

## CK2 Substrate

Synthetic peptide substrate evaluated for CK2 family assay

**Catalog # C08-58**

Lot # I278-3

### Product Description

The synthetic peptide (RRRADDSDDDDD) contains serine protein kinase phosphorylation site surrounded by acidic residues and is high selectively evaluated as a substrate for CK2 family kinases.

### Molecular Weight

The theoretical molecular weight is 1450.37

### Purity

The purity was determined to be 97.24% 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.

### Related Kinases

CK2 Substrate peptide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
CK2α1, Active	C70-10G
CK2α2, Active	C71-10G

## CK2 Substrate

Synthetic peptide substrate evaluated for CK2 family assay

Catalog Number	C08-58
Quantity	1 mg
Specific Lot Number	I278-3
Purity	97.24 %
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.**