

Plasma Kallikrein Protease Assay Service

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

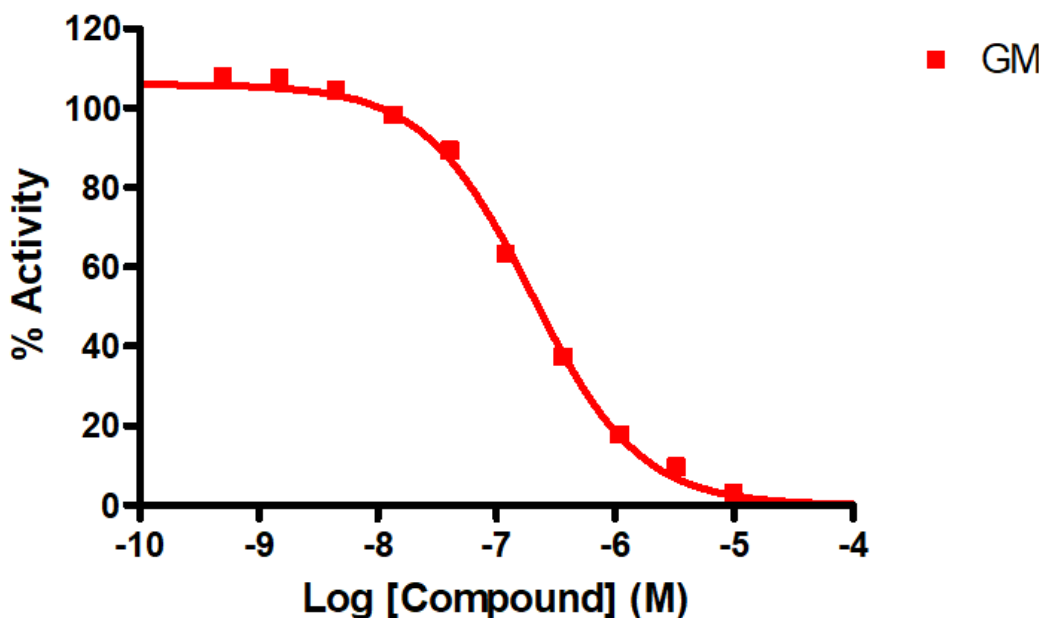
Target Overview

CAT#	Plasma Kallikrein
Enzyme	Human Plasma Kallikrein
Synonym	Fletcher factor, Kininogenin, PKK, KLK3, KLKB1
Substrates	10 μ M Fluorogenic peptide Z-FR-AMC

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>138</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Gabexate mesylate	138
Compounds	IC50 (nM)				
Gabexate mesylate	138				
Scientific Information	P03952 (https://www.uniprot.org/uniprotkb/P03952/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Plasma Kallikrein



Let's Discover Together.

© Jun 2023 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/plasma_kallikrein_protea_malvern

