



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

CERTIFICATE OF ANALYSIS

Product: Reversible Zinc Stain[™]

Catalog #: 786-32ZN

Lot #: 042210

Q.C. Tests

Tests were performed by running SDS-PAGE, using serially diluted BSA up to 10ng and G-Biosciences's in-house QC protocol. After electrophoresis, gel was stained with 5 Minute Reversible Stain[™] Zinc.

Protein Detection Sensitivity Test Passed

OK

Yes / ~~No~~

Storage: Store at RT.

This document has been produced electronically and is valid without a signature

Date of print: November 29, 2010

G-Biosciences / Genotech

92 Weldon Pkwy

St. Louis. MO 63043, USA

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

Reversible Zinc Stain C of A/Com# 3
Rev 12.16.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com