



Catalog #

Aliquot Size

D430-59-200

0.2 ml

D430-59-500

0.5 ml

DLG (Dilauroyl-sn-glycerol)

Catalog # D430-59

Lot # K4542-1

Product Description

The substrate solution used for assaying Diacylglycerol kinases.

Molecular Weight

The molecular weight of 1,2-Dilauroyl-sn-glycerol (DLG) is: 456.70.

Formulation

Frozen substrate solution containing: 10 mM of 1,2-Dilauroyl-sn-glycerol (DLG) in DMSO.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

DLG (Dilauroyl-sn-glycerol)

Catalog # D430-59

Lot # K4542-1

Concentration 10 mM

Stability 1yr at -20°C from date of shipment

Storage & Shipping Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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 - Product Identification

Product Name: DLG (Dilauroyl-sn-glycerol)**Catalog # D430-59**

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 - Hazard Identification

- **WHMIS Classification:** Not WHMIS controlled.
- **GHS classification:** Flammable Liquids (Category 4)
- **Hazard Pictograms:**



- **Signal words:** Warning
- **Hazard statements:**
 - Combustible liquid
- **Precautionary statements:**
 - Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking.
 - Wear protective gloves, protective clothing and eye protection.
 - In case of fire: Use dry chemical, CO2, water spray or regular foam to extinguish.
- **Other hazards:** None known.

Article 3 – Composition/Information on Ingredients

Chemical Characterization: Mixtures.

Description: This product consists of the substances listed below.

Common name	Chemical name	CAS-No.	Concentration
DMSO (Dimethyl sulfoxide)	1, 1'-sulfinylbis-methane	67-68-5	99.5433%
Diacylglycerol	1,2-dilauroyl-sn-glycerol	60562-15-4	0.4567%

Article 4 – First-aid Measures

- **General information:** Consult a physician by providing the SDS. Move out of dangerous area.
- **After inhalation:** Breathe in fresh air. If cannot breathe, give artificial respiration. Consult a physician if you feel unwell.
- **After skin contact:** Immediately wash with soap and plenty of water and rinse thoroughly. Remove contaminated clothing. Consult a physician.
- **After eye contact:** Rinse opened eyes thoroughly with plenty of water for at least 15 minutes. Consult a physician.
- **After swallowing:** Never give anything by mouth to an unconscious person. Rinse the mouth with plenty of water and consult a physician. Do NOT induce vomiting unless directed to do so by medical personnel.

Article 5 - Fire-fighting Measures

- **Suitable extinguishing media:** Use dry chemical, carbon dioxide, water spray or regular foam to extinguish, or other appropriate measure that is suitable to the environment.
- **Specific hazards arising from the substance or mixture:** None known.
- **Special protective equipment and precautions for fire-fighters:** Self-contained breathing apparatus if necessary.

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 6 – Accidental Release Measures

- **Personal precautions, protective equipment and emergency procedures:** Apply standard laboratory practices and personal protective equipment. Avoid breathing vapors, mist, or gas. Ensure adequate ventilation.
- **Environmental precautions:** Do not allow to enter drains.
- **Methods and materials for containment and cleaning up:** Absorb on sand or vermiculite and place in closed containers for disposal.

Article 7 - Handling and Storage

- **Precautions for safe handling:** Do not swallow. Avoid inhalation, contact with eyes, skin or clothing. Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking.
- **Conditions for safe storage:** Store according to package insert; keep container upright and tightly closed.

Article 8 - Exposure Controls/Personal Protection

- **Components with limit monitoring values at workplace:**
Dimethyl sulfoxide (DMSO); No TLV established. No PEL established.
- **Appropriate engineering controls:**
Follow standard industrial hygiene and laboratory safety practices. Wash hands before breaks and at the end of the workday.
- **Individual protection measures:**
Respiratory protection:
Not required under normal use conditions. Use appropriate respirator if there is inadequate ventilation by following the government standards.
Hand protection:
Wear gloves and use proper glove removal technique to avoid skin contact. Discard gloves after use by following the applicable laboratory regulations. Wash and dry hands.
Eye/face protection:
Safety goggles with side-shields approved under appropriate government standards.
Skin/body protection:
Use appropriate clothing, footwear and any additional protection measures to protect from splashing or contamination.

Article 9 – Physical and Chemical Properties

Appearance: Colorless fluid.	Danger of explosion: No data available.
Odour/Odour Threshold: Not determined.	Explosion limits: Not available.
pH: Not available.	Decomposition temperature: Not available.
Melting point/freezing point: Not determined.	Vapor pressure at 20 °C: Not available.
Boiling point/Boiling range: >100 °C.	Density: Not determined.
Flash point: 87°C (DMSO).	Relative density: Not determined.
Flammability (solid, gaseous): Not determined.	Vapor density: Not determined.
Ignition temperature: Not determined.	Evaporation rate: Not determined.
Auto-igniting: 215°C (DMSO).	Solubility in / Miscibility with Water: Fully miscible.

Article 10 - Stability and Reactivity

- **Reactivity:** Stable under recommended transport and storage conditions.
- **Chemical stability:** Stable under recommended transport and storage conditions.
- **Possible hazardous reactions:** No data available.
- **Conditions to avoid:** Contact with incompatible materials. Sources of ignition. Exposure to heat.
- **Incompatible materials:** Strong oxidizing agents. Perchlorates.
- **Hazardous decomposition products:** Not available.

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 11 - Toxicological Information

- **Acute toxicity:** DMSO; Oral LD₅₀ 14500 mg/kg (rat); Inhalation LD₅₀ >5330 mg/m³ (rat).
- **Skin corrosion/irritation:** DMSO; LD₅₀ dermal 40000 mg/kg.
- **Serious eye damage/eye irritation:** Not available.
- **Respiratory or skin sensitization:** Not available.
- **Germ cell mutagenicity:** Not available.
- **Carcinogenicity:** No components are listed in IARC, or NTP, or OSHA, or ACGIH.
- **Reproductive toxicity:** Not available.
- **Teratogenicity:** Not available.
- **Specific target organ toxicity - single exposure/ - repeated exposure (GHS):** Not available.
- **Aspiration hazard:** Not available.
- **Potential health effects:**
 - Inhalation: No data available
 - Ingestion: No data available
 - Skin: No data available
 - Eyes: No data available
- **Signs and Symptoms of Exposure:** No data available
- **Synergistic effects:** Not available.
- **Additional information:** To the best of our knowledge, the chemical, physical, and toxicological properties have not been thoroughly investigated.

Article 12 - Ecological Information

- **Eco-toxicity:** No data available.
- **Biodegradability:** No data available.
- **Bio-accumulative potential:** No data available.
- **Mobility in soil:** No data available.
- **PBT and vPvB assessment:** Not applicable.
- **Other adverse effects:** No data available.

Article 13 - Disposal Considerations

- **Disposal methods:** In accordance to applicable national, regional, or local laws and regulations. For additional handling information and protection of employees please refer to Article 7 and 8.
- **Contaminated packaging:** Disposal should be made in accordance to official regulations. Use water or cleansing agents to clean the area.

Article 14 - Transport Information

- **DOT:** Not dangerous goods.
- **IMDG:** Not dangerous goods.
- **IATA:** Not dangerous goods.

Article 15 - Regulatory Information

- This product has been classified in accordance with the hazard criteria of the Hazardous Products Regulations (HPR) and the SDS contains all the information required by the HPR.

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.