

Apigenin

Kinase Inhibitor

Catalog # K10-900C

Lot # S3524-5

CAS # 520-36-5

Product Description

Molecular Formula: C₁₅H₁₀O₅

Appearance: Yellow powder

Melting Point: 345-350°C

Molecular Weight: 270.2

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

Solubilization: May be dissolved in DMSO (50 mg/ml)

Alias

4',5,7-Trihydroxyflavone

Specific Activity

CK2 kinase inhibitor

Storage and Stability

Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 1 month.

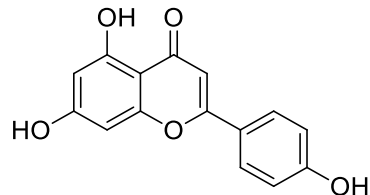
Scientific Background

Selective CK2 inhibitor. Inhibits proliferation of Wnt-1 transfected cells and suppresses phosphorylation of β-catenin. Ameliorates renal dysfunction at 20 mg/Kg in a rat model of glomerulonephritis. Cell permeable and active in vivo.

References

1. J Shen *et al.* *J. Immunol.* 2001 167:4919
2. DH Song *et al.* *J. Biol. Chem.* 2000 275:23790
3. M Yamada *et al.* *Proc. Natl. Acad. Sci. USA* 2005 102:7736

Molecular Structure



Apigenin

Kinase Inhibitor

Catalog #

K10-900C

CAS Number

520-36-5

Lot #

S3524-5

Purity

>98% (TLC); NMR (Conforms)

Format

Lyophilized powder

Solubilization

May be dissolved in DMSO (50 mg/ml)

Storage & Shipping

Store desiccated as supplied at ambient temperature 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
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: Apigenin

Catalog # K10-900C

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
Skin Corrosion/Irritation, Category 2
Serious Eye Damage/Eye Irritation, Category 1
Toxic to Reproduction, Category 2
Specific target organ toxicity – (repeated exposure) Category 2

2.2 Label Elements

Classification according to Regulation (EC) No 1272/2008



GHS Signal Word: DANGER

GHS Hazard Phrases:

H302	Harmful if swallowed
H315	Causes skin irritation.
H318	Causes serious eye damage.
H361	Suspected of damaging fertility or the unborn child
H373	May cause damage to organs through prolonged or repeated exposure.

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
P264	Wash hands thoroughly after handling.
P280	Wear protective gloves/protective clothing/eye protection/face protection.
P270	Do not eat, drink or smoke when using this product

GHS Response Phrases:

P301+312+330	IF SWALLOWED: Call a POISON CENTRE or doctor/physician if you feel unwell. Rinse Mouth
P302+352	IF ON SKIN: Wash with plenty of soap and water.
P305+P351+P338	IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
P308+313	IF exposed or concerned: Get medical attention/advice.
P332+313	If skin irritation occurs, get medical advice/attention.
P362+364	Take off contaminated clothing and wash it before reuse.

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.

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 3 – Composition/Information on Ingredients

3.1 Substance

Product Name(s): Apigenin
Chemical Name: 4',5,7-Trihydroxyflavone
Molecular Formula: C₁₅H₁₀O₅
CAS Number: 520-36-5
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: Lung, skin and eye irritant. Prolonged exposure is not known to aggravate prior medical conditions

Article 5 – Fire Fighting Measures

5.1 Extinguishing Media

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

5.2 Special Hazards

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.

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 9 – Physical and Chemical Properties

Appearance: Yellow powder
Odour: data not available
Odour Threshold: data not available
pH: data not available
Melting/Freezing Pt: 345-350°C
Boiling Pt: data not available
Flash Pt: data not available
Explosive Properties: data not available
Solubility notes: May be dissolved in DMSO (50 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 oxidizing agents
10.4 Hazardous Decomposition Products: Carbon oxides (CO, CO₂). 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: Oral TDLO (mouse): 50 mg/kg
Chronic Toxicity: Investigated as a drug.
For complete information, see Registry of Toxic Effects of Chemical Substances (RTECS) entry. Ruxolitinib RTECS Number: UQ8002700
Skin Corrosion/Irritation: data not available
Eye Damage/Irritation: data not available
Respiratory or Skin Sensitization: data not available
Germ Cell Mutagenicity: data not available
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.