



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: Silica Magnetic Beads
Catalog #: 786-915, 786-916, 786-917, 786-988
Lot#: 201510
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

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



think proteins! think G-Biosciences!

www.GBiosciences.com