

Lot specific information

Product: 6013377 Ultima Gold LLT

Mfg.Date: 19-FEB-2024

Batch#: 97-24085

Exp.Date: 01-SEP-2025

Information

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

Test	Results
Quench parameter (tSIE)	668
Quench parameter (tSIE) 50% water	278
Low Level background 50% water	2.02 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.**