



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

Certificate of Analysis

Product: Bicinchoninic Acid (BCA) Reagent A

Catalog #: 786-846

Size: 1 Liter

Lot #: 180712

Expiry: 11\2020

QC Tests:

Curve R² value (≥ 0.99)

Passed

Yes / ~~No~~

BCA Blank (< 0.08)

Passed

Yes / ~~No~~

Storage: RT

For Research Use Only

This document has been produced electronically and is valid without signature

Date of Print: October 23, 2019

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

Rev 08.25.16/MM



think proteins! think G-Biosciences!

www.GBiosciences.com