



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:** Bovine Gamma Globulin Standard; 2mg/ml

**Catalog #:** 786-007

**Lot #:** 111205

**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:** December 6, 2011

#### **G-Biosciences / Genotech**

92 Weldon Parkway  
St. Louis, MO 63043, USA  
Fax: 314-991-1504

Bovine gamma globulin Standard 2mg/ml C of A/Rev 03.31.10/IA



think proteins! think G-Biosciences!

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