

Furin Protease Assay Service

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

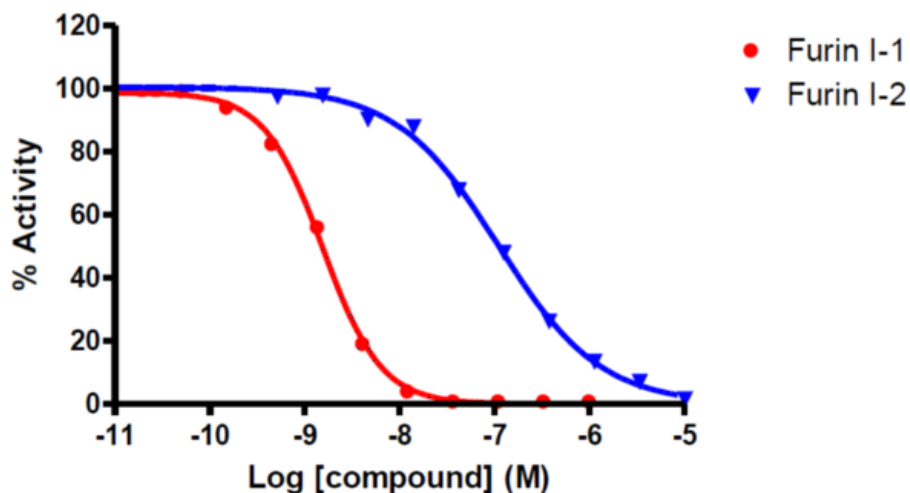
Target Overview

CAT#	Furin
Enzyme	Human Furin
Synonym	FUR, PACE, PCSK3
Substrates	Fluorogenic peptide [pERTKR -AMC], 5 μ M

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>Furin Inhibitor 1</td> <td>1.6</td> </tr> <tr> <td>Furin Inhibitor 2</td> <td>110</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Furin Inhibitor 1	1.6	Furin Inhibitor 2	110
Compounds	IC50 (nM)						
Furin Inhibitor 1	1.6						
Furin Inhibitor 2	110						
Scientific Information	P09958 (http://www.uniprot.org/uniprot/P09958)						
Screening Location	Malvern, USA						
Further information	More information can be found on our website Protease Assay Services .						

Reference compound IC50 for Furin



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

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

