

Factor XII (human, heavy chain) monoclonal antibody (10-11-37)

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-011-37-1	1mg
BPD-CAM-011-37-04	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	10-11-37
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	IgG2ak
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](#)