

Chymase Protease Assay Service

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

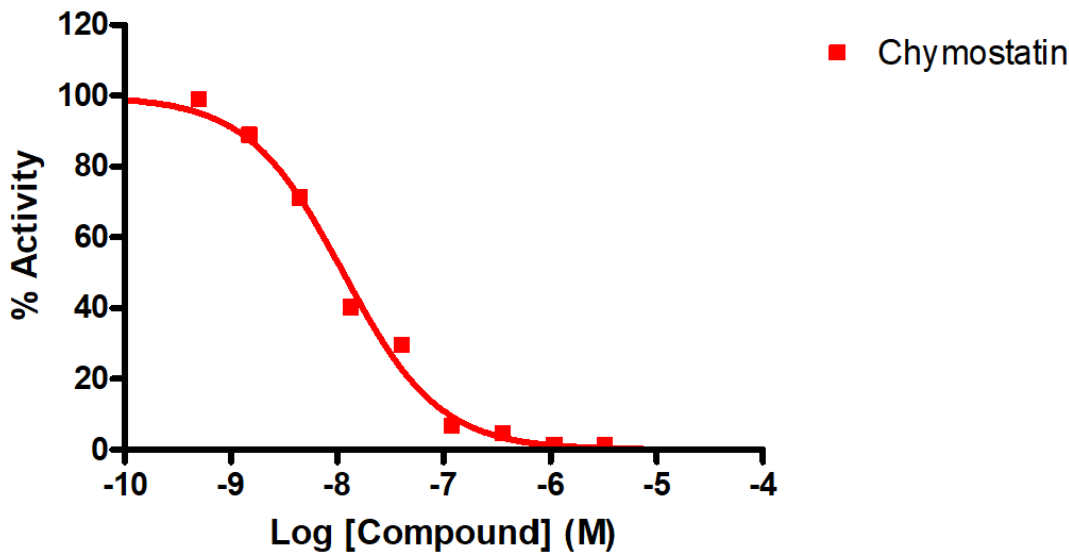
Target Overview

| | |
|------------------|---|
| CAT# | Chymase |
| Enzyme | Chymase |
| Synonym | Mast cell protease I, Skeletal muscle protease, skin chymotryptic proteinase, alpha chymase, CMA1, CYH, CYM |
| Molecular Weight | 30 kDa |
| Substrates | Fluorogenic substrate Suc-AAPF-AMC, 10 μ M |

Assay Properties

| | |
|--------------------------|---|
| Readout | Enzymatic activity |
| Reference Compound IC50s | Compounds IC50 (nM) Chymostatin 13.7 |
| Scientific Information | P23946 (https://www.uniprot.org/uniprotkb/P23946/entry) |
| Screening Location | Malvern, PA, USA |
| Further information | More information can be found on our website Protease Assay Services . |

Reference compound IC50 for Chymase



Let's Discover Together.

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

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

