

Lot specific information

Product: High Efficiency Mineral Oil Mfg. Date: 26-FEB-2024
 Lot number: 137-24095 Exp. Date: 01-OCT-2026

Information

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

Test	Temperature	Results
Appearance		Clear blue fluorescent liquid
Counting Efficiency		102.1%
Background		14 cpm

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.