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 afactor 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 IgG1ĸ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

SpecificityBinds 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

