

## Lot specific information

Product: 6013329 Ultima Gold

Mfg.Date: 13-JAN-2025

Batch#: 77-25031

Exp.Date: 01-AUG-2026

## Information

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

Test	Results
Quench parameter (tSIE)	668
Background 0-18.6 keV	15 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.