



Certificate of Analysis

Revvity Health Sciences B.V.
Rigaweg 22
9723 TH Groningen
The Netherlands

Phone: + 31-50-544 59 00
SDS.Groningen@revvity.com
www.revvity.com

Product: 6003005 Hydroxide of Hyamine 10-X

Mfg.Date: 31-JUL-2023

Batch#: 2-23311

Exp.Date: 01-FEB-2026

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

Test	Results
Chlorine content	0.28 %
Normality (mg eq OH-/ g)	1.01
Quench parameter (tSIE)	345
Quench parameter Toluene scint. (tSIE)	791

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.