

Lot specific information

Product: 6013721 Carbo-Sorb E

Mfg.Date: 01-APR-2023

Batch#: 99-24145

Exp.Date: 01-NOV-2026

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.**