

MARK2/PAR-1Ba Kinase Assay Service

Activity-based MARK2/PAR-1Ba kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

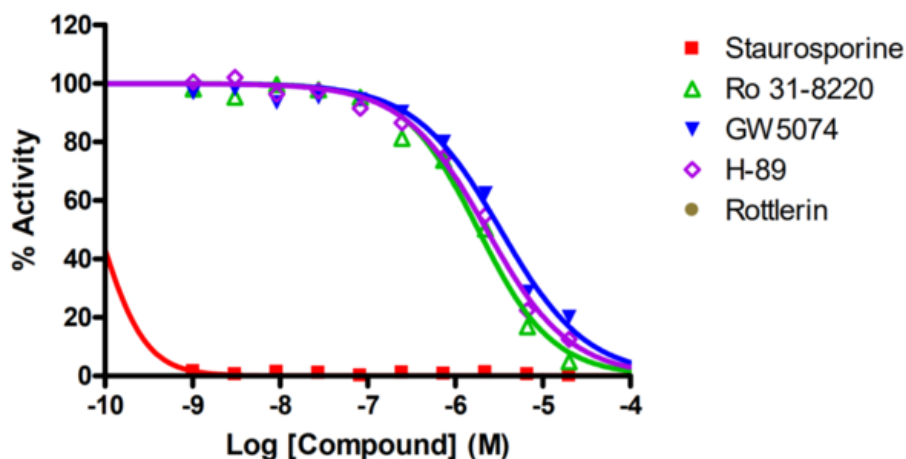
Target Overview

CAT#	MARK2/PAR-1Ba
Enzyme	Human MARK2/PAR-1Ba
HGNC	MARK2
Synonym	EMK1, MGC99619, PAR-1Ba, ELKL
Substrates	Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 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><0.5</td> </tr> <tr> <td>Ro 31-8220</td> <td>1,800</td> </tr> <tr> <td>GW 5074</td> <td>3,300</td> </tr> <tr> <td>H-89</td> <td>2,300</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	<0.5	Ro 31-8220	1,800	GW 5074	3,300	H-89	2,300
Compounds	IC50 (nM)										
Staurosporine	<0.5										
Ro 31-8220	1,800										
GW 5074	3,300										
H-89	2,300										
Scientific Information	Q7KZ17-4 (http://www.uniprot.org/uniprot/Q7KZ17-4)										
Screening Location	Malvern, USA										

Reference compound IC50 for MARK2/PAR-1Ba



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/mark2_par-1ba_kin_malvern

