

Cathepsin G Protease Assay Service

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

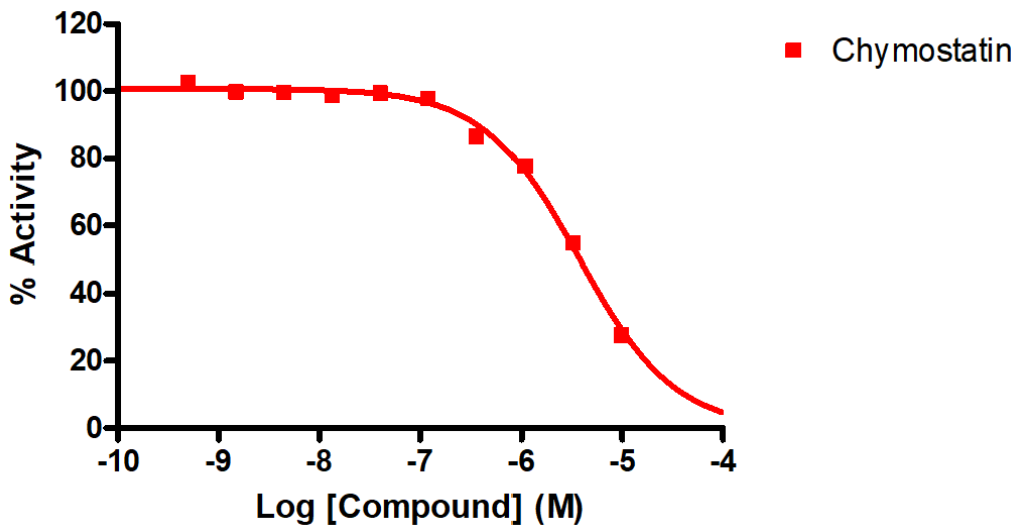
Target Overview

CAT#	Cathepsin G
Enzyme	Cathepsin G
Synonym	CTSG, CG
Substrates	Fluorogenic substrate Suc-AAPF-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (μM) Chymostatin 2.95
Scientific Information	P08311 (https://www.uniprot.org/uniprotkb/P08311/entry)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Protease Assay Services

Reference compound IC50 for Cathepsin G



Let's Discover Together.

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

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

