



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: 6013171 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.