



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 Factor Xa

Catalog #: 80112
Lot#: 121114

Description: The Human Factor Xa was activated from Human Factor X by activation with Russell's Viper Venom. This RVV-X was removed after activation. Complete activation is 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 = 46 kDa.

Concentration: 1.59 mg/ml

Aliquots: 100 units

Volume/aliquot: 0.503 ml

Specific Activity: 216.60 IU/mg

Total Protein: 0.80 mg

Buffer: 20 mM Bis Tris, 0.7 M NaCl, pH 6.0

Storage: < -70°C

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