



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com



Certificate of Analysis

Product: BLOK™ Blotto in TBS

Cat. #: 786-192

Lot #: 202007

Re-test date: 07/2021

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)

Type of membrane

PVDF

Nitrocellulose (NC)

Background Clear

Yes / ~~No~~

Yes / ~~No~~

Storage: Store at 4°C.

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

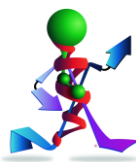
Date of print: July 30, 2020

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

Fax: 314-991-1504



think proteins! think G-Biosciences!

www.GBiosciences.com