

LOK/STK10 Kinase Assay Service

Activity-based LOK/STK10 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

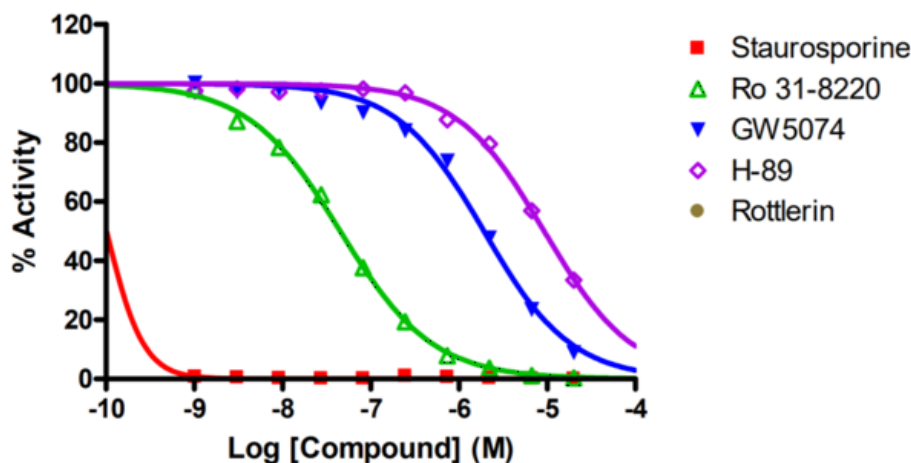
Target Overview

CAT#	LOK/STK10
Enzyme	Human LOK/STK10
HGNC	STK10
Synonym	LOK, PRO2729
Substrates	Peptide substrate, [RLGRDKYKTLRQIRQ], 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><0.5</td> </tr> <tr> <td>Ro 31-8220</td> <td>45</td> </tr> <tr> <td>GW5074</td> <td>1,900</td> </tr> <tr> <td>H-89</td> <td>9,300</td> </tr> <tr> <td>Rottlerin</td> <td>>100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	<0.5	Ro 31-8220	45	GW5074	1,900	H-89	9,300	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	<0.5												
Ro 31-8220	45												
GW5074	1,900												
H-89	9,300												
Rottlerin	>100,000												
Scientific Information	O94804 (http://www.uniprot.org/uniprot/O94804)												
Screening Location	Malvern, USA												

Reference compound IC50 for LOK/STK10



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/lok_stk10_kin_malvern

