



## 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:** 6013179 Ultima Gold F

**Mfg.Date:** 31-JUL-2023

**Batch#:** 78-23311

**Exp.Date:** 01-AUG-2024

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

Test	Results
Quench parameter (tSIE)	976
Background 0-18.6 keV	16 cpm
Quench parameter (tSIE) with Soluene-350	768
Background with Soluene-350	762 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.