

Tryptase g1 Protease Assay Service

Tryptase g1 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

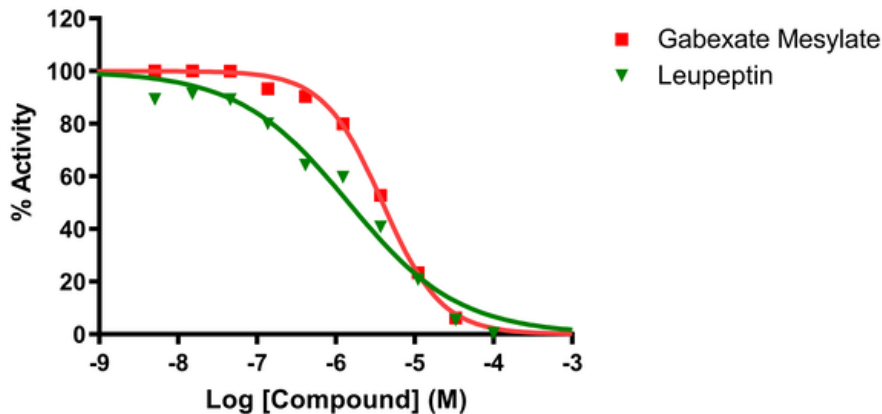
Target Overview

CAT#	Tryptase g1
Enzyme	Human Tryptase gamma 1
Synonym	TPSG1, Serine protease 31, Transmembrane tryptase
Substrates	Fluorogenic substrate Mca-RPKPVE-Nval-WRK(Dnp)-NH ₂ , 5 μ M

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (μM)
	Gabexate Mesylate	3.9
	Leupeptin	1.5
Scientific Information	Q9NRR2 (http://www.uniprot.org/uniprot/Q9NRR2)	
Screening Location	Malvern, PA, USA	
Further information	More information can be found on our website Protease Assay Services .	

Reference Compound IC50 for Tryptase g1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/tryptase_g1_protea_malvern

