



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Human alpha Thrombin

Catalog #: 80111
Lot#: 150731

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 that was tested and found negative by FDA accepted methods for Anti-HIV1/2, Anti-HTLV I & II, HBsAg, Anti-HCV, Syphilis, HBC Ab, HIV-1 p24 Ag or HIV-1 RNA, HCV RNA and HBV RNA. Donors are screened for CJD (Creutzfeldt-Jakob Disease). MW = 37 kDa.

Concentration: 3.41 mg/ml

Aliquots: 1,000 units

Volume/aliquot: 0.097 ml

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: < -70°C

Expiration Date: 2/2017

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com