

BMPR2 Kinase Assay Service

Activity-based BMPR2 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

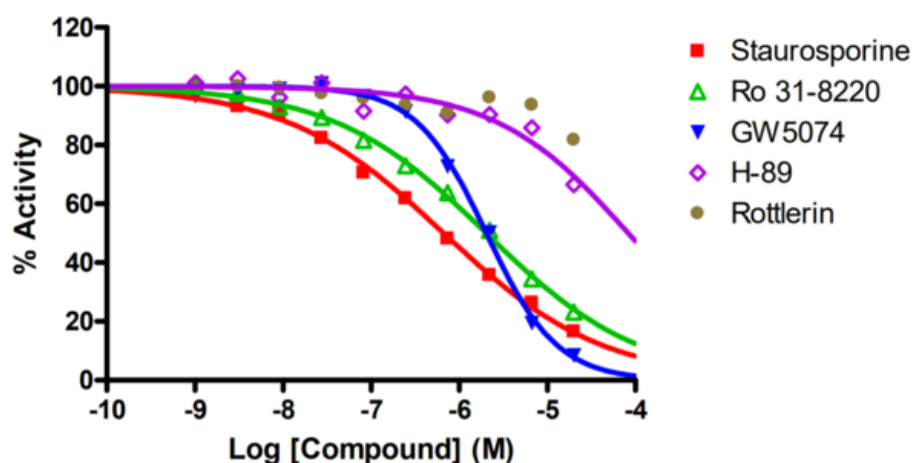
Target Overview

CAT#	BMPR2
Enzyme	Human BMPR2
HGNC	BMPR2
Synonym	BMPR-II, BMPR3, BMR2, BRK-3, PPH1, T-ALK
Substrates	MBP (Myelin Basic Protein), 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>670</td> </tr> <tr> <td>Ro 31-8220</td> <td>2,000</td> </tr> <tr> <td>GW5074</td> <td>2,000</td> </tr> <tr> <td>H-89</td> <td>>50,000</td> </tr> <tr> <td>Rottlerin</td> <td>>50,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	670	Ro 31-8220	2,000	GW5074	2,000	H-89	>50,000	Rottlerin	>50,000
Compounds	IC50 (nM)												
Staurosporine	670												
Ro 31-8220	2,000												
GW5074	2,000												
H-89	>50,000												
Rottlerin	>50,000												
Scientific Information	Q13873 (http://www.uniprot.org/uniprot/Q13873)												
Screening Location	Malvern, USA												

Reference compound IC50 for BMPR2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/bmpr2_kin_malvern

