

**Lot specific information**

**Product:** 6013621 MicroScint-20

**Mfg.Date:** 15-APR-2024

**Batch#:** 87-24165

**Exp.Date:** 01-NOV-2025

-----  
**Information**

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

| Test                           | Results |
|--------------------------------|---------|
| Counting Efficiency (TopCount) | 46.1 %  |
| 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.**