

Lot specific information

Product: 6013371 Ultima Gold LLT

Mfg.Date: 02-SEP-2024

Batch#: 97-24365

Exp.Date: 01-APR-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	277
Low Level background 50% water	2.71 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.**

