

Kallikrein 2 Protease Assay Service

Kallikrein 2 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

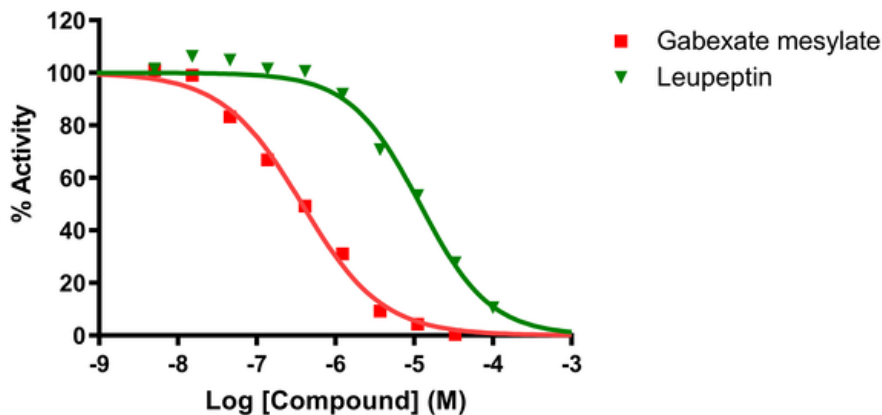
Target Overview

CAT#	Kallikrein 2
Enzyme	Kallikrein 2
Synonym	KLK2, Glandular kallikrein-1, hGK-1, Tissue kallikrein-2
Substrates	Fluorogenic substrate PFR-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (μM)
	Gabexate Mesylate	0.38
	Leupeptin	11.8
Scientific Information	P20151 (http://www.uniprot.org/uniprot/P20151)	
Screening Location	Malvern, USA	

Reference Compound IC50 for Kallikrein 2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/kallikrein_2_protea_malvern

