

## VX-702

p38 MAPK Inhibitor

### Catalog # M39-900D

Lot # J3942-25

CAS # 745833-23-2

### Product Description

Molecular Formula: C<sub>19</sub>H<sub>12</sub>F<sub>4</sub>N<sub>4</sub>O<sub>2</sub>

Appearance: White powder

Molecular Weight: 404.3

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

Solubility: May be dissolved DMSO (40 mg/ml); or ethanol (2 mg/ml)

### Alias

6-[(Aminocarbonyl)(2,6-difluorophenyl)amino]-2-(2,4-difluorophenyl)-3-pyridinecarboxamide

### Specific Activity

p38 MAPK inhibitor

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

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

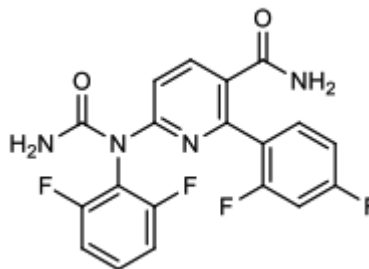
### Scientific Background

A potent and selective inhibitor of p38 MAP kinases, p38 $\alpha$  MAPK and p38 $\beta$  (K<sub>D</sub> = 3.7 and 17 nM respectively) (1). Inhibits IL-6, IL-1 $\beta$  and TNF $\alpha$  production in LPS-primed blood (1). Inhibits p38 in platelets stimulated by thrombin, collagen or thromboxane agonist U-46619 at 1  $\mu$ M (2). Provides only modest clinical efficacy in rheumatoid arthritis (3, 4). Delays platelet lesions and leads to better maintenance of stored platelets (5).

### References

1. Goldstein DM et al. J. Med. Chem. 2010 53:2345
2. Kuliopulos A et al. Thromb. Haemostasis 2004 92:1387
3. Damjanov N et al. Arthrit. Rheumat. 2009 60:1232
4. Ding C Curr. Opin. Investig. Drugs 2006 7:1020
5. Scripchenko A et al. PLoS One 2013 8(8)e:70732

### Molecular Structure



## VX-702

p38 MAPK Inhibitor

Catalog #	M39-900D
CAS #	745833-23-5
Lot #	J3942-25
Purity	>98% (HPLC); NMR (conforms)
Format	White powder
Solubility	May be dissolved in DMSO (40 mg/ml); or ethanol (2 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://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: VX-702**

**Catalog # M39-900D**

*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)

None

### 2.2 Label Elements

Classification according to Regulation (EC) No 1272/2008

None

**GHS Signal Word: None**

**GHS Hazard Phrases: None**

**GHS Precaution Phrases: None**

**GHS Response Phrases: None**

**GHS Storage and Disposal Phrases: None**

Please refer to Section 7 for Storage and Section 13 for Disposal Information

**2.3 Other hazards:** To the best of our knowledge, the toxicological properties have not been thoroughly investigated.

## Article 3 – Composition/Information on Ingredients

### 3.1 Substance

**Product Name(s):** VX-702  
**Chemical Name:** 6-[(Aminocarbonyl) (2,6-difluorophenyl)amino]-2-(2,4-difluorophenyl)-3-pyridinecarboxamide  
**Molecular Formula:** C<sub>19</sub>H<sub>12</sub>F<sub>4</sub>N<sub>4</sub>O<sub>2</sub>  
**CAS Number:** 745833-23-5  
**Concentration:** 100%

## Article 4 – First Aid

### 4.1 Description of First Aid Measures

**General:** Move out of dangerous area and contact a doctor. Provide them with this data sheet.

**If Inhaled:** Move to fresh air, and monitor breathing. Give artificial respiration if breathing stops. Get immediate medical attention.

**In Case of Skin Contact:** Remove contaminated clothing. Immediately wash with soap and water for an extended period of time. Consult a doctor if symptoms occur. Wash clothing before reuse.

**In Case of Eye Contact:** Hold eyelids apart and immediately flush eye(s) with copious amounts of water for at least 15 minutes. Have eyes examined and tested by medical personnel.

**If Swallowed:** Rinse out the mouth with water provided person is conscious. Do not give anything by mouth. Get medical attention. Do NOT induce vomiting unless directed to do so by medical personnel.

**4.2 Important Symptoms and Effects:** To the best of our knowledge, the chemical, physical and toxicological properties have not been thoroughly investigated.

## Article 5 – Fire Fighting Measures

### 5.1 Extinguishing Media

Use water spray, alcohol-resistant foam, dry chemical or carbon dioxide.

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

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

## 5.2 Special Hazards

In combustion, may emit toxic fumes.

## 5.3 Precautions for Fire Fighters

Wear suitable protection to prevent contact and respirator for fumes, and full protective gear to prevent contact with skin and eyes.

---

## Article 6 – Accidental Release Measures

---

### 6.1 Personal Precautions

Relocate personnel to a safe area and ensure ventilation is in working order. Wear protective equipment. Avoid dust formation and inhalation of fumes. As conditions warrant, wear an approved self-contained breathing apparatus, or respirator, and appropriate personal protection (rubber boots, safety goggles, and heavy rubber gloves).

### 6.2 Environmental Precautions

Do not let product enter drains, if safe to do so.

### 6.3 Methods for Containment and Clean-up

Contain spill and collect, as appropriate. Sweep up material without generating dust clouds and place in a sealed container for disposal. Clean surface where spill occurred.

---

## Article 7 – Handling and Storage

---

### 7.1 Handling Precautions

Use in a well ventilated area. Avoid formation of dust or aerosols. Keep away from sources of ignition. Wear suitable protective clothing. Avoid inhalation, and skin or eye contact.

### 7.2 Safe Storage Measures

Keep product in a sealed contained located in dry well ventilated area. 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:** Use safety glasses.

**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:** If risk is present wear a dust respirator.

---

## Article 9 – Physical and Chemical Properties

---

**Appearance:** 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 (40 mg/ml); or ethanol (2 mg/ml)

---

## Article 10 – Stability and Reactivity

---

**10.1 Stability:** Stable under recommended transport or storage conditions

**10.2 Conditions to Avoid:** Moist and hot conditions

**10.3 Incompatibility – Materials to Avoid:** Strong acids/alkalis, strong oxidizing/reducing agents

**10.4 Hazardous Decomposition Products:** Possible toxic fumes upon combustion.

---

## Article 11– Toxicological Information

---

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

**Acute Toxicity:** No data available

**Skin Corrosion/Irritation:** data not available

**Eye Damage/Irritation:** data not available

**Respiratory or Skin Sensitization:** data not available

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

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

**Germ Cell Mutagenicity:** data not available  
**Carcinogenicity:** None according to NIOSH

---

## Article 12 – Ecological Information

---

### 12.1 Toxicity

No data available.

### 12.2 Degradability

Data not available

---

## Article 13 – Disposal Considerations

---

### 13.1 Waste Treatment

Product and Contaminated Packaging: Dispose according to national legislation. Contact a licensed chemical disposal company.

---

## 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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