

STK33 Kinase Assay Service

Activity-based STK33 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

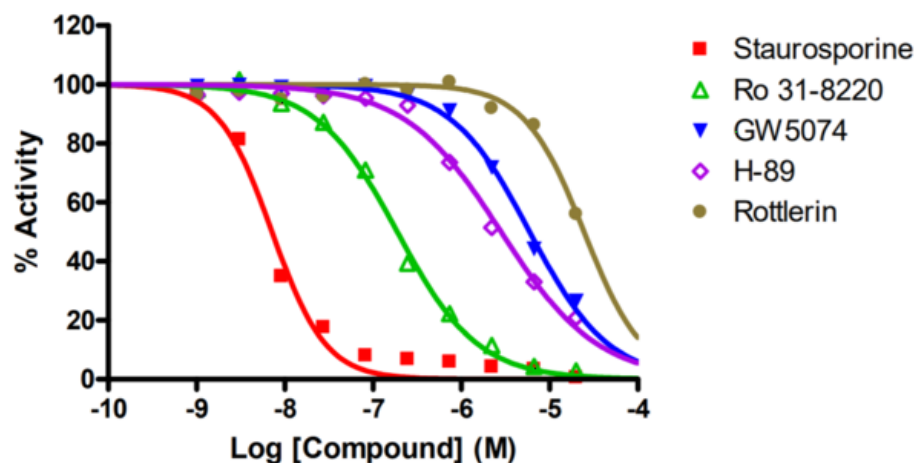
Target Overview

CAT#	STK33
Enzyme	Human STK33
HGNC	STK33
Synonym	STK33
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>7.2</td> </tr> <tr> <td>Ro 31-8220</td> <td>190</td> </tr> <tr> <td>GW5074</td> <td>6,000</td> </tr> <tr> <td>H-89</td> <td>2,800</td> </tr> <tr> <td>Rottlerin</td> <td>>20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	7.2	Ro 31-8220	190	GW5074	6,000	H-89	2,800	Rottlerin	>20,000
Compounds	IC50 (nM)												
Staurosporine	7.2												
Ro 31-8220	190												
GW5074	6,000												
H-89	2,800												
Rottlerin	>20,000												
Scientific Information	Q9BYT3 (http://www.uniprot.org/uniprot/Q9BYT3)												
Screening Location	Malvern, USA												

Reference compound IC50 for STK33



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/stk33_kin_malvern

