

**Lot specific information**

**Product:** 6013179 Ultima Gold F

**Mfg.Date:** 16-OCT-2023

**Batch#:** 78-23425

**Exp.Date:** 01-NOV-2024

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	959
Background 0-18.6 keV	16 cpm
Quench parameter (tSIE) with Soluene-350	763
Background with Soluene-350	620 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.**