

Item No. 27768 • Batch No. 0557143

Expiry date: 22APR2025

Tests	Results
Mass spec	[M+5H] ⁵⁺ : 521.2
HPLC	Purity: 95.3 %
Peptide content	62.8

more

WWW.CAYMANCHEM.COM