

Neprilysin Protease Assay Service

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

Target Overview

CAT#	Neprilysin
Enzyme	Human Neprilysin
Synonym	Atriopeptidase, Neutral endopeptidase, Skin fibroblast elastase, CALLA, MME, EPN
Substrates	10 μ M Fluorogenic peptide MCA-RPPGFSAFK(Dnp)-OH

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>Phosphoramidon</td> <td>453</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Phosphoramidon	453
Compounds	IC50 (nM)				
Phosphoramidon	453				
Scientific Information	P08473 (https://www.uniprot.org/uniprotkb/P08473/entry)				
Screening Location	Malvern, PA, USA				
Further Information	More information can be found on our website Protease Assay Services				

Reference compound IC50 for Neprilysin

