

Lot specific information

Product: 6013179 ULTIMA GOLD F / 2x5 L

Mfg.Date: 06-NOV-2023

Batch#: 78-23451

Exp.Date: 01-DEC-2024

Information

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

| Test | Results |
|--|---------|
| Quench parameter (tSIE) | 1001 |
| Background 0-18.6 keV | 14 cpm |
| Quench parameter (tSIE) with Soluene-350 | 794 |
| Background with Soluene-350 | 355 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.**

