

# PKCnu/PRKD3 Kinase Assay Service

Activity-based PKCnu/PRKD3 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

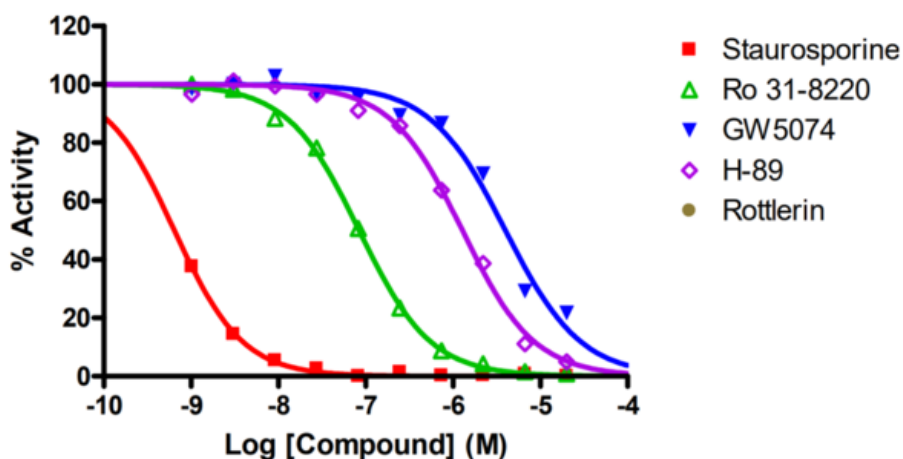
## Target Overview

<b>CAT#</b>	PKCnu/PRKD3
<b>Enzyme</b>	Human PKCnu/PRKD3
<b>HGNC</b>	PRKD3
<b>Synonym</b>	PKD3, EPK2, PKc-NU, PKD3, PRKCN
<b>Substrates</b>	Peptide substrate, [KKLNRTLSVA], 20 μM ATP 10 μM

## Assay Properties

<b>Reaction</b>	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> P-Substrate + ADP										
<b>Readout</b>	Enzymatic activity										
<b>Reference Compound IC50s</b>	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Staurosporine</td> <td>0.6</td> </tr> <tr> <td>Ro 31-8220</td> <td>84</td> </tr> <tr> <td>GW 5074</td> <td>4,000</td> </tr> <tr> <td>H-89</td> <td>1,300</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	0.6	Ro 31-8220	84	GW 5074	4,000	H-89	1,300
Compounds	IC50 (nM)										
Staurosporine	0.6										
Ro 31-8220	84										
GW 5074	4,000										
H-89	1,300										
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/O94806">O94806</a> ( <a href="http://www.uniprot.org/uniprot/O94806">http://www.uniprot.org/uniprot/O94806</a> )										
<b>Screening Location</b>	Malvern, USA										

## Reference compound IC50 for PKCnu/PRKD3



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pkcnu\\_prkd3\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/pkcnu_prkd3_kin_malvern)

