

# PKAcb Kinase Assay Service

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

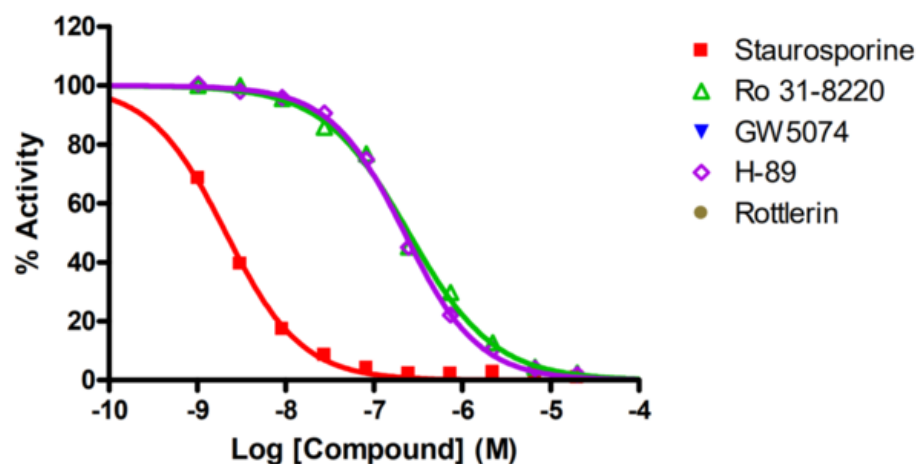
## Target Overview

<b>CAT#</b>	PKAcb
<b>Enzyme</b>	Human PKAcb
<b>HGNC</b>	PRKACB
<b>Synonym</b>	PKAcb, MGC41879, MGC9320, PKA c-beta
<b>Substrates</b>	Peptide substrate, [KEAKEKRQEIQAKRRRLSSLRASTSKSGGSQK], 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>2.1</td> </tr> <tr> <td>Ro 31-8220</td> <td>250</td> </tr> <tr> <td>H-89</td> <td>220</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	2.1	Ro 31-8220	250	H-89	220
Compounds	IC50 (nM)								
Staurosporine	2.1								
Ro 31-8220	250								
H-89	220								
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P22694">P22694</a> ( <a href="http://www.uniprot.org/uniprot/P22694">http://www.uniprot.org/uniprot/P22694</a> )								
<b>Screening Location</b>	Malvern, USA								

## Reference compound IC50 for PKAcb



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pkacb\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/pkacb_kin_malvern)

