



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

Certificate of Analysis

Product: Bovine Gamma Globulin Standard; 2mg/ml

Catalog #: 786-007

Lot #: 121408B

Appearance: Clear Liquid

Tests:

Concentration checked [2mg/ml] Passed Yes / ~~No~~

Protease Test –

Assay was performed using G-Biosciences's Protease Screening kit

No detectable protease activity found Passed Yes / ~~No~~

This document has been produced electronically and is valid without signature

Date of Print: January 12, 2017

G-Biosciences / Genotech

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

Bovine gamma globulin Standard 2mg/ml C of A/Rev 08.27.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com