

L-histidylglycyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-leucyl-L-seryl-L-lysyl-L-glutamyl-L-methionyl-L- α -glutamyl-L- α -glutamyl-L- α -glutamyl-L-alanyl-L-valyl-L-arginyl-L-leucyl-L-phenylalanyl-L-isoleucyl-L- α -glutamyl-L-tryptophyl-L-leucyl-L-lysyl-L-asparaginylglycylglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-seryl-L-lysyl-L-lysyl-L-lysyl-L-lysyl-L-lysyl-L-lysyl-L-lysineamide, acetate

Expiry date: 13JUN2028

Tests	Results
HPLC	Purity: 98.8 %
Mass spec	[M+4H]4+: 1215.3 [M+5H]5+: 972.4 [M+6H]6+: 810.5 [M+7H]7+: 695.0
Counterion	Conforms



CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM