

Factor XIa Protease Assay Service

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

Target Overview

CAT#	Factor XIa
Enzyme	Human Factor XIa
Synonym	Coagulation factor XIa, Plasma thromboplastin antecedent, PTA, F11
Substrates	Fluorogenic peptide (Boc-Glu(OBzl)-Ala-Arg)-MCA, 10 μ M

Assay Properties

Readout	Enzymatic activity				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Gabexate mesylate</td> <td>231</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Gabexate mesylate	231
Compounds	IC50 (nM)				
Gabexate mesylate	231				
Scientific Information	P03951 (https://www.uniprot.org/uniprotkb/P03951/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Factor XIa

