

Cathepsin B Protease Assay Service

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

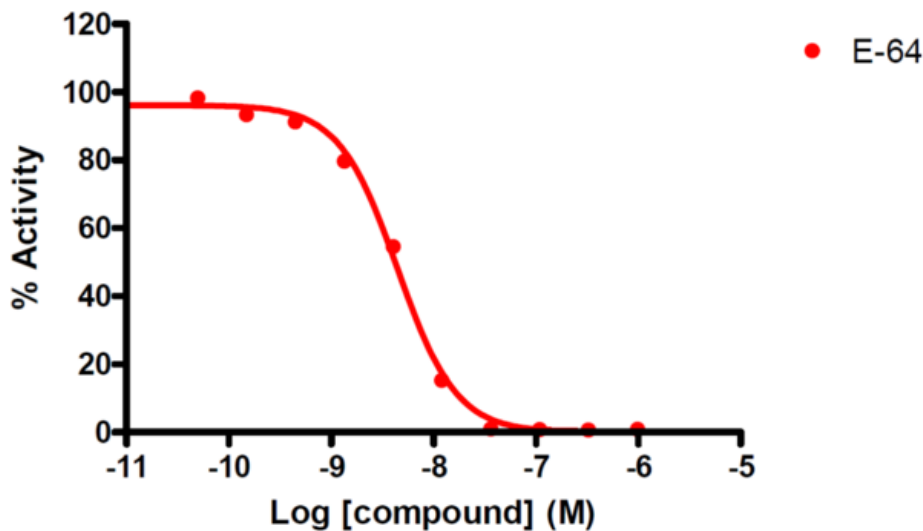
Target Overview

CAT#	Cathepsin B
Enzyme	Human Cathepsin B
Synonym	APP secretase, Cathepsin B1, CATB, CTSB, CPSB
Substrates	Fluorogenic peptide [Z-FR-AMC], 10 μ M

Assay Properties

Reference Compound IC50s	Compounds IC50 (nM) E-64 4.7
Scientific Information	P07858 (http://www.uniprot.org/uniprot/P07858)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for Cathepsin B



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

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

