

**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:** 2.16 mg/ml

**Aliquots:** 1 x 0.80 mg

**Volume/aliquot:** 0.365 ml

**Specific Activity:** 244 IU/mg

**Total Protein:** 0.80 mg

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

**Storage:** -60°C

**Expiration Date:** 05/2024