



CERTIFICATE OF ANALYSIS

Product: BLOK™ Casein in PBS

Cat. #: 786-194

Lot #: 142105

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 / ~~No~~

Yes / ~~No~~

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

Date of print: May 22, 2014

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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

BLOK Casein in PBS C of A/Com#3
Rev 05.10.12/1A



think proteins! think G-Biosciences!

www.GBiosciences.com