



CERTIFICATE OF ANALYSIS

Product: BLOK™ BSA in PBS [10X]

Cat. #: 786-195

Lot #: 172011

TESTS

Product tested with either Western blot/ Dot Blot analysis.

BLOK™ BSA in PBS protocol (Cat # 786-195) 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: July 12, 2018

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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

BLOK BSA in PBS C of A/Com#3
Rev 07.02.12/1A



think proteins! think G-Biosciences!

www.GBiosciences.com