

L- α -aspartyl-L-methionyl-L-seryl-L-seryl-L- α -aspartyl-L-leucyl-L- α -glutamyl-L-arginyl-L- α -aspartyl-L-histidyl-L-arginyl-L-prolyl-L-histidyl-L-valyl-L-seryl-L-methionyl-L-prolyl-L-glutamyl-L-asparaginyl-L-alanyl-L-asparagine, trifluoroacetate salt

Purity Specification:	≥95%			
Molecular formula:	C97H154N34O36S2 • XCF3COOH			
CAS number:	Formula weight:	2436.60	Expiration date:	20SEP2029

Tests	Results
HPLC	Purity: 99.1 %
Mass Spec	[M+3H] ³⁺ : 813.2 [M+2H] ²⁺ : 1219.1

Shin

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM