



Product Data Sheet

Product Name: SAMS Peptide
Catalog Number: AS-60167-5 (5 mg) **Lot Number:** See label on vial

Sequence: H-His-Met-Arg-Ser-Ala-Met-Ser-Gly-Leu-His-Leu-Val-Lys-Arg-Arg-NH₂
(3-letter code)
HMRSAMSGHLHLVKRR-NH₂ (1-letter code)

Molecular Weight: 1778.2

% Peak Area by HPLC: ≥95

Appearance: Lyophilized white powder

Peptide Reconstitution: Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.

Storage: Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.

Description: A synthetic peptide substrate specific for AMP-activated protein kinase (AMPK).
Ref: Davies, SP. et al. *Eur. J. Biochem.* **186**, 123 (1989).

Additional Information: *Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.*

AMPK was immunoprecipitated from 5 µg of cell lysate by incubation with 1 µg of anti- 1 and 1 µg of anti-α2 polyclonal antibodies, and AMPK activity in the IP complexes was determined by phosphorylation of peptide HMRSAMSGHLHLVKRR (SAMS). Synthetic peptides encompassing Ser77 (71QGTVNWSVDDIVKGI85), Ser105 (98QAARKLLSREKQ PPID113), and Ser179 (172ASPHAHISEQAVWALG187) were obtained from AnaSpec, Inc. (San Jose, CA) and used in phosphorylation assays employing AMPK as kinase, following the procedures described above for the SAMS peptide. Phosphorylated substrates were measured by scintillation counting-[Wang, W. et al. *J Biol Chem* **279**, 48376 \(2004\).](#)

Published Citations:

[Wang, W. et al. *J Biol Chem* **279**, 48376 \(2004\).](#)

For Research Use Only