

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-014-22-1	1mg
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](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-
fr@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)