



CERTIFICATE OF ANALYSIS

Product: BLOK™ Casein in PBS

Cat. #: 786-194

Lot #: 072702

TESTS

Product tested with either Western blot/ Dot Blot analysis.

BLOK™ Casein in PBS protocol (Cat. # 786-194) was followed.

Product tested with both PVDF and Nitrocellulose (NC) membranes.

METHODS

G-Bioscience Lab procedure

SDS-PAGE & Western Blot or Dot Blot Method

Detection method, G-Bioscience Chemiluminescence kit (*femto*-LUCENT, Cat. # 786-10)

Critical Agents: Multiple tissue lysates / Western Blot / Dot Blot

Probe- actin antibodies (primary) polyclonal, using HRP-labeled antibodies (secondary)

Type of membrane

PVDF

Nitrocellulose (NC)

Background Clear

Yes / ~~Ne~~

Yes / ~~Ne~~

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

Date of print: June 13, 2007

G-Biosciences / Genotech

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

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

BLOK Casein in PBS CofA/Com#3 Rev 12.18.08-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com