

Item No. 36922 * Batch No. 0678457

CAS number: Formula weight: 1796.00 Expiration date: 12NOV2029

Tests	Results
HPLC	Purity: 100.0 %
Mass Spec	[M-2H]2-: 896.6 [M-3H]3-: 597.5

Reviewed and approved by: Libby Perry

Gibby Perry

WWW.CAYMANCHEM.COM