

Product Name:	RS domain derived peptide	
Catalog Number:	AS-61722 (1 mg)	Lot Number: See label on vial
Sequence:	H-Gly-Arg-Ser-Arg-Ser-Arg-Ser-Arg-Ser-Arg-OH (3-letter code) GRSRSRSRSR (1-letter code)	
Molecular Weight:	1205.3	
% Peak Area by HPLC: ≥ 95		
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Using H_2O , reconstitute by adding 100 μ I to 1 mg RS domain derived peptide.

Storage: RS domain derived 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.

Description: This peptide is a substrate for Clk/Sty and is phosphorylated by Clk/Sty protein kinase (Km=102 microM). Zahler, AM. et al. *Genes Dev.* **6**, 837 (1992).

For Research Use Only