

Lot specific information

Product: 6013321 Ultima Gold

Mfg.Date: 22-JAN-2024

Batch#: 77-24045

Exp.Date: 01-AUG-2025

Information

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

| Test | Results |
|-------------------------|---------|
| Quench parameter (tSIE) | 673 |
| 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.**