

# Kallikrein 7 Protease Assay Service

Kallikrein 7 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

## Target Overview

CAT#	Kallikrein 7
Enzyme	Human Kallikrein 7
Synonym	Serine protease 6, PRSS6, SCCE, KLK7
Substrates	Fluorogenic peptide MCA-RPLPVE-Nval-WRK(Dup)-NH <sub>2</sub> , 10 μM

## 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>29.9</td> </tr> </tbody> </table>	Compounds	IC50 (μM)	Gabexate mesylate	29.9
Compounds	IC50 (μM)				
Gabexate mesylate	29.9				
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P49862/entry">P49862</a> ( <a href="https://www.uniprot.org/uniprotkb/P49862/entry">https://www.uniprot.org/uniprotkb/P49862/entry</a> )				
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 7

