

**Lot specific information**

**Product:** 6013721 Carbo-Sorb E

**Mfg.Date:** 10-JUN-2024

**Batch#:** 99-24245

**Exp.Date:** 01-JAN-2027

-----  
**Information**

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

Test	Results
Background	30 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.**