

**Lot specific information**

**Product:** 6013329 Ultima Gold

**Mfg.Date:** 04-DEC-2023

**Batch#:** 77-23495

**Exp.Date:** 01-JUL-2025

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	671
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.**