

Item No. 38850 * Batch No. 0693115

CAS number: Formula weight: 1461.70 Expiration date: 06FEB2030

Tests

Results

Purity: 99.7 %

[M+2H]²⁺: 731.5

Conforms

Reviewed and approved by: Erik Guetschow

E. Johnson

WWW.CAYMANCHEM.COM