

Everolimus

mTOR Inhibitor

Catalog # F17-901D

Lot # O3936-9

CAS # 159351-69-6

Product Description

Molecular Formula: C₅₃H₈₃NO₁₄

Appearance: White powder

Molecular Weight: 956.2

Purity: ≥98% (HPLC); NMR (Conforms)

Solubility: Soluble in DMSO (100 mg/ml); or ethanol (100 mg/ml)

Alias

40-O-(2-hydroxyethyl)rapamycin

Specific Activity

mTOR 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.

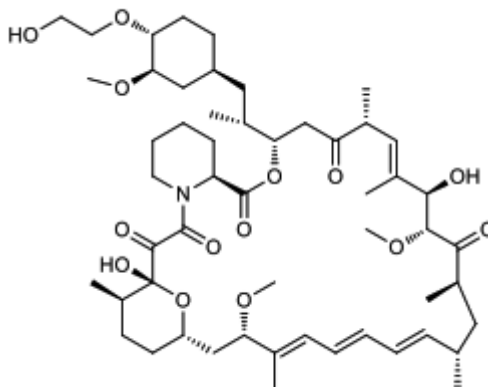
Scientific Background

Semisynthetic analog of rapamycin with potent immunosuppressive activity (1). Inhibitors mTOR (2). Clinically useful immunosuppressant (3).

References

1. HJ Schuurman et al. Transplantation 1997 64:32.
2. P Neuhaus et al. Liver Transpl. 2001 7: 473
3. BD Kahan et al. Clin. Ther. 2002 24:330

Molecular Structure



Everolimus

mTOR Inhibitor

Catalog #	F17-901D
CAS #	159351-69-6
Lot #	O3936-9
Purity	≥98% (HPLC); NMR (Conforms)
Format	White powder
Solubility	Soluble in DMSO (100 mg/ml); or ethanol (100 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 1-866-9KINASE (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: Everolimus

Catalog # F17-901D

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)

Repeated oral exposure: Category 1

2.2 Label Elements

Classification according to Regulation (EC) No 1272/2008



GHS Signal Word: DANGER

GHS Hazard Phrases:

H372	Causes damage to organs through prolonged or repeated exposure if swallowed.
------	--

GHS Precaution Phrases:

P260	Do not breathe dust/fumes/mist/vapours/sprays.
P264	Wash face, hands and any exposed skin thoroughly after handling.
P270	Do not eat, drink, or smoke while using this product.
P273	Avoid release to the environment.
P314	Get medical advice/attention if you feel unwell.
P501	Dispose of any 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: Exercise caution.

Article 3 – Composition/Information on Ingredients

3.1 Substance

Product Name(s): Everolimus
Chemical Name: 40-O-(2-hydroxyethyl)rapamycin
Molecular Formula: C₅₃H₈₃NO₁₄
CAS Number: 159351-69-6
Concentration: 100%

Article 4 – First Aid

4.1 Description of First Aid Measures

General: If symptoms persist, call a physician.

If Inhaled: Remove from exposure. Move to fresh air. If breathing is difficult, administer oxygen. If breathing stops, administer artificial respiration.

In Case of Skin Contact: Wash with soap and water for 15 minutes. Remove contaminated clothing and shoes.

In Case of Eye Contact: Immediately flush with water. After initial flushing remove any contact lenses and continue flushing for at least 15 minutes separating eyelids.

If Swallowed: Rinse mouth. Do NOT induce vomiting unless directed to do so by medical personnel.

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.

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 extinguishing measures that are appropriate to local circumstances.

5.2 Special Hazards

Combustion products are carbon monoxide, carbon dioxide, nitrogen oxide.

5.3 Precautions for Fire Fighters

Use self-contained breathing apparatus and protective clothing to prevent contact with skin and eyes.

Article 6 – Accidental Release Measures

6.1 Personal Precautions

Use personal protective equipment. Avoid contact with skin, eyes, and clothing. Do not eat, drink or smoke when using this product. Remove and wash contaminated clothing before reuse.

6.2 Environmental Precautions

Prevent material from entering drains or water courses.

6.3 Methods for Containment and Clean-up

Prevent further leakage or spillage if safe to do so. Cover powder with wet paper towels. Minimize spreading. Pick up and transfer to properly labeled containers for disposal.

Article 7 – Handling and Storage

7.1 Handling Precautions

Wear personal protective equipment. Avoid contact with skin, eyes, and clothing. Use only in an area approved for laboratory research.

7.2 Safe Storage Measures

This product should be stored at -20°C. It should be kept in a tightly closed container. This product should be handled only by qualified experienced professionals. Wash thoroughly after handling this material.

Article 8 – Exposure Controls/Personal Protection

8.1 Control Parameters

This product does not contain any hazardous materials with occupational exposure limits established by region-specific regulatory bodies.

8.2 Specific Engineering Controls

Showers, eyewash stations, ventilation systems

8.3 Personal Protection

Eye/Face Protection: Tightly fitting safety goggles. Face shield.

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: Wear NIOSH / MSHA-approved respirator if required.

Personal Hygiene Measures: When using this product, do not eat, drink or smoke. Remove and wash contaminated clothing before re-use.

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: soluble in DMSO (100 mg/ml); or ethanol (100 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: None known

10.4 Hazardous Decomposition Products: Carbon monoxide (CO); carbon dioxide (CO₂)

Article 11– Toxicological Information

11.1 Toxicological Effects: The toxicological effects of this product have not been thoroughly studied. Oral Rat LD50: >2000 mg/kg

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

Cyprinus carpio (Carp) LC50 > 18.4 mg/l – 96h

12.2 Degradability

Not readily biodegradable

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 1-866-9KINASE (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.