



METTL3/METTL14 Substrate

Synthetic oligonucleotide as METTL3/METTL14 complex substrate

Catalog # M323-58

Lot# U4760-1

Product Description

The synthetic single-stranded 27 base oligoribonucleotide is used as a substrate for the METTL3/METTL14 methyltransferase complex assay.

Application

A methyltransferase METTL3/METTL14 complex assay.

Molecular Weight

The size of single-stranded oligoribonucleotide is 27 bases. The molecular weight is 8553.2.

Format

White to off-white lyophilized powder.

Reconstitution Protocol

2.5 nmol of lyophilized single-stranded oligoribonucleotide is supplied. Rehydrate in 25µl of nuclease-free water or 20mM Tris-HCl, pH 7.5, to a final concentration of 100µM. Prior to use, add 1 unit/µl RNase inhibitor.

Storage and Stability

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

Related METTL3/METTL14 complex

METTL3/METTL14 Substrate can be utilized as a substrate for the active METTL3/METTL14 methyltransferase.

Product Name	Catalog Number
METTL3/METTL14, Active	M323-380G

METTL3/METTL14 Substrate

Synthetic oligonucleotide as METTL3/METTL14 complex substrate

Catalog #	M323-58
Quantity	2.5 nmol
Lot #	U4760-1
Format	Lyophilized powder
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.

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: METTL3/METTL14 Substrate

Catalog # M323-58

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.

Supplier'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 - Hazards Identification

Emergency Overview: The product contains no substances which at their given concentration, are considered to be hazardous to health.
 WHMIS Classification: Not WHMIS controlled
 GHS Classification: Not a dangerous substance according to GHS.

Labeling: Observe the general safety regulations when handling chemicals.
 This substance is not subject to identification regulations.

Safety Phrases: The material and its container must be disposed of in a safe way.

Article 3 - Composition/Information on Ingredients

Common Name	Chemical Name/Formula	CAS-No.	Concentration
ssRNA	-	NA	≤100%

Article 4 - First-aid Measures

- **General information:** No special measures required.
- **After inhalation:** Supply fresh air; consult a physician in case of complaint.
- **After skin contact:** Wash with soap and plenty of water and rinse thoroughly. Generally, the product does not irritate skin.
- **After eye contact:** Rinse opened eyes with plenty of water for at least 15 minutes. Consult a physician.
- **After swallowing:** Rinse the mouth with plenty of water and consult a physician.

Article 5 - Fire-fighting Measures

- **Suitable extinguishing media:** Use water spray, extinguishing powder, carbon dioxide, 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.

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:** No special measures required.
- **Methods and materials for containment and cleaning up:** Dispose contaminated material as waste according to item 13.

Article 7 - Handling and Storage

- **Precautions for safe handling:** Apply standard laboratory practices and personal protective equipment.
- **Conditions for safe storage:** Store according to product specification.

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 8 - Exposure Controls/Personal Protection

- **Appropriate engineering controls:**
Follow standard laboratory practices.
 - **Individual protection measures:**
 - Respiratory protection:**
Ensure adequate ventilation.
 - Skin/body protection:**
Follow standard laboratory practices and use standard personal protective equipment.
 - Eye/face protection:**
Safety glasses approved under appropriate government standards.
 - General hygiene considerations:**
Handle in accordance with good industrial hygiene and safety practice. Wash and dry hands after use.
-

Article 9 – Physical and Chemical Properties

Appearance: Solid.	Danger of explosion: No data available.
Odour/Odour Threshold: Not determined.	Explosion limits: No data available.
pH: No data available.	Decomposition temperature: No data available.
Melting point/freezing point: Not determined.	Vapor pressure at 20 °C: No data available.
Boiling point/Boiling range: No data available .	Density: No data available.
Flash point: No data available .	Relative density: No data available.
Flammability (solid, gaseous): Not determined.	Vapor density: No data available.
Ignition temperature: No data available.	Evaporation rate: No data available.
Auto-igniting: No data available .	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 dangerous reactions known.
 - **Conditions to avoid:** No data available.
 - **Incompatible materials:** No data available.
 - **Hazardous decomposition products:** No data available.
-

Article 11 - Toxicological Information

- **Acute toxicity:** Product does not present an acute toxicity hazard based on known or supplied information.
- **LD/LC50:** Not available.
- **Skin corrosion/irritation:** Not available.
- **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 irritant effect.
 - Ingestion:** No known harmful effects.
 - Skin:** No irritant effect.
 - Eyes:** No irritant effect.
- **Signs and Symptoms of Exposure:** Not available.
- **Synergistic effects:** 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 12 - Ecological Information

- **Eco-toxicity:** Not applicable.
- **Biodegradability:** Not applicable.
- **Bio-accumulative potential:** Not applicable.
- **Mobility in soil:** Not applicable.
- **PBT and vPvB assessment:** Not applicable.
- **Other adverse effects:** Not applicable.

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

Safety, health, and environmental regulations/legislation specific for the substance or mixture.

Canadian substance listings:

- **WHMIS Classification:** Non-hazardous.
- **GHS label elements:** Not applicable.
- **Signal word:** Not applicable.
- **Hazard statements:** Not applicable.

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.