

tPA Protease Assay Service

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

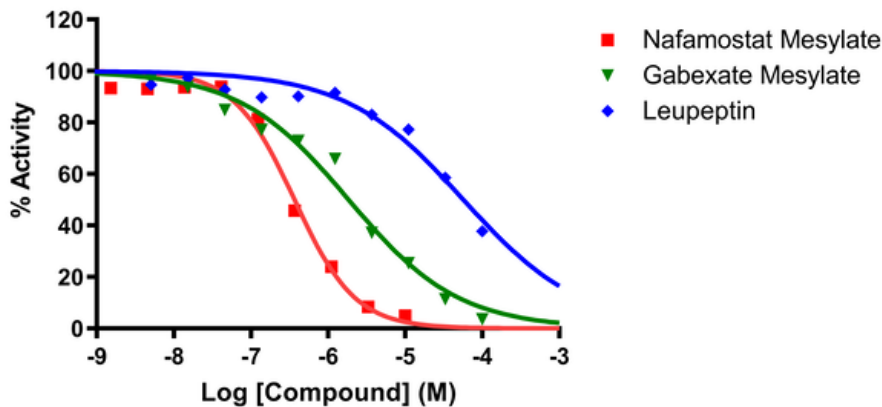
Target Overview

CAT#	tPA
Enzyme	Tissue Plasminogen Activator
Synonym	Tissue-type plasminogen activator, tPA, Alteplase, Reteplase
Substrates	Fluorogenic substrate Z-GPR-AMC, 50 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (μM)
	Nafamostat Mesylate	0.37
	Gabexate Mesylate	1.85
	Leupeptin	57.99
Scientific Information	P00750 (http://www.uniprot.org/uniprot/P00750)	
Screening Location	Malvern, USA	

Reference Compound IC50 for tPA



Let's Discover Together.

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

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

