



## 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:** 6013329 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.