Catalog #: AS-60298 (5mg)

AnaSpec, Inc.

Certificate of Analysis

Peptide Name: PKG Substrate 2

Run Number: 32707

> H-Arg-Lys-Arg-Ser-Arg-Ala-Glu-OH Sequence:

903.03 Theoretical Mass (M+H⁺):

> 902.75 Mass Found (M+H⁺):

> > Dissolve 1mg of peptide in 1ml Water Solubility:

Appearance: White crystal

HPLC Purity: > 95 %

Amount Packed: net mg

Amino Acid Analysis

	Peptide Content:		76.1	%				
	CALC	FOUND	C	CALC	FOUND		CALC	FOUND
Asx			Ala	1	0.89	lle		
Ser	1	0.84	Pro			Leu		
Glx	1	1.05	Cys			Phe		
Gly			Tyr			Trp		
His			Val			NorLeu		
Arg	3	3.00	Met			AHX		
Thr			Lys	1	1.22	Other		
Note:						Other		

Remarks: Not for Human Use. Research Purposes Only.

Released by: Date: _____ Quality Control