

Papain Protease Assay Service

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

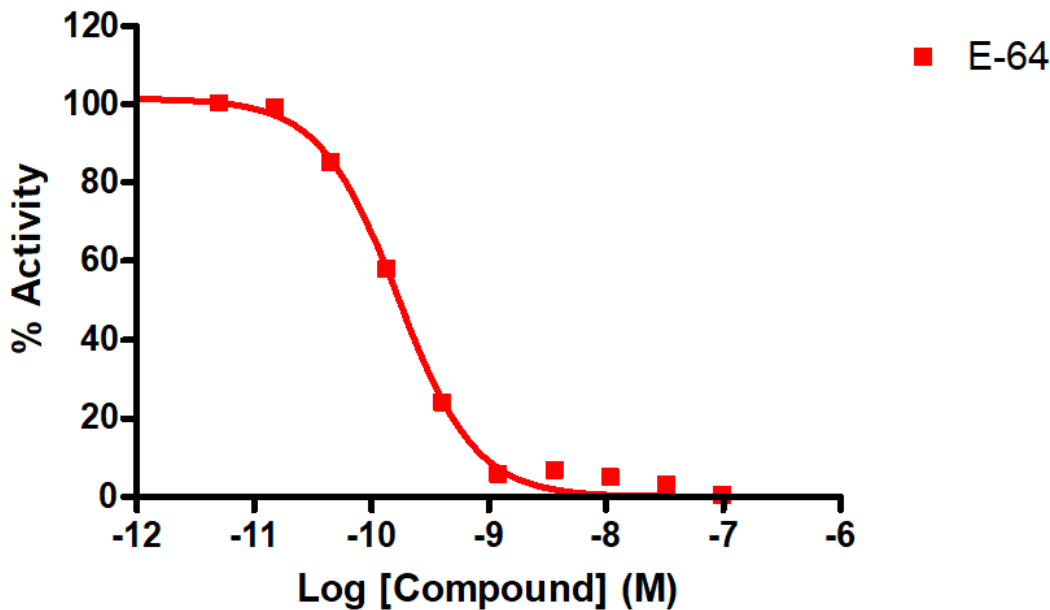
Target Overview

CAT#	Papain
Enzyme	Papaya Papain
Synonym	Papaya proteinase I
Substrates	10 μ M Fluorogenic peptide Z-FR-AMC

Assay Properties

Readout	Enzymatic activity				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Gabexate mesylate</td> <td>0.14</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Gabexate mesylate	0.14
Compounds	IC50 (nM)				
Gabexate mesylate	0.14				
Scientific Information	P00784 (https://www.uniprot.org/uniprotkb/P00784/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Papain (Cysteine Protease)



Let's Discover Together.

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

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

