

## PDKtide

Synthetic peptide substrate derived from Akt1 and PKN2

### Catalog # P10-58

Lot # E357-5

### Product Description

The 39 amino acids of PDKtide peptide sequence (KTFCGTPEYLAPEVRREPRILSEEEQEMFRDFDYIADWC) is derived from two human proteins: residues 1-14 are based on AKT1 (307-320) and residues 16-39 are based on PKN2/PRK2 (961-984).

### Molecular Weight

The theoretical molecular weight is 4771.36

### Purity

The purity was determined to be more than 96.37% 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 40mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

### Related Kinases

PDKtide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
PDK1, Active	P14-10H
CDC7/DBF4, Active	C26-10G
CDK9/CyclinK, Active	C40-10G

## PDKtide

Synthetic peptide substrate derived from AKT1 and PKN2

Catalog Number P10-58

Quantity 1 mg

Specific Lot Number E357-5

Purity >96.37%

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.

# MATERIAL SAFETY DATA SHEET

---

## Article 1 - Product Identification and Use

---

**Product Name: PDKtide**

**Catalog # P10-58**

*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: 550-5600 Parkwood Way  
City, Prov. Postal Code: Richmond, BC, V6V 2M2  
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

---

This product consists of purified protein in Tris-HCl buffer shipped on dry ice. 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.**