



### Certificate of Analysis

**Product:** Potassium Iodide  
**Catalog #:** RC-151  
**Size:** 500g  
**Lot#:** 211207  
**Re-test date:** 07/2026  
**Mol Formula:** KI  
**MW:** 166.00  
**CAS#:** 7681-11-0

	Min	Max	Actual	Units	Disposition
Purity	99.0		100.00	%	PASS
pH (5% in Water) at 25°C	6.0	9.2	6.61		PASS
Loss on Drying		0.2	0.022	%	PASS
Insoluble(s)		0.005	<0.0005	%	PASS
Iodate		0.0003	<0.0003	%	PASS
Chloride & Bromide		0.01	0.0075	%	PASS
Phosphate		0.001	0.0002	%	PASS
Sulfate		0.005	0.0025	%	PASS
Barium		0.002	<0.0001	%	PASS
Heavy Metals (as Pb)		0.0005	0.0001	%	PASS
Iron		0.0003	0.0001	%	PASS
Sodium		0.005	0.0039	%	PASS
Calcium		0.002	0.0004	%	PASS
Magnesium		0.001	<0.0001	%	PASS

**Storage Temp:** Store at Room Temp. Protect from Moisture

This document has been produced electronically and is valid without signature

**Date of Print:** November 9, 2021

**G-Biosciences / Genotech**  
 9800 Page Avenue, St. Louis, MO 63132, USA  
 Ph: 314-991-6034, Fax: 314-991-1504

Potassium Iodide C of A/Com#3  
Rev 02.25.13/MM

