

Lot specific information

Product: 6003005 Hydroxide of Hyamine 10-X

Mfg.Date: 21-MAY-2024

Batch#: 2-24215

Exp.Date: 01-DEC-2026

Information

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

| Test | Results |
|--|---------|
| Chlorine content | 0.32 % |
| Normality (mg eq OH-/ g) | 1.05 |
| Quench parameter (tSIE) | 340 |
| Quench parameter Toluene scint. (tSIE) | 790 |

**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.**

