
F7 Polyclonal Antibody, HRP Conjugated

(Catalog #A62531)

Background

Initiates the extrinsic pathway of blood coagulation. Serine protease that circulates in the blood in a zymogen form. Factor VII is converted to factor VIIa by factor Xa, factor XIIa, factor IXa, or thrombin by minor proteolysis. In the presence of tissue factor and calcium ions, factor VIIa then converts factor X to factor Xa by limited proteolysis. Factor VIIa will also convert factor IX to factor IXa in the presence of tissue factor and calcium.

Description

F7 Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P08709

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Coagulation factor VII protein (61-212AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Coagulation factor VII, Proconvertin, Serum prothrombin conversion accelerator, SPCA, INN: Eptacog alfa, F7

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
