

AKT (PKB) Substrate

Synthetic peptide substrate derived from GSK3

Catalog # A05-58

Lot # H103-7

Product Description

The AKT (PKB) Substrate peptide sequence (CKRPRAASFAE) is based on the N-terminus of GSK3.

Molecular Weight

The theoretical molecular weight is 1235.43

Purity

The purity was determined to be 95.71% by HPLC analysis.

Formulation

1 mg 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

AKT (PKB) Substrate can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
AKT1, Active	A16-10G
AKT2, Active	A17-10G
AKT3, Active	A18-10G
SGK1, Active	S06-11G
SGK2, Active	S07-10G
SGK3, Active	S08-10G
PAK4, Active	P05-10G
PAK7, Active	P07-10G

AKT (PKB) Substrate

Synthetic peptide substrate derived from GSK3

Catalog Number	A05-58
Quantity	1 mg
Specific Lot Number	H103-7
Purity	95.71%
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
www.signalchem.com

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