

NEK1 Kinase Assay Service

Activity-based NEK1 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

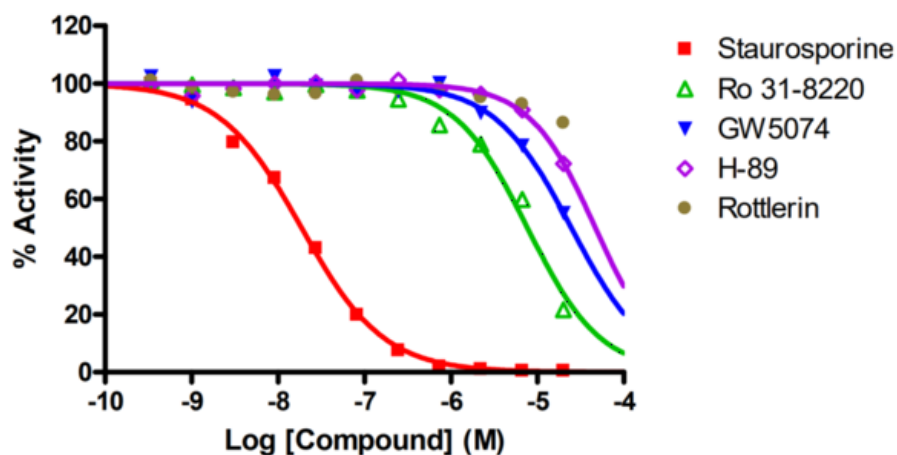
Target Overview

CAT#	NEK1
Enzyme	Human NEK1
HGNC	NEK1
Synonym	KIAA1901, NY-REN-55,
Substrates	MBP (Myelin Basic Protein), 10 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>18.7</td> </tr> <tr> <td>Ro 31-8220</td> <td>7,700</td> </tr> <tr> <td>GW 5074</td> <td>24,700</td> </tr> <tr> <td>H-89</td> <td>47,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	18.7	Ro 31-8220	7,700	GW 5074	24,700	H-89	47,000
Compounds	IC50 (nM)										
Staurosporine	18.7										
Ro 31-8220	7,700										
GW 5074	24,700										
H-89	47,000										
Scientific Information	Q96PY6 (http://www.uniprot.org/uniprot/Q96PY6)										
Screening Location	Malvern, USA										

Reference compound IC50 for NEK1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/nek1_kin_malvern

