

Caspase 2 Protease Assay Service

Caspase 2 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

Target Overview

CAT#	Caspase 2
Enzyme	Human Caspase 2
Synonym	Neural precursor cell expressed developmentally down-regulated protein 2, Protease ICH-1, CASP2, ICH1, NEDD2
Substrates	Fluorogenic peptide [Ac-LEHD-AMC], 5 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (μM) IETD-CHO 1.71
Scientific Information	P42575 (http://www.uniprot.org/uniprot/P42575)
Screening Location	Malvern, USA

Reference compound IC50 for Caspase 2

