

LRRK2 (G2019S) Kinase Assay Service

Activity-based LRRK2 (G2019S) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

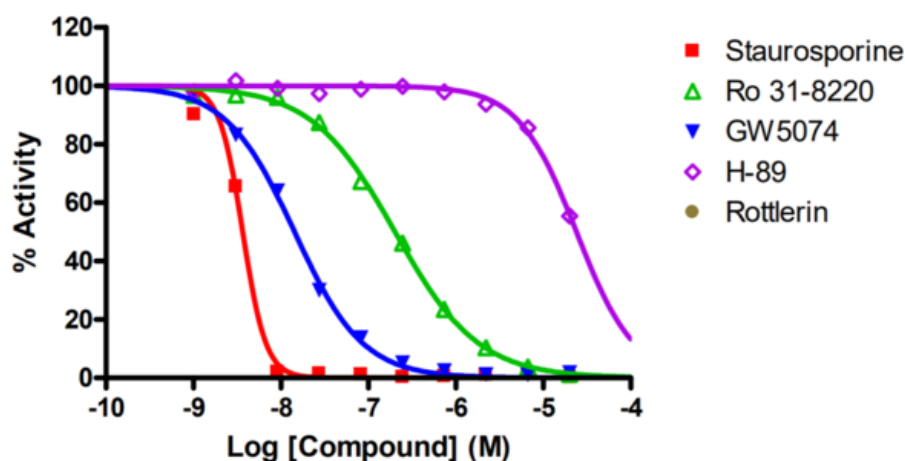
Target Overview

CAT#	LRRK2 (G2019S)
Enzyme	Human LRRK2 (G2019S)
HGNC	LRRK2 (G2019S)
Synonym	PARK8, ROCO2, FLJ45829, DKFZp434H2111
Substrates	Peptide substrate, [RLGRDKYKTLRQIRQ], 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>3.7</td> </tr> <tr> <td>Ro 31-8220</td> <td>202</td> </tr> <tr> <td>GW5074</td> <td>14</td> </tr> <tr> <td>H-89</td> <td>24,000</td> </tr> <tr> <td>Rottlerin</td> <td>>100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	3.7	Ro 31-8220	202	GW5074	14	H-89	24,000	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	3.7												
Ro 31-8220	202												
GW5074	14												
H-89	24,000												
Rottlerin	>100,000												
Scientific Information	Q5S007 (http://www.uniprot.org/uniprot/Q5S007)												
Screening Location	Malvern, USA												

Reference compound IC50 for LRRK2 (G2019S)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/lrrk2_g2019s_kin_malvern

