

# Kallikrein 13 Protease Assay Service

Kallikrein 13 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

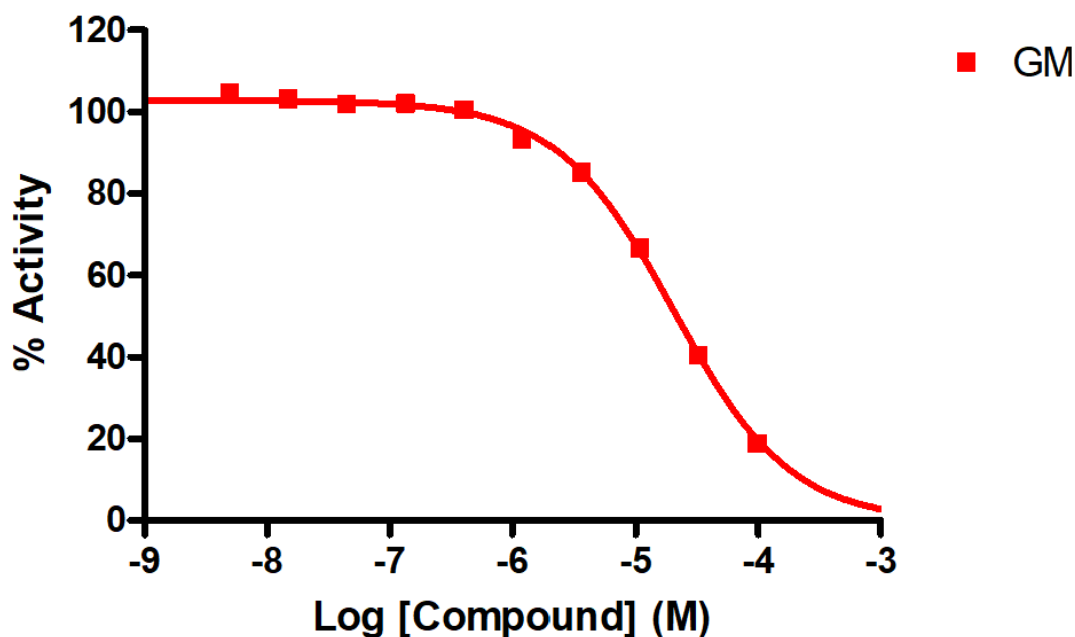
## Target Overview

CAT#	Kallikrein 13
Enzyme	Human Kallikrein 13
Synonym	Kallikrein-like protein 4, KLK13
Substrates	Fluorogenic peptide VPR-AMC, 10 $\mu$ M

## Assay Properties

Readout	Enzymatic activity				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (<math>\mu</math>M)</th> </tr> </thead> <tbody> <tr> <td>Gabexate mesylate</td> <td>12.6</td> </tr> </tbody> </table>	Compounds	IC50 ( $\mu$ M)	Gabexate mesylate	12.6
Compounds	IC50 ( $\mu$ M)				
Gabexate mesylate	12.6				
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/Q9UKR3/entry">Q9UKR3</a> (https://www.uniprot.org/uniprotkb/Q9UKR3/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website <a href="#">Protease Assay Services</a>				

## Reference compound IC50 for Kallikrein 13



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/kallikrein\\_13\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/kallikrein_13_protea_malvern)

