

Lot specific information

Product: 6013327 ULTIMA GOLD / 4x2.5 L

Mfg.Date: 04-DEC-2023

Batch#: 77-23495

Exp.Date: 01-JUL-2025

Information

The product has been analyzed and is certified to have met the specifications.

Test	Results
Quench parameter (tSIE)	671
Background 0-18.6 keV	17 cpm

**This Certificate of Analysis has been produced electronically and is valid without a signature.
For questions regarding this Certificate of Analysis, please contact your customer service representative.**