



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: [2N] NaOH; Sodium hydroxide

Catalog #: R007

Lot #: 210202

Size: 100ml

Tests:

| | Specification | Result | Disposition |
|-----------------------|---------------|-----------|-------------|
| Appearance - | Clear Liquid | Clear Liq | Passed |
| Specific gravity test | 0.99-1.2g/mL | 1.09 | Passed |
| pH Test | >13 | 13.9 | Passed |

Storage: Store it at RT.

This document has been produced electronically and is valid without signature

Date of Print: February 19, 2021

Date of Expiry: 02-02-2027

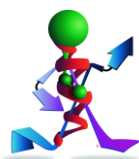
G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

NaOH ; Sodium hydroxide 2N C of A/Com#3
Rev 04.04.13/MM



think proteins! think G-Biosciences!

www.GBiosciences.com