

# MINK/MINK1 Kinase Assay Service

Activity-based MINK/MINK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

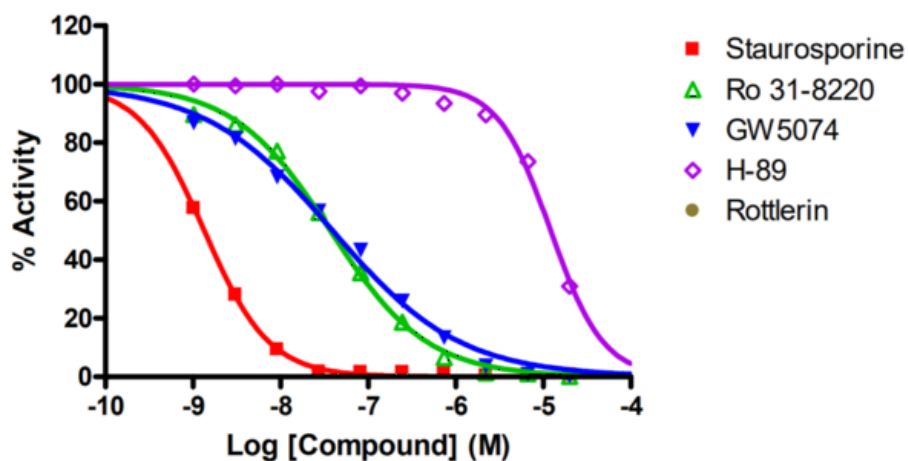
## Target Overview

<b>CAT#</b>	MINK/MINK1
<b>Enzyme</b>	Human MINK/MINK1
<b>HGNC</b>	MINK1
<b>Synonym</b>	B55, YSK2, MAP4K6, ZC3
<b>Substrates</b>	MBP (Myelin Basic Protein), 20 μM ATP 10 μM

## Assay Properties

<b>Reaction</b>	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> P-Substrate + ADP										
<b>Readout</b>	Enzymatic activity										
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>1.3</td> </tr> <tr> <td>Ro 31-8220</td> <td>37</td> </tr> <tr> <td>GW 5074</td> <td>39</td> </tr> <tr> <td>H-89</td> <td>12,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	1.3	Ro 31-8220	37	GW 5074	39	H-89	12,000
Compounds	IC50 (nM)										
Staurosporine	1.3										
Ro 31-8220	37										
GW 5074	39										
H-89	12,000										
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/Q8N4C8">Q8N4C8</a> ( <a href="http://www.uniprot.org/uniprot/Q8N4C8">http://www.uniprot.org/uniprot/Q8N4C8</a> )										
<b>Screening Location</b>	Malvern, USA										

## Reference compound IC50 for MINK/MINK1



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mink\\_mink1\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/mink_mink1_kin_malvern)

