

BRAF-KIAA1549 (Kex16Bex9) Kinase Assay Service

Activity-based BRAF-KIAA1549 (Kex16Bex9) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

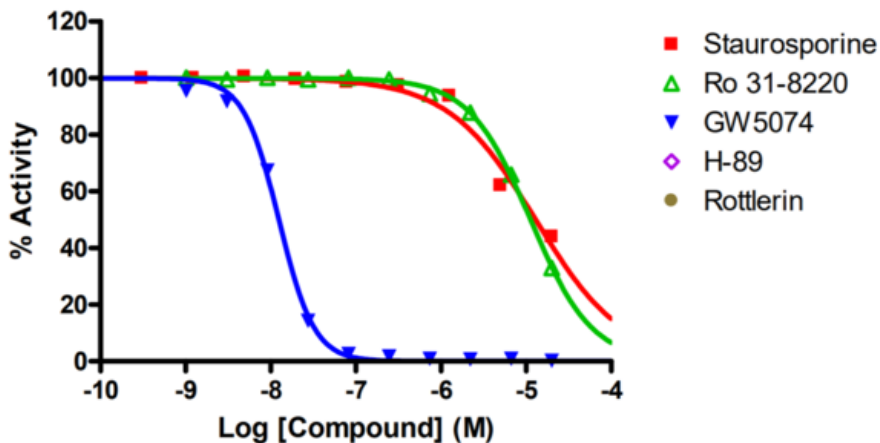
Target Overview

CAT#	BRAF, KIAA1549-
Enzyme	Human KIAA1549-BRAF fusion
HGNC	BRAF
Synonym	B-Raf, B-raf-1, BRAF1, RAFB1
Substrates	MEK1 (K97R), 5 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>13,000</td> </tr> <tr> <td>Ro 31-8220</td> <td>11,300</td> </tr> <tr> <td>GW5074</td> <td>13</td> </tr> <tr> <td>H-89</td> <td>>100,000</td> </tr> <tr> <td>Rottlerin</td> <td>>100,000</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Staurosporine	13,000	Ro 31-8220	11,300	GW5074	13	H-89	>100,000	Rottlerin	>100,000
Compounds	IC50 (nM)												
Staurosporine	13,000												
Ro 31-8220	11,300												
GW5074	13												
H-89	>100,000												
Rottlerin	>100,000												
Scientific Information	KIAA1549: Q9HCM3 (http://www.uniprot.org/uniprot/Q9HCM3) BRAF: P15056 (http://www.uniprot.org/uniprot/P15056)												
Screening Location	Malvern, USA												

Reference compound IC50 for BRAF KIAA1549 (Kex16Bex9)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/braf-kiaa1549_kex16bex9_kin_malvern

