

Lot specific information

Product: 6003054 Combustaid

Mfg.Date: 18-DEC-2023

Batch#: 25-23515

Exp.Date: 01-JUL-2026

Information

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

Test	Results
Refraction	1.4041

**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.**

