

Lot specific information

Product: 6013179 Ultima Gold F

Mfg.Date: 19-AUG-2024

Batch#: 78-24345

Exp.Date: 01-SEP-2025

Information

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

Test	Results
Quench parameter (tSIE)	970
Background 0-18.6 keV	15 cpm
Quench parameter (tSIE) with Soluene-350	756
Background with Soluene-350	2104 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.**