

## LRRKtide

Synthetic peptide substrate derived from ezrin/moesin/radixin

**Catalog # L10-58**

Lot # E3174-7

### Product Description

The LRRKtide peptide sequence (RLGRDKYKTLRQIRQ) is derived from human ezrin (amino acids 561-573), moesin (amino acids 539-553) and radixin (amino acids 558-570) and is suitable for use as a substrate for LRRK kinases.

### Molecular Weight

The theoretical molecular weight is 1931.29

### Purity

The purity was determined to be 97.36% by HPLC analysis.

### Formulation

1mg of peptide supplied as a lyophilized powder.

### 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.

### Reconstitution Protocol

Dilute peptide in 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

## LRRKtide

Synthetic peptide substrate derived from ezrin/moesin/radixin

Catalog Number L10-58

Quantity 1 mg

Specific Lot Number E3174-7

Purity 97.36%

Format 1 mg 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-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**

Use of this product in assays for LRRK2 inhibitors falls within the scope of the claims of the Medical Research Council's granted US patent 7,947,468 and international patent application WO2008122789 and consequently any commercial use of this product in such assays may require a licence from the Medical Research Council. For enquiries about commercial licences to this patent family, please contact Dr. Mary Canning at MRC Technology. Email: [mary.canning@tech.mrc.ac.uk](mailto:mary.canning@tech.mrc.ac.uk) Telephone: +44(0)20 7391 2801

# MATERIAL SAFETY DATA SHEET

## Article 1 - Product Identification and Use

**Product Name: LRRKtide**

**Catalog # L10-58**

*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.*

Manufacturer's Name: SignalChem Pharmaceuticals 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 Ingredients

NOT AVAILABLE. We are not aware of any hazards associated with this product or its ingredients, but the chemical, physical, and toxicological properties of this product have not been investigated thoroughly. Observe normal laboratory precautions.

## Article 3 - Physical Data

The physical properties of this product have not been investigated thoroughly.

## Article 4 - Fire and Explosion Hazard

NOT APPLICABLE

## Article 5 - Reactivity Data

NOT APPLICABLE

## Article 6 - Toxicologically Data

May be harmful by inhalation, ingestion, or skin absorption. The toxicological properties of this product have not been investigated thoroughly. Exercise due caution.

## Article 7 - Preventative Measures

Wear chemical safety goggles and compatible chemical-resistant gloves. Avoid inhalation, contact with eyes, skin or clothing.

\*\*\*\*\*MULTIPLE COMPONENT SPILL OR LEAK PROCEDURES\*\*\*\*\*

- Wear protective equipment.
- Absorb on sand or vermiculite and place in closed containers for disposal.
- Observe all federal, state and local environmental regulations.

## Article 8 - First Aid Measures

- If swallowed, wash out mouth with water, provided person is conscious. Call a physician.
- In case of skin contact, flush with copious amounts of water for at least 15 minutes. Remove contaminated clothing and shoes. If a rash or other irritation develops, call a physician.
- If inhaled, remove to fresh air. If breathing becomes difficult, call a physician.
- In case of eye contact, flush with copious amounts of water for at least 15 minutes while separating the eyelids with fingers. Call a physician.

## Article 9 - Preparation

Prepared By: Jun Yan

Phone #: 1-866-954-6273

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 Catalogue 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**

Use of this product in assays for LRRK2 inhibitors falls within the scope of the claims of the Medical Research Council's granted US patent 7,947,468 and international patent application WO2008122789 and consequently any commercial use of this product in such assays may require a licence from the Medical Research Council. For enquiries about commercial licences to this patent family, please contact Dr. Mary Canning at MRC Technology. Email: [mary.canning@tech.mrc.ac.uk](mailto:mary.canning@tech.mrc.ac.uk) Telephone: +44(0)20 7391 2801