

**Lot specific information**

**Product:** 6013377 Ultima Gold LLT

**Mfg.Date:** 21-OCT-2024

**Batch#:** 97-24435

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

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	661
Quench parameter (tSIE) 50% water	278
Low Level background 50% water	2.25 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.**