Factor Xa (human plasma), (purified)

Factor Xa (FXa) is a serine endopeptidase comprising a light and a heavy chain, the latter being the catalytic subunit. As part of the prothrombinase complex in the clotting cascade, FXa converts prothrombin to thrombin and is important in cardiovascular research and drug development.

Ordering Information

Order Online »

BML-SE362-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

MW 46 kDa

Source Isolated from human plasma.

P00742 **UniProt ID**

Last modified: May 29, 2024



info-

eu@enzolifesciences.com