

Lot specific information

Product: 6013171 Ultima Gold F

Mfg.Date: 08-JAN-2024

Batch#: 78-24025

Exp.Date: 01-FEB-2025

Information

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

Test	Results
Quench parameter (tSIE)	991
Background 0-18.6 keV	13 cpm
Quench parameter (tSIE) with Soluene-350	783
Background with Soluene-350	274 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.**