

Lot specific information

Product: 6013611 MicroScint-O

Mfg.Date: 29-JAN-2024

Batch#: 86-24055

Exp.Date: 01-FEB-2025

Information

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

Test	Results
Quench parameter (tSIE)	875
Counting Efficiency (TopCount)	59.8 %
Background (TopCount)	4 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.**