



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** BLOK™ Casein in TBS

**Cat. #:** 786-196

**Lot #:** 202301

**Re-test date:** 02/2021

### TESTS

Product tested with either Western blot/ Dot Blot analysis.

BLOK™ Casein in TBS protocol 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)

<u>Type of membrane</u>	<u>Background Clear</u>	
PVDF		Yes / <del>No</del>
Nitrocellulose (NC)		Yes / <del>No</del>

**Storage:** Store at 4°C.

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

**Date of print:** March 4, 2020

### G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

Fax: 314-991-1504



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)