



SC-79

AKT Activator

Catalog # A16-901

Lot # E4297-1

CAS # 305834-79-1

Catalog #

Aliquot Size

A16-901-5

5 mg

A16-901-25

25 mg

Product Description

Molecular Formula: $C_{17}H_{17}ClN_2O_5$

Appearance: White powder

Molecular Weight: 364.80

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

Solubilization: May be dissolved in DMSO (100 mg/ml) or ethanol (40 mg/ml, warm)

Alias

2-Amino-6-chloro- α -cyano-3-(ethoxycarbonyl)-4H-1-benzopyran-4-acetic acid, ethyl ester

Specific Activity

AKT Activator

Storage and Stability

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

Store solutions at -20°C for up to 2 months.

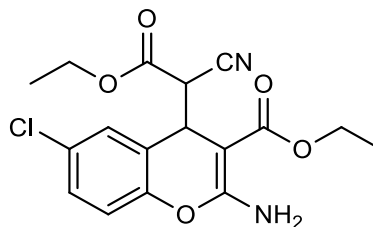
Scientific Background

Akt activator. Binds to the pleckstrin homology domain of Akt preventing membrane translocation but activating it in the cytosol. Enhances Akt phosphorylation by upstream protein kinases. Suppresses excitotoxicity-induced neuronal death *in vitro* and *in vivo* (1). A useful tool for studying PI3K/Akt axis (2,3).

References

1. H Jo *et al.* *Proc. Natl. Acad. Sci. USA* 2012 109:10581
2. W. Yang *et al.* *Cell. Physiol. Biochem.* 2016 38:2366
3. YX Chen *et al.* *Oncotarget* 2017 8:31065

Molecular Structure



SC-79

AKT Activator

Catalog #

A16-901

CAS #

305834-79-1

Lot #

E4297-1

Purity

98% (TLC); NMR (Conforms)

Format

Lyophilized white powder

Solubilization

May be dissolved in DMSO (100 mg/ml)
Ethanol (40 mg/ml, warm)

Storage & Shipping

Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 2 months.

To place your order, please contact us by phone 1-866-9KINASE (54-6273), 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: SC-79

Catalog # A16-901

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)

The substance is not classified, according to the Globally Harmonized System (GHS).

2.2 Label Elements

None

GHS Signal Word: Not applicable

GHS Hazard Phrases: Not applicable

GHS Precaution Phrases: Not applicable

GHS Response Phrases: Not applicable

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): SC-79
Chemical Name: 2-Amino-6-chloro-a-cyano-3-(ethoxycarbonyl)-4H-1-benzopyran-4-acetic acid, ethyl ester
Molecular Formula: C₁₇H₁₇ClN₂O₅
CAS Number: 305834-79-1
Concentration: 100%

Article 4 – First Aid

4.1 Description of First Aid Measures

General: No special measures required.

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 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 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.

To place your order, please contact us by phone 1-866-9KINASE (54-6273), 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 6 – Accidental Release Measures

6.1 Personal Precautions

Not required.

6.2 Environmental Precautions

Do not allow to enter sewers/surface or ground water.

6.3 Methods for Containment and Clean-up

Pick up mechanically.

Article 7 – Handling and Storage

7.1 Handling Precautions

No special measures required.

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: Not required.

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 (100 mg/ml) or ethanol (40 mg/ml, warm)

Article 10 – Stability and Reactivity

10.1 Stability: No further relevant information available.

10.2 Conditions to Avoid: No further relevant information available.

10.3 Incompatibility – Materials to Avoid: Strong oxidizing agents

10.4 Hazardous Decomposition Products: Carbon oxides (CO, CO₂), hydrogen chloride, nitrogen oxides.

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: No irritating effect

Eye Damage/Irritation: No irritating effect.

Respiratory or Skin Sensitization: No sensitizing effects known.

Germ Cell Mutagenicity: data not available

Carcinogenicity: None according to NIOSH

To place your order, please contact us by phone 1-866-9KINASE (54-6273), 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 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

Smaller quantities can be disposed of with household waste. 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 1-866-9KINASE (54-6273), 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.