

# Cathepsin S Protease Assay Service

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

## Target Overview

CAT#	Cathepsin S
Enzyme	Cathepsin S
Synonym	CTSS, CATS
Substrates	Fluorogenic substrate Z-FR-AMC, 10 $\mu$ M

## Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> E-64      1.67
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P25774/entry">P25774</a> ( <a href="https://www.uniprot.org/uniprotkb/P25774/entry">https://www.uniprot.org/uniprotkb/P25774/entry</a> )
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 S

