

Hepsin Protease Assay Service

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

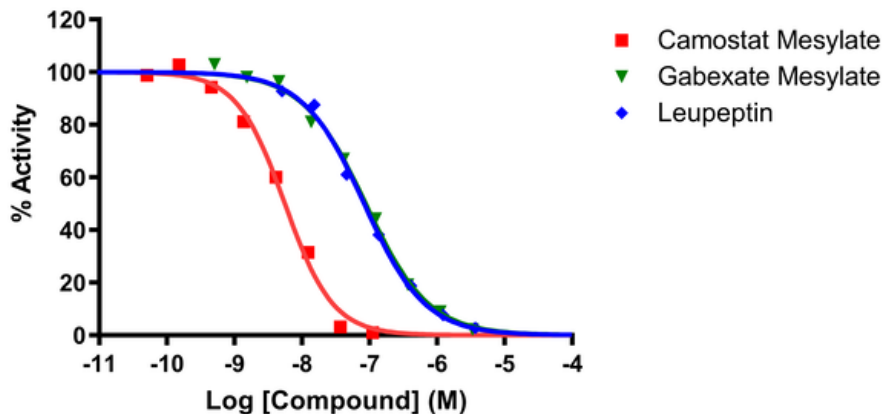
Target Overview

CAT#	Hepsin
Enzyme	Human Serine Protease Hepsin
Synonym	Transmembrane protease serine 1, TMPRSS1, HPN
Substrates	Fluorogenic substrate Boc-QRR-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (nM)
	Camostat Mesylate	5.57
	Gabexate Mesylate	87.25
	Leupeptin	82.99
Scientific Information	P05981 (http://www.uniprot.org/uniprot/P05981)	
Screening Location	Malvern, USA	

Reference Compound IC50 for Hepsin



Let's Discover Together.

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

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

