

Lot specific information

Product: 6013721 Carbo-Sorb E

Mfg.Date: 18-SEP-2023

Batch#: 99-23385

Exp.Date: 01-APR-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.**

