

Lot specific information

Product: 6013621 MicroScint-20

Mfg.Date: 13-MAR-2023

Batch#: 87-23115

Exp.Date: 01-OCT-2024

Information

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

Test	Results
Counting Efficiency (TopCount)	45.7 %
Background (TopCount)	13 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.**