



6405 Mira Mesa Blvd. Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Human α Thrombin

Catalog #: 80111
Lot#: 230321

Description: The Human Thrombin was activated from homogeneous human prothrombin by activation with Factor Xa, Factor Va, and phospholipid. These activating 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.

Concentration: 3.41 mg/ml

Aliquots: 1,000 units

MW: 37 kDa

Specific Activity: 3027 NIH units/mg

Total Protein: 0.33 mg

Buffer: 50 mM sodium citrate, 0.2 M NaCl, 0.1% PEG-8000, pH 6.5

Storage: < -60°C

Expiration Date: If properly stored and handled, this enzyme is expected to maintain activity for at least 6 months after shipping.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com