

# Cathepsin C Protease Assay Service

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

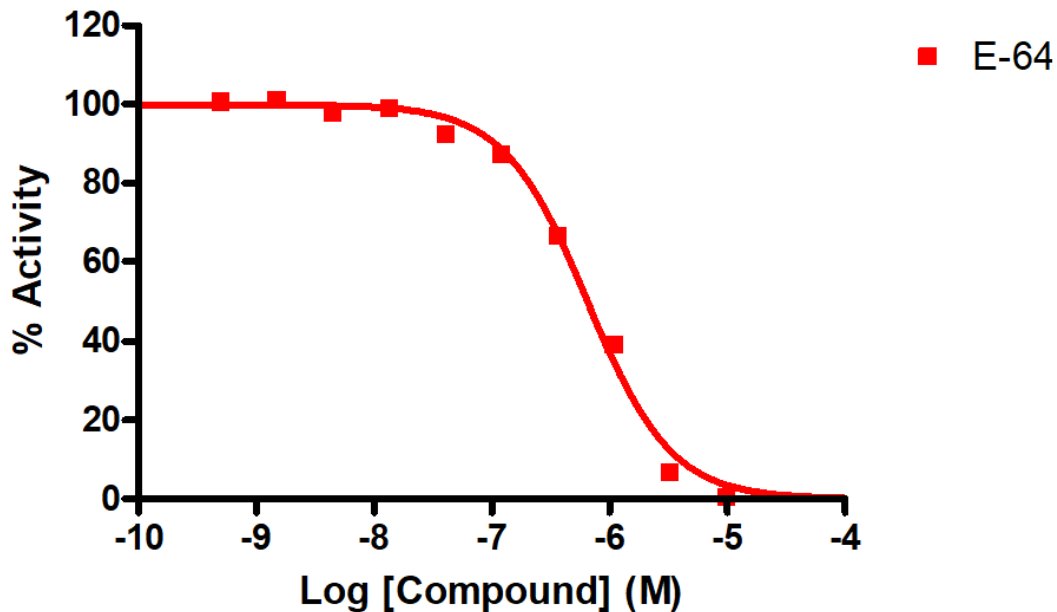
## Target Overview

CAT#	Cathepsin C
Enzyme	Cathepsin C
Synonym	Dipeptidyl-peptidase (DPPI), Cathepsin J, Dipeptidyl transferase, CTSC, CPPI, CATC
Substrates	Fluorogenic substrate H-GR-AMC, 10 $\mu$ M

## Assay Properties

Reference Compound IC50s	<b>Compounds IC50 (nM)</b> E-64      771
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P53634/entry">P53634</a> (https://www.uniprot.org/uniprotkb/P53634/entry)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website <a href="#">Protease Assay Services</a> .

## Reference compound IC50 for Cathepsin C



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/cathepsin\\_c\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/cathepsin_c_protea_malvern)

