

Product Data Sheet

Product Name: [Ser25]-PKC (19-31), biotinylated

Catalog Number: AS-23018 (1 mg) Lot Number: See label on vial

Sequence: H-Lys(Biotin)-Arg-Phe-Ala-Arg-Lys-Gly-Ser-Leu-Arg-Gln-Lys-Asn-Val-OH

(3-letter code)

K(Biotin)-RFARKGSLRQKNV (1-letter code)

Molecular Weight: 1914.3 % 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.

For Research Use Only