

Chymotrypsin Protease Assay Service

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

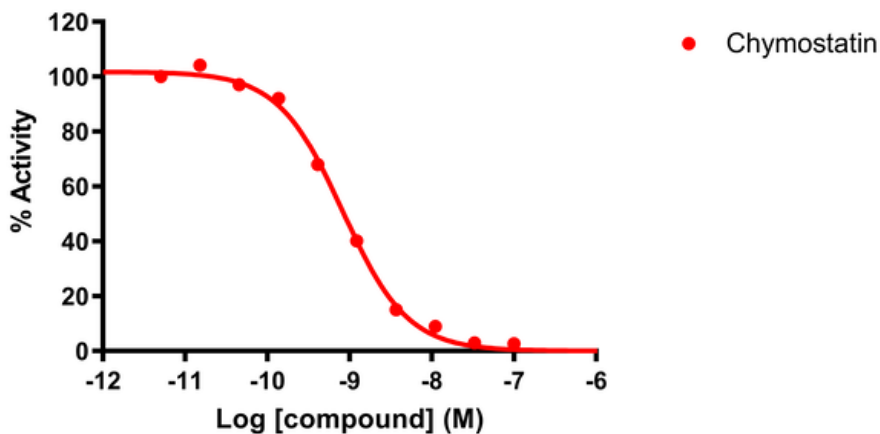
Target Overview

CAT#	Chymotrypsin
Enzyme	Human Chymotrypsin
Synonym	Alpha-chymotrypsin
Substrates	Fluorogenic peptide [Suc-AAPF-AMC], 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) Chymostatin 0.8
Scientific Information	Q99895 (https://www.uniprot.org/uniprot/Q99895)
Screening Location	Malvern, USA
Further information	More information can be found on our website Protease Assay Services .

Reference Compound IC50 for Chymotrypsin



Let's Discover Together.

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

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

