



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: 6013721 Carbo-Sorb E

Mfg.Date: 14-AUG-2023

Batch#: 99-23331

Exp.Date: 01-MAR-2026

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

Test	Results
Background	25 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.