

Lot specific information

Product: 6013611 MicroScint-O

Mfg.Date: 09-SEP-2024

Batch#: 86-24375

Exp.Date: 01-OCT-2025

Information

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

| Test | Results |
|--------------------------------|---------|
| Quench parameter (tSIE) | 854 |
| Counting Efficiency (TopCount) | 58.4 % |
| Background (TopCount) | 3 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.**