

## **Product Data Sheet**

Product Name: UOM9, PKC Substrate, phosphorylated

Catalog Number: AS-20294 (1 mg) Lot Number: See label on vial

Sequence: H-Lys-Arg-Pro-pSer-Gln-Arg-His-Gly-Ser-Lys-Tyr-NH2 (3-letter code)

KRP-pS-QRHGSKY-NH2 (1-letter code)

Molecular Weight: 1422.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: UOM9, PKC peptide is freely soluble in water.

Storage: UOM9, PKC peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Reference: Ref: Rusnak, F. et al. J. Biomol. Tech. 13, 228 (2002).

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Anti-PKC- $\alpha$  and anti-PKC- $\beta$ I antibodies were preincubated with a ten-fold weight excess of the corresponding peptide provided by the manufacturer of the antibody (Santa Cruz Biotechnology) at +4C for 18 hr. Anti-phospho-PKC substrate antibodies were preincubated similarly with a phosphorylated peptide that corresponds to the phosphorylation pattern and amino acid sequence that this antibody detects (phosphorylated PKC substrate 3; H-Lys-Arg-ProphosphoSer-Gln-Arg-His-Gly-Ser-Lys-Tyr-NH2) (AnaSpec, Inc.; San Jose, CA)- Aaltonen, V. et al. J. Histochem. Cytochem. 54, 795 (2006).

## **Published Citations:**

Aaltonen, V. et al. J. Histochem. Cytochem. 54, 795 (2006).

## Related Products:

Name Cat # Size

Phosphopeptide Mass Spec Standards AS-61145 1 kit

For Research Use Only