

# MASTL Kinase Assay Service

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

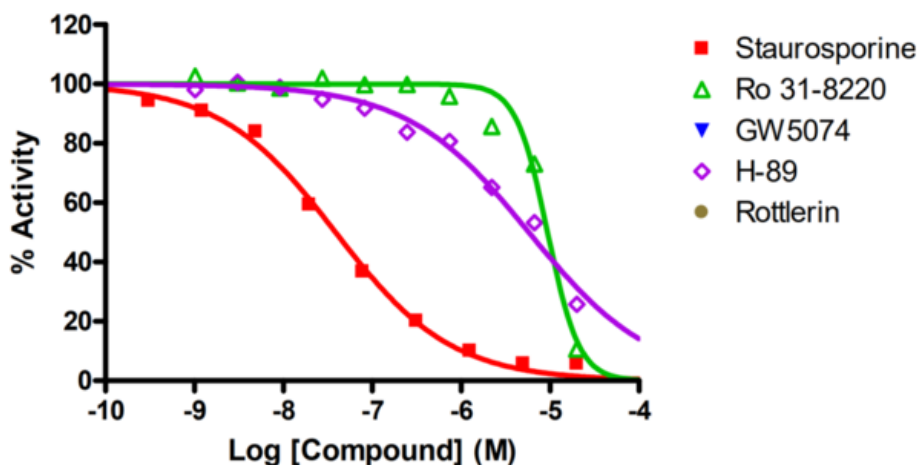
## Target Overview

CAT#	MASTL
Enzyme	Human MASTL
HGNC	MASTL
Synonym	GLW, GREATWALL, GWL, hGWL, THC2
Substrates	ARPP19, 5 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>39</td> </tr> <tr> <td>Ro 31-8220</td> <td>9,400</td> </tr> <tr> <td>GW 5074</td> <td>&gt;100,000</td> </tr> <tr> <td>H-89</td> <td>5,800</td> </tr> <tr> <td>Rottlerin</td> <td>&gt;100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	39	Ro 31-8220	9,400	GW 5074	>100,000	H-89	5,800	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	39												
Ro 31-8220	9,400												
GW 5074	>100,000												
H-89	5,800												
Rottlerin	>100,000												
Scientific Information	<a href="http://www.uniprot.org/uniprot/Q96GX5">Q96GX5</a> ( <a href="http://www.uniprot.org/uniprot/Q96GX5">http://www.uniprot.org/uniprot/Q96GX5</a> )												
Screening Location	Malvern, USA												

## Reference compound IC50 for MASTL



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mastl\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/mastl_kin_malvern)

