

Item No. 27425 * Batch No. 0820044

CAS number: Formula weight: 1827.26 Expiration date: 29MAY2030

Tests	Results
Mass Spec	[M+2H+Na]2+: 936.1
HPLC	Purity: 95.2 %

Reviewed and approved by: Michelle Riley

MORLEY

Copyright Cayman Chemical Company, 10/12/2018

WWW.CAYMANCHEM.COM