

**Lot specific information**

**Product:** 6013179 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.**