

Dabrafenib

B-Raf(mutant V600E) inhibitor

Catalog # B08-901

Lot # T4298-6

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.

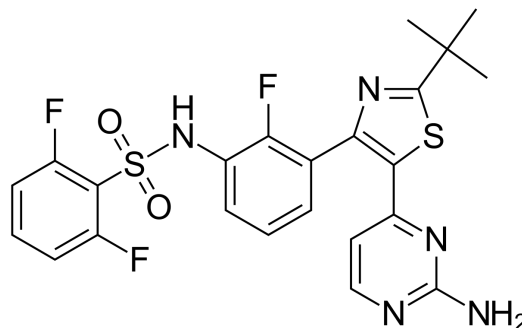
Scientific Background

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 #	T4298-6
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.

SAFETY DATA SHEET

Article 1 – Identification of the Substance

Product Name: Dabrafenib

Catalog # B08-901

This product is sold only for research use by qualified laboratory personnel, and is not to be used as a drug, medical device, food additive, cosmetic, nor household chemical. It is not to be used in diagnostic, therapeutic, consumer, agricultural, nor pesticidal applications.

Manufacturer's Name: SignalChem Biotech Inc.
Street Address: 110-13120 Vanier Place
City, Prov. Postal Code: Richmond, BC, V6V 2J2
Fax: 604-232-4601
EMERGENCY PHONE: 604-232-4600

Article 2 - Hazardous Identification

2.1 Classification of Substance (GHS)

Repr. Category 2
Aquatic Acute, Category 1
Aquatic Chronic, Category 2

2.2 Label Elements

Classification according to Regulation (EC) No 1272/2008



GHS09



GHS08

GHS Signal Word: WARNING

GHS Hazard Phrases:

H361	Suspected of damaging fertility to the unborn child.
H373	May cause damage to the hearth through prolonged or repeated exposure.
H400	Very toxic to aquatic life.
H411	Toxic to aquatic life with long-lasting effects.

GHS Precaution Phrases:

P201	Obtain special instructions before use.
P202	Do not handle until all safety precautions have been read and understood.
P260	Do not breathe dust/fume/gas/mist/vapors/spray
P273	Avoid release to the environment.
P280	Wear protective gloves/protective clothing/eye protection/face protection.

GHS Response Phrases:

P308+313	IF exposed or concerned: Get medical attention/advice.
P391	Collect spillage.
P405	Store locked up.

2.3 Other hazards

Data not available

Article 3 – Composition/Information on Ingredients

3.1 Substance

Product Name(s): Dabrafenib
Chemical Name: N-[3-[5-(2-amino-4-pyrimidinyl)-2-(1,1-dimethylethyl)-4-thiazolyl]-2-fluorophenyl]-2,6-difluorobenzenesulfonamide; GSK2118436
Molecular Formula: C₂₃H₂₀F₃N₅O₂S₂
CAS Number: 1195765-45-7

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.

Article 4 – First Aid

4.1 Description of First Aid Measures

General: Symptoms of poisoning may even occur after several hours; therefore, medical observation for at least 48 hours after the accident.

If Inhaled: Supply fresh air; consult doctor in case of complaints.

In Case of Skin Contact: Generally, the product does not irritate the skin.

In Case of Eye Contact: Rinse opened eye for several minutes under running water.

If Swallowed: If symptoms persist, consult a doctor.

4.2 Important Symptoms and Effects: May cause anemia, cough, CNS depression, drowsiness, headache, heart damage, lassitude (weakness, exhaustion), liver damage, narcosis, reproductive effects, teratogenic effects. No further relevant information available.

Article 5 – Fire Fighting Measures

5.1 Extinguishing Media

Use fire fighting measures that suit the environment. A solid water stream may be inefficient.

5.2 Special Hazards

No further relevant information available.

5.3 Precautions for Fire Fighters

No special measures required.

Article 6 – Accidental Release Measures

6.1 Personal Precautions

Not required.

6.2 Environmental Precautions

Inform respective authorities in case of seepage into water course or sewage system. Do not allow to enter sewers/surface or ground water.

6.3 Methods for Containment and Clean-up

Dispose contaminated material as waste according to Article 13.

Article 7 – Handling and Storage

7.1 Handling Precautions

No special precautions are necessary if used correctly. Avoid breathing dust/fume/gas/mist/vapours/spray. Avoid prolonged or repeated exposure. Keep away from sources of ignition. Take precautionary measures against static charge.

7.2 Safe Storage Measures

Keep container tightly closed. Store in accordance with information listed on the product insert.

Article 8 – Exposure Controls/Personal Protection

8.1 Control Parameters

Contains no substances with occupational exposure limit values.

8.2 Specific Engineering Controls

Keep airborne levels low by using a fumehood or local exhaust.

8.3 Personal Protection

Eye/Face Protection: Tightly sealed goggles

Skin Protection: Use chemical resistant gloves. Remove and dispose of gloves properly after use. Wash hands thoroughly after use.

Body Protection: Wear a lab coat.

Respiratory Protection: In case of brief exposure or low pollution, used respiratory filter device. In case of intensive or longer exposure, use respiratory protective device that is independent of circulating air.

Article 9 – Physical and Chemical Properties

Appearance: Off White powder

Odour: data not available

Odour Threshold: data not available

pH: data not available

Melting/Freezing Pt: Data not available

Boiling Pt: data not available

Flash Pt: data not available

Explosive Properties: data not available

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

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.

Article 10 – Stability and Reactivity

10.1 Stability: No decomposition if used according to specifications.

10.2 Conditions to Avoid: No relevant information available.

10.3 Incompatibility – Materials to Avoid: Strong oxidizing agents.

10.4 Hazardous Decomposition Products: Carbon oxides, hydrogen fluoride, hydrogen sulfide, nitrogen oxides.

Article 11– Toxicological Information

11.1 Toxicological Effects: The toxicological effects of this product have not been thoroughly studied.

Acute Toxicity: None known.

Chronic Toxicity: None known.

Skin Corrosion/Irritation: No irritant effect.

Eye Damage/Irritation: No irritant effect.

Respiratory or Skin Sensitization: No sensitizing effects known.

Germ Cell Mutagenicity: data not available

Carcinogenicity: None according to NIOSH

Article 12 – Ecological Information

12.1 Toxicity

Avoid release into the environment. Very toxic for fish.

12.2 Degradability

Data not available

Article 13 – Disposal Considerations

13.1 Waste Treatment

Must not be disposed of together with household garbage. Do not allow product to reach sewage system. Disposal must be made according to official regulations.

Article 14 – Transport Information

14.1 DOT

Not dangerous goods

14.2 IMDG

Not dangerous goods

14.2 IATA

Not dangerous goods

Article 15 – Regulatory Information

This product has been classified in accordance with the hazard criteria of the Controlled Products Regulations and the SDS contains all the information required by the Controlled Products Regulations.

Article 16 – Other Information

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. SignalChem shall not be held liable for any damage resulting from handling or from contact with the above product. See the Technical Specification, Packing Slip, Invoice, and Product Catalog for additional terms and conditions of sale.

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.