

MRCKa/CDC42BPA Kinase Assay Service

Activity-based MRCKa/CDC42BPA kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

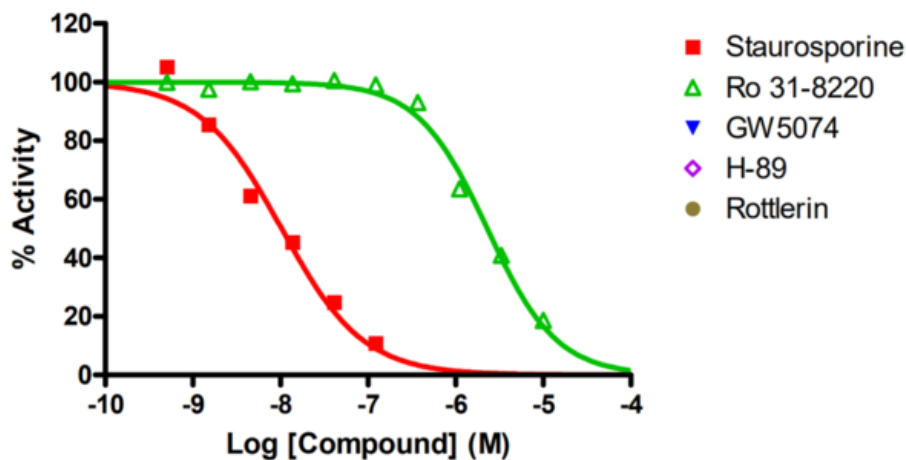
Target Overview

CAT#	MRCKa/CDC42BPA
Enzyme	Human MRCKa/CDC42BPA
HGNC	CDC42BPA
Synonym	MRCK-alpha
Substrates	Peptide substrate, [KEAKEKRQEIQAKRRRLSSLRASTSKSGGSQK], 20 μM ATP 10 μM

Assay Properties

Reaction	Substrate + [³² P]-ATP = ³² P-Substrate + ADP						
Readout	Enzymatic activity						
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>10.3</td> </tr> <tr> <td>Ro 31-8220</td> <td>2,300</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	10.3	Ro 31-8220	2,300
Compounds	IC50 (nM)						
Staurosporine	10.3						
Ro 31-8220	2,300						
Scientific Information	Q5VT25 (http://www.uniprot.org/uniprot/Q5VT25)						
Screening Location	Malvern, USA						

Reference compound IC50 for MRCKa/CDC42BPA



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/mrcka_cdc42bpa_kin_malvern

