

Kallikrein 5 Protease Assay Service

Kallikrein 5 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

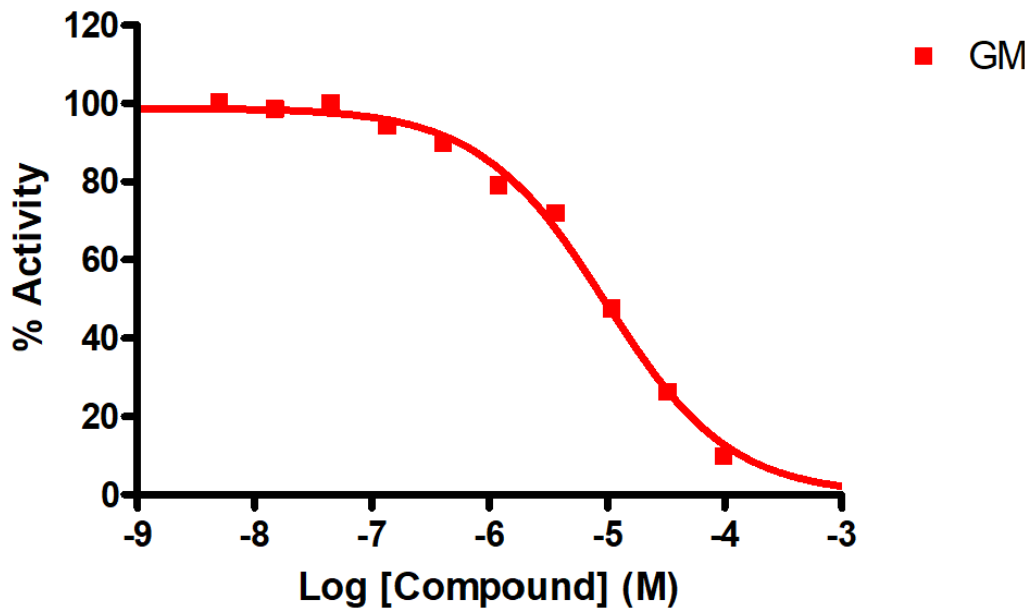
Target Overview

CAT#	Kallikrein 5
Enzyme	Human Kallikrein 5
Synonym	KLK5, KLK-L2, SCTE
Substrates	Fluorogenic peptide Z-VVR-AMC, 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>9.07</td> </tr> </tbody> </table>	Compounds	IC50 (μ M)	Gabexate mesylate	9.07
Compounds	IC50 (μ M)				
Gabexate mesylate	9.07				
Scientific Information	Q9Y337 (https://www.uniprot.org/uniprotkb/Q9Y337/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Kallikrein 5



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/kallikrein_5_protea_malvern

