

Lot specific information

Product: 6013329 Ultima Gold

Mfg.Date: 18-MAR-2024

Batch#: 77-24125

Exp.Date: 01-OCT-2025

Information

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

Test	Results
Quench parameter (tSIE)	665
Background 0-18.6 keV	13 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.**