

## CK1tide

Synthetic peptide evaluated as a substrate for CK1 family

### Catalog # C07-58

Lot # K1768-2

### Product Description

The CK1tide synthetic peptide [HAAIGDDDDAYSITA-NH<sub>2</sub>] is routinely evaluated as a substrate for CK1 family kinases, such as CK1 and CK1 delta.

### Molecular Weight

The theoretical molecular weight is 1533.58.

### Purity

The purity was determined to be 99.72% 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.

## CK1tide

Synthetic peptide evaluated as a substrate for CK1 family

Catalog #	C07-58
Quantity	1 mg
Lot #	K1768-2
Purity	99.72%
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.