

Flavopiridol HCl

CDK Inhibitor

Catalog # C22-900

Lot # J3942-18

CAS # 131740-09-5

Product Description

Molecular Formula: C₂₁H₂₀ClNO₅ · HCl

Appearance: Pale-yellow powder

Molecular Weight: 438.3

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

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

Alias

2-(2-Chlorophenyl)-5,7-dihydroxy-8-[(3S, 4R)-3-hydroxy-1-methyl-4-piperidiny]-4H-1-benzopyran-4-one, hydrochloride; L 86-8275; Alvocidib; NSC-649890

Specific Activity

CDK Inhibitor

Storage and Stability

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

Store solutions at -20°C for up to 1 month.

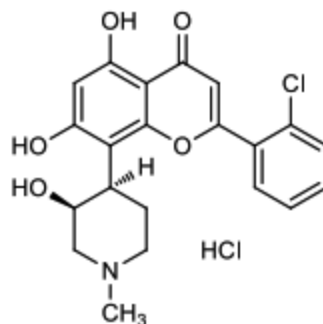
Scientific Background

Cyclin-dependent kinase (CDK) inhibitor that causes cell cycle arrest at G₁ and G₂ phase (1). Potentiates (100-200 nM) PMA-induced apoptosis in HL-60 and U937 cells (2). Potentiates the effects of topoisomerase I inhibitors by suppressing Rad51 expression in a p53-dependent manner. Displays potent antiviral activity against HIV-1 and HSV-1 *in vitro* (4).

References

1. Kaur G et al. J. Natl. Cancer Inst. 1992 84:1736
2. Cartee L et al. Mol. Pharmacol. 2002 61:1313
3. Ambrosini G et al. Cancer Res. 2008 68:2312
4. Schang LM et al. Biochim. Biophys. Acta 2004 1697:197

Molecular Structure



Flavopiridol HCl

CDK Inhibitor

Catalog #	C22-900
CAS #	131740-09-5
Lot #	J3942-18
Purity	>98% (HPLC); NMR (Conforms)
Format	Pale-yellow powder
Solubility	May be dissolved DMSO (40 mg/ml); or water (40 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

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: Flavopiridol HCl

Catalog # C22-900

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)

Acute toxicity, Oral, Category 4

2.2 Label Elements

Classification according to Regulation (EC) No 1272/2008



GHS Signal Word: WARNING

GHS Hazard Phrases:

H302	Harmful if swallowed
------	----------------------

GHS Precaution Phrases:

P264	Wash hands thoroughly after handling.
P270	Do not eat, drink or smoke when using this product.

GHS Response Phrases:

P301+P312	IF SWALLOWED: Call a POISON CENTRE/doctor if you feel unwell
P330	Rinse mouth.
P501	Dispose of contents/container to an approved waste disposal plant.

GHS Storage and Disposal Phrases:

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): Flavopiridol HCl
Chemical Name: 2-(2-Chlorophenyl)-5,7-dihydroxy-8-[(3S, 4R)-3-hydroxy-1-methyl-4-piperidinyl]-4H-1-benzopyran-4-one, hydrochloride
Molecular Formula: C₂₁H₂₀ClNO₅ · HCl
CAS Number: 131740-09-5
Concentration: 100%

Article 4 – First Aid

4.1 Description of First Aid Measures

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. Get medical attention. Do NOT induce vomiting unless directed to do so by medical personnel.

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.

4.2 Important Symptoms and Effects: The most important known symptoms and effects are described in section 2.2.

Article 5 – Fire Fighting Measures

5.1 Extinguishing Media

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

5.2 Special Hazards

During combustion, may emit irritant 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: Pale-yellow 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 DMSO (40 mg/ml); or water (40 mg/ml)

Article 10 – Stability and Reactivity

10.1 Stability: Stable under recommended transport or storage conditions

10.2 Conditions to Avoid: No data available

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

10.4 Hazardous Decomposition Products: Possible toxic fumes upon combustion.

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 11– Toxicological Information

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

Carcinogenicity: None according to NIOSH

Article 12 – Ecological Information

12.1 Toxicity

Avoid release into the environment.

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
www.signalchem.com

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