

PRODUCT DATA SHEET



PEPTIDE NAME: CaM kinase II (290-309), calmodulin antagonist
SEQUENCE: Leu-Lys-Lys-Phe-Asn-Ala-Arg-Arg-Lys-Leu-Lys-Gly-Ala-Ile-Leu-Thr-Thr-Met-Leu-Ala

SEQUENCE DESCRIPTION: Calcium/calmodulin-dependent protein kinase II pseudosubstrate site.
[115044-69-4]

CATALOG NO.: P-200

LOT NO.: K1918

MOLECULAR WEIGHT: 2274.2

PURITY: >97% by HPLC analysis.

AMINO ACID ANALYSIS AND IDENTITY: Confirms.

PEPTIDE CONTENT: 80.3%

SUPPLIED AS: Trifluoroacetate salt; 0.5 mg net peptide/vial.

PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: H₂O

STORAGE: -20°C

OTHER: See attached Guide to Handling and Storing Peptides.

LITERATURE REFERENCES:

M.E. Payne *et al.* *J. Biol. Chem.* 1988 **263** 7190