

Item No. 24984 * Batch No. 0527843

CAS number: Formula weight: 1455.68 Expiration date: 25JAN2029

Tests	Results
Mass Spec	[M+3H]3+: 486.0
HPLC	Purity: 98.1 %
Peptide content	0.00
Peptide Co	77.00

Reviewed and approved by: Timothy Schrot

Tim Sauer

Copyright Cayman Chemical Company, 10/12/2018

WWW.CAYMANCHEM.COM