Certificate of Analysis

	ide Name:	Amiı	Aminopeptidase N Ligand(CD13)							
Sequence:			Н-Су	H-Cys-Asn-Gly-Arg-Cys-Gly-OH						
Theoretical Mass (M+H⁺):			609.7	609.7						
Ма	d (M+H⁺):	609.	609.5							
Solubility:			Diss	Dissolve 1mg of peptide in 1ml Water						
Appearance:			White Powder							
	HPI	_C Purity:	> 95	; o	6					
		-		-						
Amount Packed: 5 mg net peptide										
Amin	o Acid A	nalysis								
	Peptide	e Content:	71.0	%						
	CALC	FOUND	C	ALC	FOUND		CALC	FOUND		
Asx	1	1.06	Ala			lle				
Ser			Pro			Leu				
Glx			Cys	2	1.12	Phe				
Gly	2	1.73	Tyr			Trp				
His			Val			NorLeu				
Arg	1	1.11	Met			AHX				
Thr			Lys			Other				
Note:			-			Other				
		Rem	arks: Not	for Hu	man Use. Res	earch Purpo	ses Only	<i>.</i> .		
Re	leased b	oy:	Quality Cor		Da	ate:				