

Lot specific information

Product: 6013389 Emulsifier-Safe

Mfg.Date: 19-AUG-2024

Batch#: 76-24345

Exp.Date: 01-MAR-2027

Information

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

Test	Results
Quench parameter (tSIE)	474

**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.**

