

Item No. 42401 * Batch No. 0724949

CAS number: Formula weight: 1737.16 Expiration date: 16JAN2029

Tests	Results
HPLC	Purity: 98.9 %
Mass Spec	[M+3H]3+: 579.6 [M+4H]4+: 435.4
19F-NMR	Conforms

Gibby Perry

WWW.CAYMANCHEM.COM