

Factor XII (human, heavy chain) monoclonal antibody (10-11-5)

Factor XII is a component in the coagulation cascade. It is a plasma glycoprotein that participates in the initiation of blood coagulation, fibrinolysis, and the generation of bradykinin and angiotensin. Prekallikrein is cleaved by factor XII to form kallikrein, which then cleaves factor XII to a-factor XIIa, which in turn activates factor XI to factor XIa.

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

Ordering Information

[Order Online »](#)

BPD-CAM-011-05-1	1mg
BPD-CAM-011-05-04	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 Coagulation factor XII, Hageman factor

Application ELISA, WB

Clone 10-11-5

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Affinity purified human factor XII.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Binds the heavy chain of human factor XII, including factor XII derivatives such as α-factor XIIa and α-factor XIIa C1-inhibitor complexes.

UniProt ID P00748

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)