

**Lot specific information**

**Product:** 6003005 Hydroxide of Hyamine 10-X

**Mfg.Date:** 07-OCT-2024

**Batch#:** 2-24415

**Exp.Date:** 01-MAY-2027

-----  
**Information**

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

| Test                                   | Results |
|--|---------|
| Chlorine content                       | 0.30 %  |
| Normality (mg eq OH-/ g)               | 1.05    |
| Quench parameter (tSIE)                | 345     |
| Quench parameter Toluene scint. (tSIE) | 794     |

**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.**

