrombin (TAT) complexes |
| Isotype | IgG1κ |
| Purity Detail | Protein A/G-purified. |
| Source | From culture supernatant. |
| Species Reactivity | Human |
| Specificity | Specific for human Thrombin-antithrombin (TAT) complexes |
| UniProt ID | P00734 (Prothrombin) |

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024