

**Lot specific information**

**Product:** 6013389 Emulsifier-Safe

**Mfg.Date:** 25-MAR-2024

**Batch#:** 76-24135

**Exp.Date:** 01-OCT-2026

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	468

**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.**

