



Certificate of Analysis

Product: Bovine Gamma Globulin Standard; 2mg/ml

Catalog #: 786-007

Lot #: 092007

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: November 30, 2010

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