

GSK3b Kinase Assay Service

Activity-based GSK3b kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

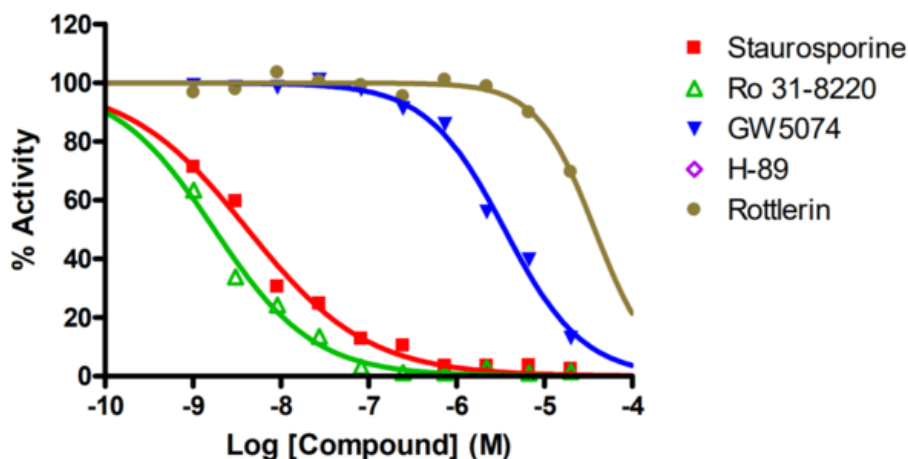
Target Overview

| | |
|------------|--|
| CAT# | GSK3b |
| Enzyme | Human GSK3b |
| HGNC | GSK3B |
| Synonym | GSK3beta |
| Substrates | Peptide substrate, [YRRAAVPPSPSLSRHSSPHQ(pS)EDEEE], 20 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>4.2</td> </tr> <tr> <td>Ro 31-8220</td> <td>1.8</td> </tr> <tr> <td>GW 5074</td> <td>3,500</td> </tr> <tr> <td>H-89</td> <td>>100,000</td> </tr> <tr> <td>Rottlerin</td> <td>37,400</td> </tr> </tbody> </table> | Compounds | IC50 (nM) | Staurosporine | 4.2 | Ro 31-8220 | 1.8 | GW 5074 | 3,500 | H-89 | >100,000 | Rottlerin | 37,400 |
| Compounds | IC50 (nM) | | | | | | | | | | | | |
| Staurosporine | 4.2 | | | | | | | | | | | | |
| Ro 31-8220 | 1.8 | | | | | | | | | | | | |
| GW 5074 | 3,500 | | | | | | | | | | | | |
| H-89 | >100,000 | | | | | | | | | | | | |
| Rottlerin | 37,400 | | | | | | | | | | | | |
| Scientific Information | P49841 (http://www.uniprot.org/uniprot/P49841) | | | | | | | | | | | | |
| Screening Location | Malvern, USA | | | | | | | | | | | | |

Reference compound IC50 for GSK3b



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/gsk3b_kin_malvern

