

Factor VIIa (human plasma), (purified)

FVIIa is a serine endopeptidase comprising a light and a heavy chain, the latter being the catalytic subunit. When complexed with tissue factor (TF), FVIIa converts factors X and IX to factors Xa and IXa as part of the clotting cascade, making it important in cardiovascular research and drug development.

Ordering Information

[Order Online »](#)

BML-SE438-0010	10µg
-----------------------	------

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

Formulation Liquid. In 20mM TRIS-HCl, pH 7.4, containing 0.1M NaCl.

MW 50 kDa

Source Isolated from human plasma and activated.

UniProt ID P08709

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)