



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:** Silica Magnetic Beads  
**Catalog #:** 786-915, 786-916, 786-917, 786-988  
**Lot#:** 201012  
**Re-test date:** 09/2022

### Tests:

Appearance (color)  
Concentration  
Particle size  
DNA binding test  
(>10ug in 50ug/ml salmon sperm DNA)

### Disposition

Black to brown  
50.1mg beads/ml  
3.19um  
20.1ug DNA/ml beads

This document has been produced electronically and is valid without signature

**Date of Print:** March 16, 2020

### **Genotech/GBiosciences**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504  
[www.GBiosciences.com](http://www.GBiosciences.com)

Silica magnetic beads C of A/Com#3  
Created 6/5/18 MN



think proteins! think G-Biosciences!

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