

Factor Xa Protease Assay Service

Factor Xa [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

Target Overview

CAT#	Factor Xa
Enzyme	Human Factor Xa
Synonym	Coagulation factor Xa, Stuart factor, F10
Substrates	Fluorogenic peptide CH3SO2-D-CHA-Gly-Arg-AMC-AcOH, 10 μ M

Assay Properties

Readout	Enzymatic activity				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (μM)</th> </tr> </thead> <tbody> <tr> <td>Gabexate mesylate</td> <td>1.92</td> </tr> </tbody> </table>	Compounds	IC50 (μ M)	Gabexate mesylate	1.92
Compounds	IC50 (μ M)				
Gabexate mesylate	1.92				
Scientific Information	P00742 (https://www.uniprot.org/uniprotkb/P00742/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Factor Xa

