



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: 07-AUG-2023

Batch#: 78-23321

Exp.Date: 01-SEP-2024

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

Test	Results
Quench parameter (tSIE)	968
Background 0-18.6 keV	13 cpm
Quench parameter (tSIE) with Soluene-350	780
Background with Soluene-350	597 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.