

**Lot specific information**

**Product:** 6013329 Ultima Gold

**Mfg.Date:** 02-JUL-2024

**Batch#:** 77-24275

**Exp.Date:** 01-JAN-2026

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	663
Background 0-18.6 keV	15 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.**