

Activated Protein C Protease Assay Service

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

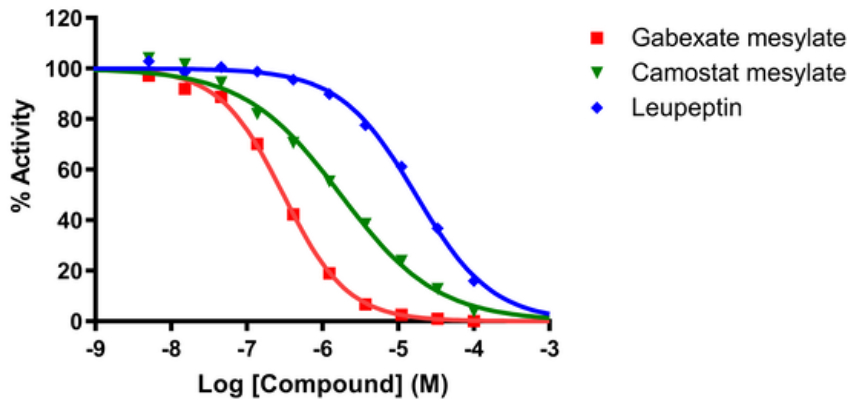
Target Overview

CAT#	Activated Protein C
Enzyme	Human Activated Protein C
Synonym	Vitamin K-dependent Protein C, APC, PROC
Substrates	Fluorogenic substrate Z-VLR-AMC, 50 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (μM)
	Gabexate mesylate	0.3
	Camostat mesylate	1.7
	Leupeptin	17.2
Scientific Information	P04070 (http://www.uniprot.org/uniprot/P04070)	
Screening Location	Malvern, USA	

Reference Compound IC50 for Activated Protein C



Let's Discover Together.

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

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

