

## SAMStide

Synthetic peptide substrate derived from the mouse ACACA

**Catalog # S07-58**

Lot # R339-3

### Product Description

The SAMStide peptide sequence (HMRSAMSGHLVKKRR) 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 95.9% by HPLC analysis.

### Formulation

1mg of peptide supplied as a lyophilized powder.

### Storage and Stability

Store product at  $-20^{\circ}\text{C}$ . 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 R339-3

Purity 95.9%

Format 1 mg lyophilized powder

Stability 1yr at  $-20^{\circ}\text{C}$  from date of shipment

Storage & Shipping Store product at  $-20^{\circ}\text{C}$ . 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.