



## Certificate of Analysis

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

**Catalog #:** 786-007

**Lot #:** 052701

**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:** July 8<sup>th</sup>, 2009

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

92 Weldon Parkway, St. Louis, MO 63043, USA

Fax: 314-991-1504

Bovine gamma globulin Standard 2mg/ml CofA/Com3 Rev 07.08.09-SA/MM/IA



**think proteins! think G-Biosciences!**

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