

# MST4 Kinase Assay Service

Activity-based MST4 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

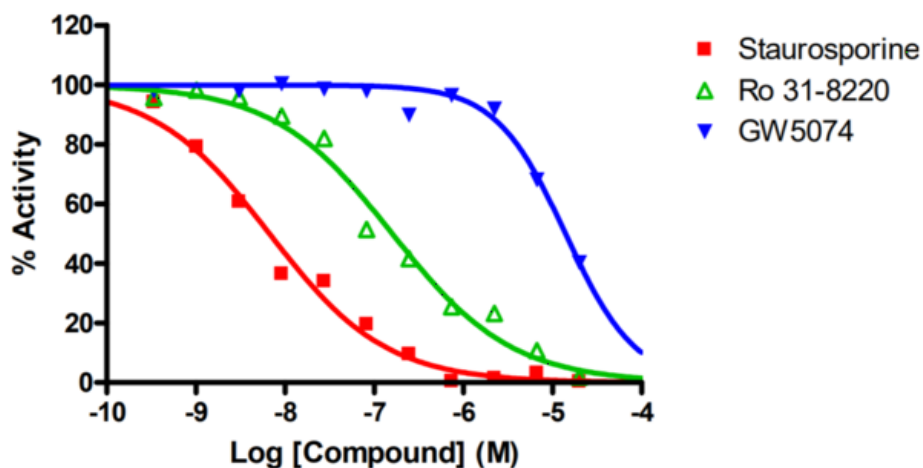
## Target Overview

CAT#	MST4
Enzyme	Human MST4
HGNC	STK26
Synonym	MASK
Substrates	MBP (Myelin Basic Protein), 10 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>6.7</td> </tr> <tr> <td>Ro 31-8220</td> <td>160</td> </tr> <tr> <td>GW 5074</td> <td>14,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	6.7	Ro 31-8220	160	GW 5074	14,000
Compounds	IC50 (nM)								
Staurosporine	6.7								
Ro 31-8220	160								
GW 5074	14,000								
Scientific Information	<a href="http://www.uniprot.org/uniprot/Q9P289">Q9P289</a> ( <a href="http://www.uniprot.org/uniprot/Q9P289">http://www.uniprot.org/uniprot/Q9P289</a> )								
Screening Location	Malvern, USA								

## Reference compound IC50 for MST4



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mst4\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/mst4_kin_malvern)

