

**Lot specific information**

**Product:** 6013171 Ultima Gold F

**Mfg.Date:** 04-NOV-2024

**Batch#:** 78-24455

**Exp.Date:** 01-DEC-2025

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	962
Background 0-18.6 keV	13 cpm
Quench parameter (tSIE) with Soluene-350	731
Background with Soluene-350	5087 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.**

