

Kallikrein 14 Protease Assay Service

Kallikrein 14 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

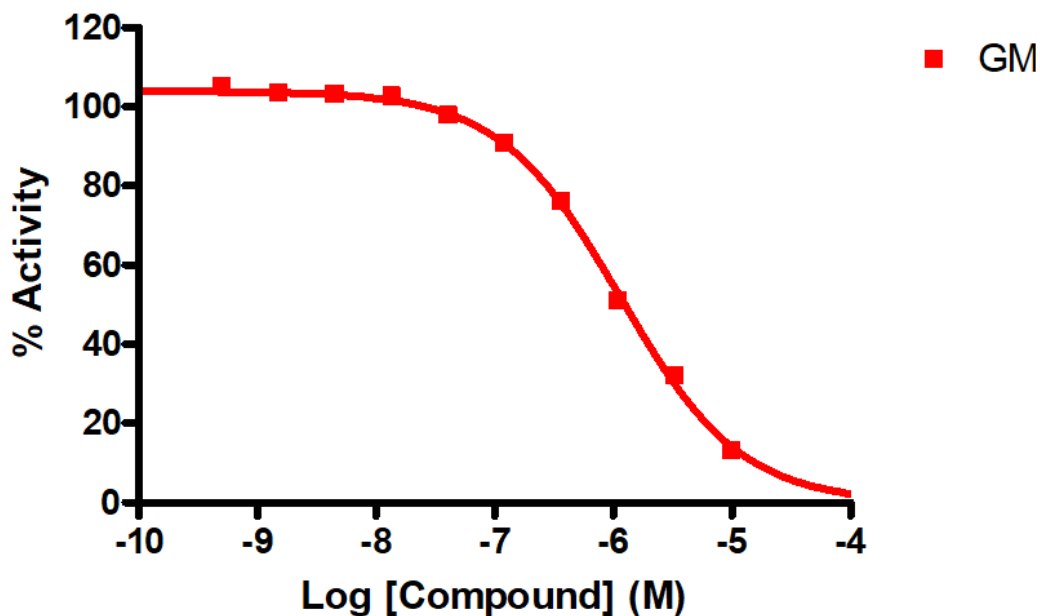
Target Overview

CAT#	Kallikrein 14
Enzyme	Human Kallikrein 14
Synonym	Kallikrein-like protein 6, KLK14
Substrates	Fluorogenic peptide VPR-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	Gabexate mesylate	678
Scientific Information	Q9P0G3 (https://www.uniprot.org/uniprotkb/Q9P0G3/entry)	
Screening Location	Malvern, PA, USA	
Further Information	More information can be found on our website Protease Assay Services	

Reference compound IC50 for Kallikrein 14



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/kallikrein_14_protea_malvern

