

# P38d/MAPK13 Kinase Assay Service

Activity-based P38d/MAPK13 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

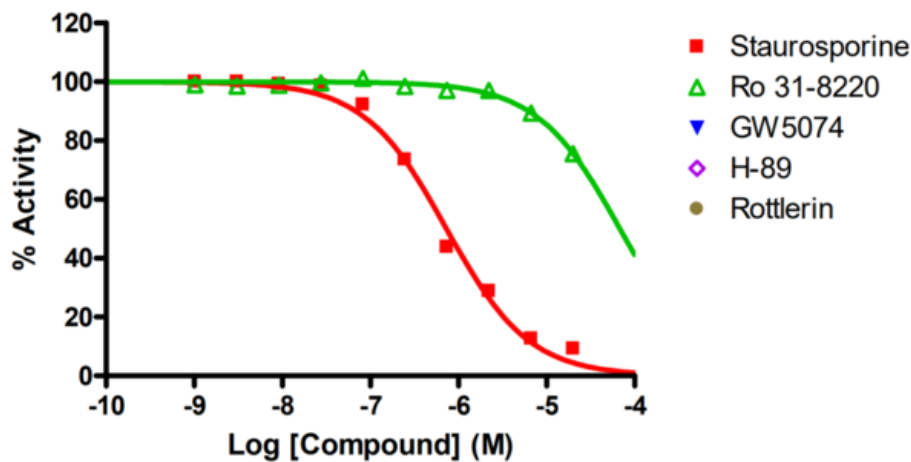
## Target Overview

CAT#	P38d/MAPK13
Enzyme	Human P38d/MAPK13
HGNC	MAPK13
Synonym	p38-delta, PRKM13, SAPK4
Substrates	MBP (Myelin Basic Protein), 20 μM ATP 10 μM

## Assay Properties

Reaction	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> 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>730</td> </tr> <tr> <td>Ro 31-8220</td> <td>&gt;20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	730	Ro 31-8220	>20,000
Compounds	IC50 (nM)						
Staurosporine	730						
Ro 31-8220	>20,000						
Scientific Information	<a href="http://www.uniprot.org/uniprot/O15264">O15264</a> ( <a href="http://www.uniprot.org/uniprot/O15264">http://www.uniprot.org/uniprot/O15264</a> )						
Screening Location	Malvern, USA						

## Reference compound IC50 for P38d/MAPK13



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/p38d\\_mapk13\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/p38d_mapk13_kin_malvern)

