

# SAMS peptide (phosphorylated) (biotinylated)

## AMPK substrate

This peptide, biotin-HMRSAMpSGLHLVKRR, contains a biotin group at its amino terminus and a phospho-serine. It is derived from rat acetyl-CoA carboxylase. It may be used as a control peptide for unphosphorylated SAMS Peptide (Prod. No. BML-P231 and BML-P232). The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P233-0001	1mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**MW** 2199

**Purity** ≥95%

**Sequence** biotin-His-Met-Arg-Ser-Ala-Met-pSer-Gly-Leu-His-Leu-Val-Lys-Arg-Arg-OH



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)