

CDK16/cyclin Y Kinase Assay Service

Activity-based CDK16/cyclin Y kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

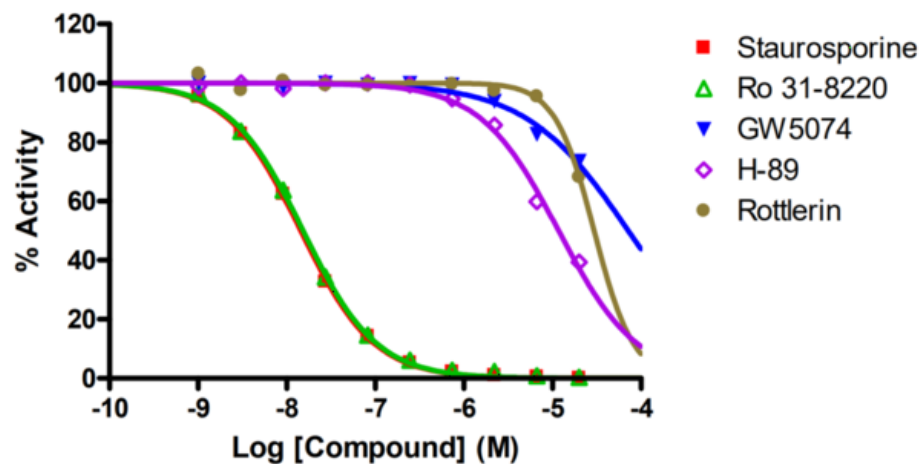
Target Overview

CAT#	CDK16/Cyclin Y
Enzyme	Human CDK16/Cyclin Y
HGNC	CDK16/CCNY
Synonym	PCTAIRE1, PCTAIRE, PCTGAIRE, PCKT1
Substrates	RB protein, 3 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>14</td> </tr> <tr> <td>Ro 31-8220</td> <td>15.3</td> </tr> <tr> <td>GW5074</td> <td>72,000</td> </tr> <tr> <td>H-89</td> <td>11,800</td> </tr> <tr> <td>Rottlerin</td> <td>30,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	14	Ro 31-8220	15.3	GW5074	72,000	H-89	11,800	Rottlerin	30,000
Compounds	IC50 (nM)												
Staurosporine	14												
Ro 31-8220	15.3												
GW5074	72,000												
H-89	11,800												
Rottlerin	30,000												
Scientific Information	CDK16: Q00536 (http://www.uniprot.org/uniprot/Q00536) Cyclin Y: Q8ND76 (http://www.uniprot.org/uniprot/Q8ND76)												
Screening Location	Malvern, USA												

Reference compound IC50 for CDK16/Cyclin Y



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/cdk16_cyclin_y_kin_malvern

