

**Lot specific information**

**Product:** 6013371 Ultima Gold LLT

**Mfg.Date:** 13-NOV-2023

**Batch#:** 97-23461

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

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	663
Quench parameter (tSIE) 50% water	279
Low Level background 50% water	1.91 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.**

