

MLCK/MYLK Kinase Assay Service

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

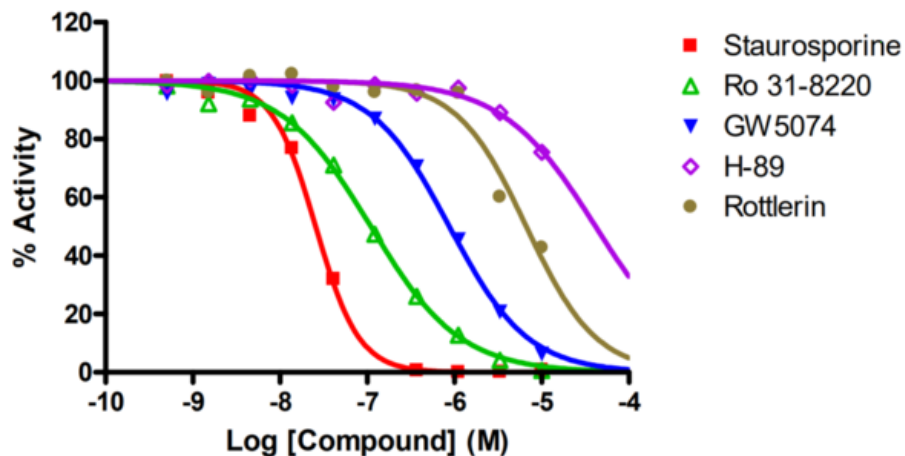
Target Overview

CAT#	MLCK
Enzyme	Human MLCK
HGNC	MYLK
Synonym	smMLCK, FLJ12216, KRP, MLCK, MLCK108, MLCK210
Substrates	Peptide substrate, [KKLNRTLFAEPG], 20 μM Ca ²⁺ -Calmodulin, 1 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>25.6</td> </tr> <tr> <td>Ro 31-8220</td> <td>110</td> </tr> <tr> <td>GW 5074</td> <td>876</td> </tr> <tr> <td>H-89</td> <td>41,600</td> </tr> <tr> <td>Rottlerin</td> <td>7,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	25.6	Ro 31-8220	110	GW 5074	876	H-89	41,600	Rottlerin	7,000
Compounds	IC50 (nM)												
Staurosporine	25.6												
Ro 31-8220	110												
GW 5074	876												
H-89	41,600												
Rottlerin	7,000												
Scientific Information	Q15746 (http://www.uniprot.org/uniprot/Q15746)												
Screening Location	Malvern, USA												

Reference compound IC50 for MLCK/MYLK



Let's Discover Together.

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

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

