



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

Certificate of Analysis

Product: Blue Dextran Loading Dye

Catalog #: 786-557

Lot #: 101410

Size: 3 x 1.5ml

Tests:

Specific gravity test	Passed	Yes / No
Dye Run Test	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
Ph: 314-991-6034, Fax: 314-991-1504

Blue Dextran Loading Dye
Rev 10.15.10-IA



think proteins! think G-Biosciences!

www.GBiosciences.com