

# ERN1 (R727A) Kinase Assay Service

Activity-based ERN1 (R727A) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

## Target Overview

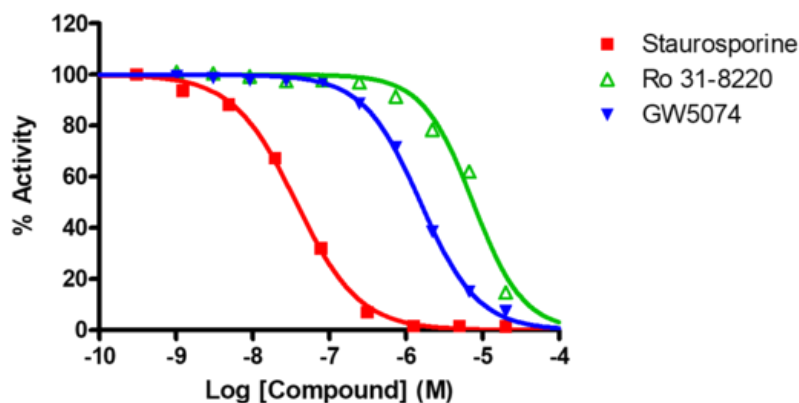
<b>CAT#</b>	ERN1/IRE1 (R727A)
<b>Enzyme</b>	Human ERN1/IRE1 (R727A)
<b>HGNC</b>	ERN1 (R727A)
<b>Synonym</b>	Serine/threonine-protein kinase/endoribonuclease IRE1, Endoplasmic reticulum-to-nucleus signaling 1, Inositol-requiring protein 1
<b>Substrates</b>	MBP (Myelin Basic Protein), 20 μM ATP 10 μM

## Assay Properties

<b>Reaction</b>	Substrate + [ <sup>33</sup> P]-ATP = <sup>33</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>37</td> </tr> <tr> <td>Ro 31-8220</td> <td>7,500</td> </tr> <tr> <td>GW5074</td> <td>1,600</td> </tr> <tr> <td>H-89</td> <td>&gt;100,000</td> </tr> <tr> <td>Rottlerin</td> <td>&gt;100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	37	Ro 31-8220	7,500	GW5074	1,600	H-89	>100,000	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	37												
Ro 31-8220	7,500												
GW5074	1,600												
H-89	>100,000												
Rottlerin	>100,000												
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/O75460">O75460</a> ( <a href="http://www.uniprot.org/uniprot/O75460">http://www.uniprot.org/uniprot/O75460</a> )												
<b>Screening Location</b>	Malvern, USA												

## Reference compound IC50 for ERN1/IRE1 (R727A)

Reference compound IC50 for ERN1/IRE1 (R727A)



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/ern1\\_r727a\\_kin\\_malvern](https://www.reactionbiology.com/datasheet/ern1_r727a_kin_malvern)

