

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 »](#)

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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