

Lot specific information

Product: 6013171 Ultima Gold F

Mfg.Date: 04-DEC-2023

Batch#: 78-23495

Exp.Date:

Information

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

Test	Results
Quench parameter (tSIE)	977
Background 0-18.6 keV	11 cpm
Quench parameter (tSIE) with Soluene-350	773
Background with Soluene-350	458 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.**