

DYRK3 Kinase Assay Service

Activity-based DYRK3 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

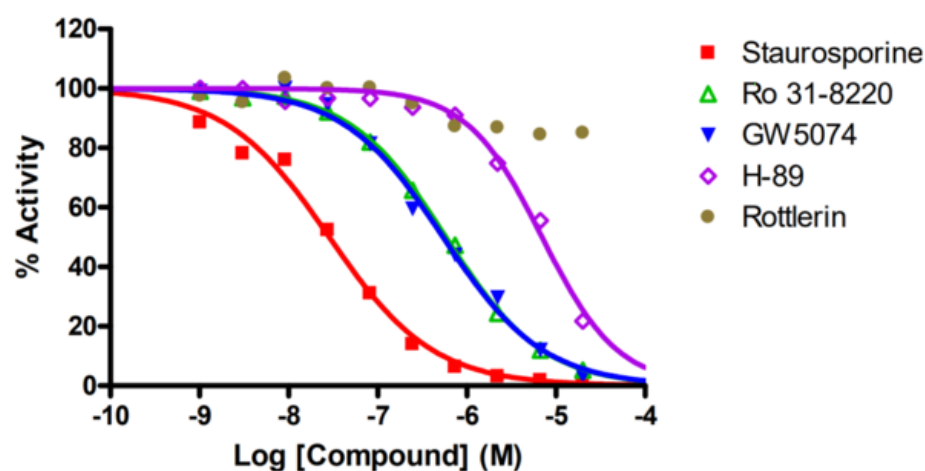
Target Overview

CAT#	DYRK3
Enzyme	Human DYRK3
HGNC	DYRK3
Synonym	DYRK5, RED, REDK
Substrates	Peptide substrate, [RRRFRPASPLRGPPK], 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>580</td> </tr> <tr> <td>GW 5074</td> <td>540</td> </tr> <tr> <td>H-89</td> <td>7,000</td> </tr> <tr> <td>Rottlerin</td> <td>>20,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	29	Ro 31-8220	580	GW 5074	540	H-89	7,000	Rottlerin	>20,000
Compounds	IC50 (nM)												
Staurosporine	29												
Ro 31-8220	580												
GW 5074	540												
H-89	7,000												
Rottlerin	>20,000												
Scientific Information	O43781 (http://www.uniprot.org/uniprot/O43781)												
Screening Location	Malvern, USA												

Reference compound IC50 for DYRK3



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/dyrk3_kin_malvern

