

Cathepsin L Protease Assay Service

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

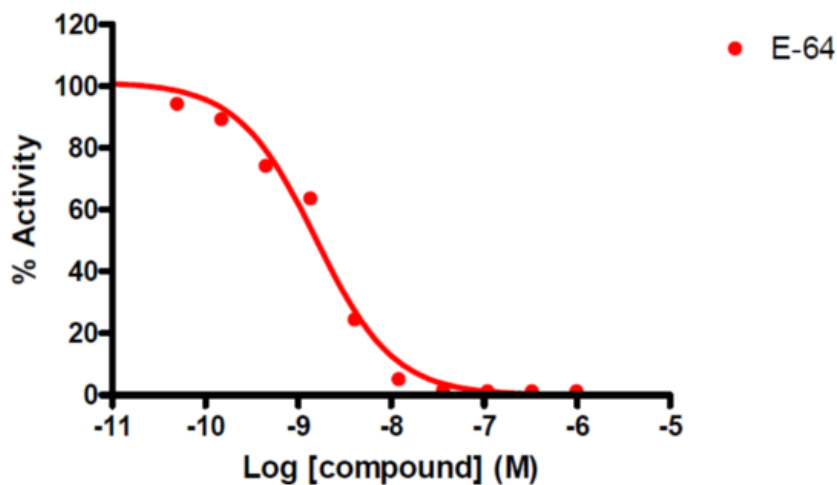
Target Overview

CAT#	Cathepsin L
Enzyme	Human Cathepsin L
Synonym	CTSL, CTSL1, MEP
Substrates	Fluorogenic peptide [Z-FR-AMC], 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) E-64 1.6
Scientific Information	P07711 (http://www.uniprot.org/uniprot/P07711)
Screening Location	Malvern, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for Cathepsin L



Let's Discover Together.

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

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

