

Dabrafenib

B-Raf(mutant V600E) inhibitor

Catalog # B08-901

Lot # D3370-15

CAS # 1195765-45-7

Product Description

Molecular Formula: C₂₃H₂₀F₃N₅O₂S₂

Appearance: Off-White powder

Molecular Weight: 519.6

Purity: >98% (TLC); NMR (conforms)

Solubilization: May be dissolved in DMSO (30 mg/ml); or ethanol (1 mg/ml, warm)

Alias

N-[3-[5-(2-amino-4-pyrimidinyl)-2-(1,1-dimethylethyl)-4-thiazolyl]-2-fluorophenyl]-2,6-difluorobenzenesulfonamide; GSK2118436

Specific Activity

B-Raf(mutant V600E) inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 1 year.

Store solutions at -20°C for up to 3 months.

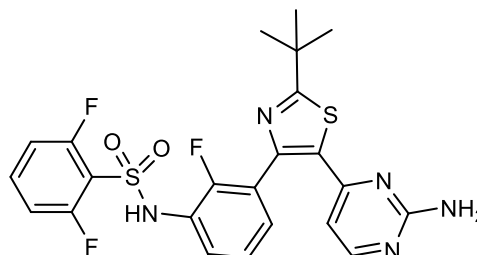
Scientific Background

Selective inhibitor of mutant B-Raf^{V600E} (IC₅₀ = 0.8 nM), with 4- and 6-fold reduced potency against wild type B-Raf and c-Raf (IC₅₀ = 3.2 and 5 nM, respectively).¹ Has been in clinical trials in patients with B-Raf^{V600E} metastatic melanoma and other solid tumors

References

1. J Huang *et al.* *J. Hematol. Oncol.* 2013 6:1
2. C Ji *et al.* *Drug Des. Dev. Ther.* 2016 10:2491
3. R Herr *et al.* *Cancer Res.* 2015 75:216

Molecular Structure



Dabrafenib

B-Raf(mutant V600E) inhibitor

Catalog #	B08-901
CAS #	1195765-45-7
Lot #	D3370-15
Purity	>98% (TLC); NMR (conforms)
Format	Off-White powder
Solubilization	May be dissolved in DMSO (30 mg/ml); or Ethanol (1 mg/ml) warm
Storage & Shipping	Store desiccated as supplied at -20°C for up to 1 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.