



## Certificate of Analysis

**Product:** Cesium chloride  
**Catalog #:** RC-034  
**Quantity:** 100g  
**Lot#:** 121007  
**Molecular Formula:** CsCl  
**Molecular Weight:** 168.36  
**Appearance:** Fine white powder  
**Purity:** > 99.9%

### Tests:

UV absorbance (50% in Water, equilibrated overnight)

260nm	<0.02
280nm	<0.02
405nm	<0.01

Chemical Analysis, ppm

Rb	32	K	<1	Na	5	Li	<1
Insol.	<100	Al	<1	Ba	<1	Ca	<1
Cr	<1	Cu	<1	Fe	<1	Mg	<1
Mn	<1	Ni	<1	Pb	<1	Si	<1

**Storage:** Store at room temperature. Protect from moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** July 19, 2017

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

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

Cesium chloride 500g, RC-035 C of A/Com#3  
Rev 07.11.12/IA

