

TSSK3/STK22C Kinase Assay Service

Activity-based TSSK3/STK22C kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

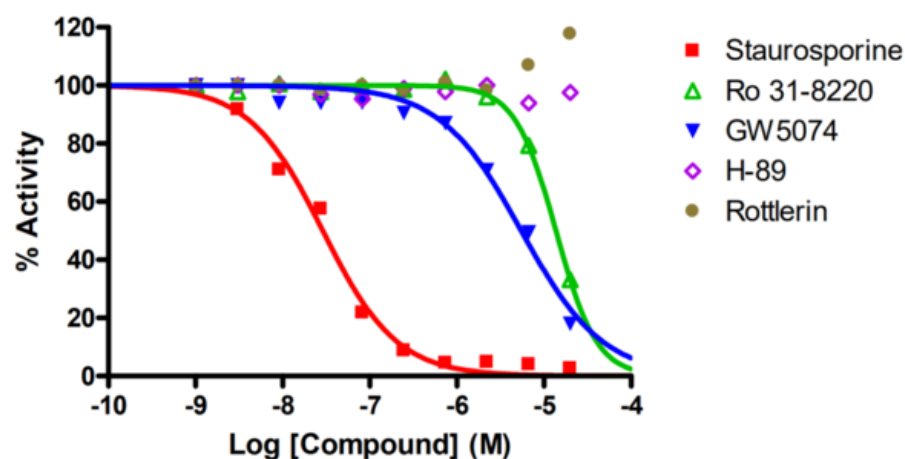
Target Overview

CAT#	TSSK3/STK22C
Enzyme	Human TSSK3/STK22C
HGNC	TSSK3
Synonym	STK22C, STK22D, SPOGA3
Substrates	Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 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>29</td> </tr> <tr> <td>Ro 31-8220</td> <td>14,000</td> </tr> <tr> <td>GW 5074</td> <td>5,500</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	29	Ro 31-8220	14,000	GW 5074	5,500	H-89	>50,000	Rottlerin	>50,000
Compounds	IC50 (nM)												
Staurosporine	29												
Ro 31-8220	14,000												
GW 5074	5,500												
H-89	>50,000												
Rottlerin	>50,000												
Scientific Information	Q96PN8 (http://www.uniprot.org/uniprot/Q96PN8)												
Screening Location	Malvern, USA												

Reference compound IC50 for TSSK3/STK22C



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/tssk3_stk22c_kin_malvern

