



Certificate of Analysis

Revvity Health Sciences B.V.
Rigaweg 22
9723 TH Groningen
The Netherlands

Phone: + 31-50-544 59 00
SDS.Groningen@revvity.com
www.revvity.com

Product: 6013327 Ultima Gold

Mfg.Date: 14-AUG-2023

Batch#: 77-23331

Exp.Date: 01-MAR-2025

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

Test	Results
Quench parameter (tSIE)	661
Background 0-18.6 keV	14 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.