

BRAF (V599E) Kinase Assay Service

Activity-based BRAF (V599E) kinase assay for compound screening and profiling via [radiometric HotSpot™ kinase assay](#).

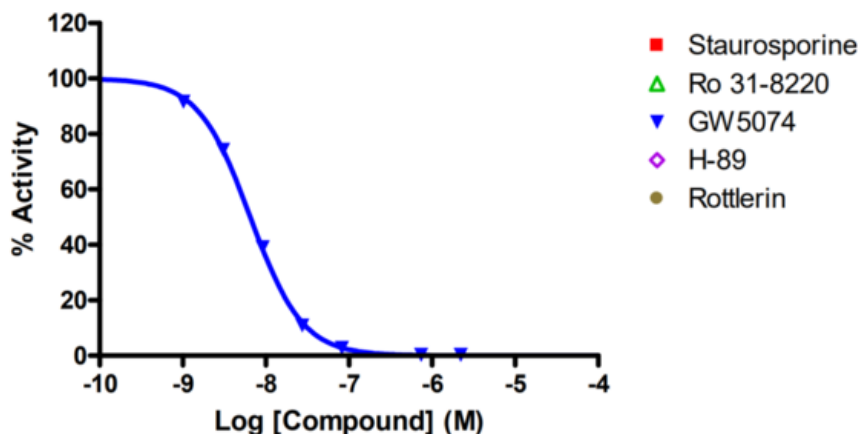
Target Overview

CAT#	BRAF (V599E)
Enzyme	Human BRAF (V599E)
HGNC	BRAF (V600E)
Synonym	B-Raf, B-raf-1, BRAF1, RAFB1
Substrates	MEK1 (K97R), 1 μM ATP 10 μM

Assay Properties

Reaction	Substrate + [³³ P]-ATP = ³³ P-Substrate + ADP
Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) GW 5074 8.5
Scientific Information	P15056 (http://www.uniprot.org/uniprot/P15056)
Screening Location	Malvern, USA
Further information	More information can be found on our website Kinase Assays .

Reference compound IC50 for BRAF (V599E)



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/braf_v599e_kin_malvern

