

Lot specific information

Product: 6NE9100 SOLVABLE

Mfg.Date: 15-JUL-2024

Batch#: 140-24295

Exp.Date: 01-FEB-2027

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	12 cpm
Background in Ultima Gold after 2 hrs	9 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.**