

**Lot specific information**

**Product:** 6013721 Carbo-Sorb E

**Mfg.Date:** 09-OCT-2023

**Batch#:** 99-23411

**Exp.Date:** 01-MAY-2026

-----  
**Information**

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

Test	Results
Background	16 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.**