



## CERTIFICATE OF ANALYSIS

**Product:** BLOK™ Casein in PBS

**Cat. #:** 786-194

**Lot #:** 202011

**Re-test date:** 11/2023

### 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:** December 2, 2020

**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

