



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Potassium hydroxide [2N]

**Catalog #:** R008

**Lot #:** 202610

**Size:** 100ml

### Tests:

Appearance - Clear Liquid	Passed	Yes / <del>No</del>
Specific gravity test	Passed	Yes / <del>No</del>
Potassium acetate 2N solution (tolerance $\pm 0.001$ )	Passed	Yes / <del>No</del>

**Storage:** Store at room temperature

*For Research Use Only*

This document has been produced electronically and is valid without signature

**Date of Print:** November 6, 2020

### **G-Biosciences / Genotech**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

KOH 2N C of A/Com#3  
Rev 10.11.13/MM



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)