



CERTIFICATE OF ANALYSIS

Product Name: **BLOT-Fast Stain™**

Catalog #: **786-34**

Lot #: **131604**

Q.C. Tests

Tests performed using protein (BSA) dot-blot. Serially diluted BSA up to 0.5ng/μl in Protein dotMetric dilution buffer was applied on a nitrocellulose membrane strip, using G-Biosciences's application device (Cat.# 786-63). The Test Strip was developed as per G-Biosciences's protocol.

Spot developed in 2-3 minutes – O.K.

Detection sensitivity (up to 0.5ng BSA) – O.K.

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

Date of print: April 24, 2013

G-Biosciences / Genotech

9800 Page Avenue, St. Louis. MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1405

BlotFast Stain /C of A/Com#3 Rev 04.24.13/MM

