



Product Data Sheet

Product Name: Myristolated PKC Zeta, Pseudosubstrate (ZIP)
Catalog Number: AS-63361 (1 mg) Lot Number: See label on vial
Sequence: Myr-Ser-Ile-Tyr-Arg-Arg-Gly-Ala-Arg-Arg-Trp-Arg-Lys-Leu-OH (3-letter code)
Myr-SIYRRGARRWRKL (1-letter code)

Molecular Weight: 1928.5

% 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.

Published Citations:

[Migues, P. et al. *Nat Neurosci* 13, 630 \(2010\).](#)

For Research Use Only