

Lot specific information

Product: 6013389 Emulsifier-Safe

Mfg.Date: 29-JAN-2024

Batch#: 76-24055

Exp.Date: 01-AUG-2026

Information

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

| Test | Results |
|-------------------------|---------|
| Quench parameter (tSIE) | 470 |

**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.**