

Lot specific information

Product: 6003005 Hydroxide of Hyamine 10-X

Mfg.Date: 13-NOV-2023

Batch#: 2-23461

Exp.Date: 01-JUN-2026

Information

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

Test	Results
Chlorine content	0.40 %
Normality (mg eq OH-/ g)	1.03
Quench parameter (tSIE)	338
Quench parameter Toluene scint. (tSIE)	786

**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.**