

Cathepsin V Protease Assay Service

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

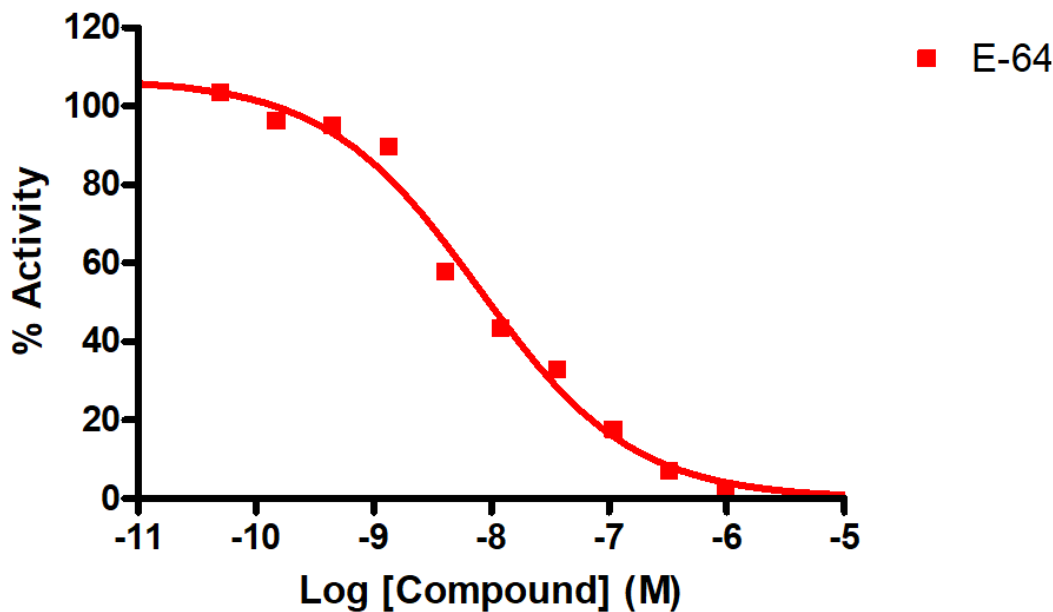
Target Overview

CAT#	Cathepsin V
Enzyme	Cathepsin V
Synonym	Cathepsin L2, CTSL2, CATL2, CTSU, CTSV
Substrates	Fluorogenic substrate Z-FR-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) E-64 6.45
Scientific Information	O60911 (https://www.uniprot.org/uniprotkb/O60911/entry)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for Cathepsin V



Let's Discover Together.

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

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

