

**Lot specific information**

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

**Mfg.Date:** 16-OCT-2023

**Batch#:** 2-23425

**Exp.Date:** 01-MAY-2026

-----  
**Information**

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

Test	Results
Chlorine content	0.29 %
Normality (mg eq OH-/ g)	1.02
Quench parameter (tSIE)	348
Quench parameter Toluene scint. (tSIE)	796

**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.**