

Certificate of Analysis

Peptide Name: Biotin-LC-Angiotensin II
Run Number: 31997
Sequence: Biotin-LC-Asp-Arg-Val-Tyr-Ile-His-Pro-Phe-OH

Theoretical Mass (M+H⁺): 1386.7

Mass Found (M+H⁺): 1386.8

Solubility: Dissolve 1mg of peptide in 1ml water

Appearance: White powder

HPLC Purity: > 95 %

Amount Packed: 5 mg net peptide

Amino Acid Analysis

Peptide Content: 70.4 %

	CALC	FOUND		CALC	FOUND		CALC	FOUND
Asx	1	1.11	Ala			Ile	1	0.97
Ser			Pro	1	1.02	Leu		
Glx			Cys			Phe	1	0.99
Gly			Tyr	1	0.73	Trp		
His	1	0.93	Val	1	1.12	NorLeu		
Arg	1	1.13	Met			AHX		
Thr			Lys			Other		
Note:						Other		

Remarks: Not for Human Use. Research Purposes Only.

Released by: _____ **Date:** _____

Quality Control