

**Lot specific information**

**Product:** 6013119 Ultima Gold XR

**Mfg.Date:** 02-SEP-2024

**Batch#:** 79-24365

**Exp.Date:** 01-APR-2026

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	530
Background 0-18.6 keV	15 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.**