



Certificate of Analysis

Product: Sodium Acetate

Catalog #: RC-090

Quantity: 1kg

Lot#: 211002

Mol Formula: CH₃COONa

MW: 82.03

Tests:

	<u>Min</u>	<u>Max</u>
Insolubles		0.01%
Loss on Drying		1.0%
pH (5%, water) @25C	7.0	9.2
Chloride		0.002%
Phosphate		0.001%
Sulfate		0.003%
Calcium		0.005%
Heavy Metals (as Pb)		0.001%
Iron		0.001%
Purity	99.0%	
Magnesium		0.002%

Storage Temp: Store at room temperature, protect from Moisture.

This document has been produced electronically and is valid without signature

Date of Print: February 19, 2021

G-Biosciences / Genotech

9800 Page Avenue, St. Louis, MO 63132, USA

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

Sodium Acetate C of A/Common# 3/Rev 09.25.12-IA

