

Thrombin (human plasma), (purified)

Thrombin (factor IIa) is a serine endopeptidase comprising a light and a heavy chain, the latter being the catalytic subunit. Thrombin is a necessary component of the clotting cascade, and is important in cardiovascular research and drug development, and may play a role in Alzheimer's disease as well.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE363-1000	1000U
----------------	-------

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 37 kDa

Source The human Thrombin was activated from homogeneous human prothrombin by activation with Factor Xa, Factor Va, and phospholipid. These activation enzymes were removed after activation. Complete activation was observed on SDS-PAGE. The above protein was purified from Human plasma units that test negative for antibodies to HBc, HCV, HIV 1/2, HTLV-I/II, and non-reactive for HBsAg, HIV-1 RNA, HCV RNA, HBV DNA and Syphilis by FDA approved tests. Donors are screened for CJD.

UniProt ID P00734



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)