

L-histidyl-2-methylalanyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-valyl-L-seryl-L-seryl-L-tyrosyl-L-leucyl-L- α -glutamylglycyl-L-glutaminy-L-alanyl-d4-L-alanyl-d4-N6-[N-(17-carboxy-1-oxoheptadecyl)-L- γ -glutamyl-2-[2-(2-aminoethoxy)ethoxy]acetyl-2-[2-(2-aminoethoxy)ethoxy]acetyl]-L-lysyl-L- α -glutamyl-L-phenylalanyl-L-isoleucyl-L-alanyl-L-tryptophyl-L-leucyl-L-valyl-L-arginylglycyl-L-arginyl-glycine

Purity Specification:	≥90%			
Molecular formula:	C187H283D8N45O59			
CAS number:	Formula weight:	4121.69	Expiration date:	27JUL2030

Tests	Results
Mass Spec	[M+4H]4+: 1031.0
HPLC	Purity: 91.5 %
NMR	Conforms

Reviewed and approved by: Michelle Riley

MSK

WWW.CAYMANCHEM.COM