

Lot specific information

Product: 6013377 Ultima Gold LLT

Mfg.Date: 16-OCT-2023

Batch#: 97-23421

Exp.Date: 01-MAY-2025

Information

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

| Test | Results |
|-----------------------------------|----------|
| Quench parameter (tSIE) | 667 |
| Quench parameter (tSIE) 50% water | 281 |
| Low Level background 50% water | 1.83 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.**