

Lot specific information

Product: 6013179 Ultima Gold F

Mfg.Date: 25-MAR-2024

Batch#: 78-24135

Exp.Date: 01-MAY-2025

Information

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

Test	Results
Quench parameter (tSIE)	1007
Background 0-18.6 keV	14 cpm
Quench parameter (tSIE) with Soluene-350	811
Background with Soluene-350	201 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.**