

Item No. 43858 * Batch No. 0817701

Expiration date: 02FEB2029

Results

Conforms

Reviewed and approved by: Sara Rozny

Jan 5. 1888

WWW.CAYMANCHEM.COM