

SAMStide

Synthetic peptide substrate derived from the mouse ACACA

Catalog # S07-58

Lot # D303-3

Product Description

The SAMStide peptide sequence (HMRSAMSGHLVKRR) is based on the mouse acetyl-Coenzyme A carboxylase alpha (amino acid 73-85).

Molecular Weight

The theoretical molecular weight is 1779.18

Purity

The purity was determined to be 95% 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

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

Product Name	Catalog Number
AMPK (A1/B1G1), Active	P47-10H
AMPK (A2/B1G1), Active	P48-10H
AMPK (A1/B1G2), Active	P55-10H
AMPK (A1/B1G3), Active	P56-10H
AMPK (A1/B2G1), Active	P50-10H

SAMStide

Synthetic peptide substrate derived from the mouse ACACA

Catalog Number S07-58

Quantity 1 mg

Specific Lot Number D303-3

Purity 95%

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.