



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: 6013326 Ultima Gold

Mfg.Date: 04-SEP-2023

Batch#: 77-23365

Exp.Date: 01-APR-2025

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

| Test | Results |
|-------------------------|---------|
| Quench parameter (tSIE) | 665 |
| 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.