

Lot specific information

Product: 6NE9100 SOLVABLE

Mfg.Date: 27-NOV-2023

Batch#: 140-23485

Exp.Date: 01-JUN-2026

Information

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

Test	Results
Normality (OH-)	0.41
Background in Ultima Gold after 30 min	14 cpm
Background in Ultima Gold after 2 hrs	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.**