

Reversine

A3 adenosine receptor antagonist; Aurora A/B/C inhibitor

Catalog # A28-900

Lot # J3942-24

CAS # 656820-32-5

Product Description

Molecular Formula: C₂₁H₂₇N₇O

Appearance: Off-white powder

Molecular Weight: 393.5

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

Solubility: May be dissolved DMSO (7 mg/ml)

Alias

2-(4-Morpholinoanilino)-6-cyclohexylaminopurine

Specific Activity

A3 adenosine receptor antagonist

Storage and Stability

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

Scientific Background

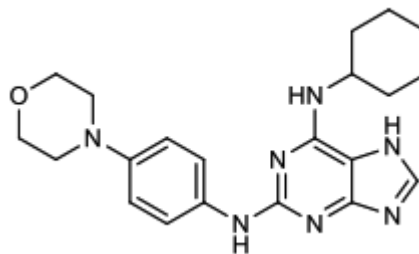
Induces differentiated myogenic-lineage-committed cells to become multipotent mesenchymal progenitor cells (1, 2). Acts as a selective human A₃ adenosine receptor antagonist (K_i=0.66 μM) (3). Cell permeable.

Reversine is also a known ATP-competitive Aurora kinase inhibitor with nanomolar IC₅₀s for Aurora A, Aurora B and Aurora C, as well as a monopolar spindle 1 (Mps1) kinase inhibitor and mitotic spindle assembly checkpoint (SAC) inhibitor.

References

1. Chen S et al. J. Am. Chem. Soc. 2004 126:410
2. Chen S et al. Proc. Natl. Acad. Sci USA 2007 104:10482
3. Pereira M et al. J. Med. Chem. 2005 48:4910
4. D'Alise AM et al. Mol. Cancer Ther. 2008 7:1140.
5. Qin HX et al. Cytotechnology 2013 65:643-53.

Molecular Structure



Reversine

A3 adenosine receptor antagonist; Aurora A/B/C inhibitor

| | |
|--------------------|---|
| Catalog # | A28-900 |
| CAS # | 656820-32-5 |
| Lot # | J3942-24 |
| Purity | >98% (HPLC, TLC); NMR (conforms) |
| Format | Off-white powder |
| Solubility | May be dissolved in DMSO (7 mg/ml) |
| Storage & Shipping | Store desiccated as supplied at -20°C. 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: Reversine

Catalog # A28-900

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)

Not a hazardous substance.

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): Reversine
Chemical Name: 2-(4-Morpholinoanilino)-6-cyclohexylaminopurine
Molecular Formula: C₂₁H₂₇N₇O
CAS Number: 656820-32-5
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.

Article 5 – Fire Fighting Measures

5.1 Extinguishing Media

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

5.2 Special Hazards

Carbon oxides, Nitrogen oxides (NOx).

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.

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 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 container 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: 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 (7 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 oxidizing agents

10.4 Hazardous Decomposition Products: Hazardous decomposition products formed upon combustion: Carbon oxides (CO, CO₂), Nitrogen oxides (NO_x).

Article 11– Toxicological Information

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

Acute Toxicity: data not available.

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

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 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.