

Item No. 24992 * Batch No. 0527851

CAS number: Formula weight: 1528.69 Expiration date: 22JUN2030

Tests	Results
Mass Spec	[M+2H]2+: 765.3
HPLC	Purity: 95.4 %
Peptide content	0.00
Peptide Co	77.00

Reviewed and approved by: Timothy Schrot

Tim Sauer

Copyright Cayman Chemical Company, 10/12/2018

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM