

Lot specific information

Product: 6013329 Ultima Gold

Mfg.Date: 03-MAR-2025

Batch#: 77-25101

Exp.Date: 01-OCT-2026

Information

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

Test	Results
Quench parameter (tSIE)	672
Background 0-18.6 keV	14 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.