



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: 6013377 Ultima Gold LLT

Mfg.Date: 24-JUL-2023

Batch#: 97-23301

Exp.Date: 01-FEB-2025

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

Test	Results
Quench parameter (tSIE)	686
Quench parameter (tSIE) 50% water	281
Low Level background 50% water	2.36 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.