



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:** Carboxyl Magnetic Beads

**Catalog #:** 786-908

**Lot#:** 180901

**Re-test date:** 01/2023

**Tests:**

Appearance (color)  
Concentration  
Particle size

**Disposition**

Black to brown  
50.7mg beads/ml  
3.64um

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)



think proteins! think G-Biosciences!

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