

# PAK6 Kinase Assay Service

Activity-based PAK6 kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

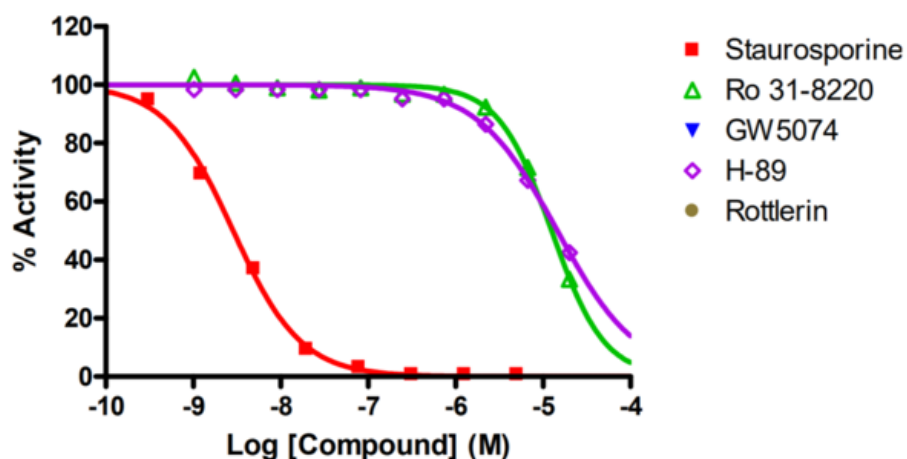
## Target Overview

CAT#	PAK6
Enzyme	Human PAK6
HGNC	PAK6
Synonym	PAK6, PAK5
Substrates	Peptide substrate, [RRRLSFAEPG], 20 μM ATP 10 μM

## Assay Properties

Reaction	Substrate + [ <sup>32</sup> P]-ATP = <sup>32</sup> 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>2.9</td> </tr> <tr> <td>Ro 31-8220</td> <td>12,500</td> </tr> <tr> <td>GW5074</td> <td>&gt;100,000</td> </tr> <tr> <td>H-89</td> <td>14,600</td> </tr> <tr> <td>Rottlerin</td> <td>&gt;100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	2.9	Ro 31-8220	12,500	GW5074	>100,000	H-89	14,600	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	2.9												
Ro 31-8220	12,500												
GW5074	>100,000												
H-89	14,600												
Rottlerin	>100,000												
Scientific Information	<a href="http://www.uniprot.org/uniprot/Q9NQU5">Q9NQU5</a> ( <a href="http://www.uniprot.org/uniprot/Q9NQU5">http://www.uniprot.org/uniprot/Q9NQU5</a> )												
Screening Location	Malvern, USA												

## Reference compound IC50 for PAK6



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/pak6\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/pak6_kin_malvern)

