

Lot specific information

Product: 6013721 Carbo-Sorb E

Mfg.Date: 02-JAN-2024

Batch#: 99-24015

Exp.Date: 01-AUG-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.**