

Lot specific information

Product: Emulsifier-Safe Mfg. Date: 20-NOV-2023
 Lot number: 76-23475 Exp. Date: 01-JUN-2026

Information

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

Test	Temperature	Results
Appearance		Clear blue fluorescent liquid
Quench parameter		75.7%
Water Holding Capacity	5 °C	≥ 1.1 ml clear
0.01 M PBS Holding Capacity	27 °C	≥ 1.9 ml clear
	20 °C	≥ 2.6 ml clear
Appearance 0.9 ml 1.0 M NaCl	27 °C	Clear within 1.5 minute

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.