



Certificate of Analysis

Product: BLOK™ Casein in TBS

Cat. #: 786-196

Lot #: 181903

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 / No
Nitrocellulose (NC)		Yes / No

Storage: Store at 4⁰C.

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

Date of print: May 4, 2018

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Fax: 314-991-1504

