



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

### CERTIFICATE OF ANALYSIS / DATASHEET

**Product:** Protein Free Blocking Buffer PBST

**Catalog #:** 786-665

**Size:** 500ml

**Lot #:** 203010

#### TESTS

- Product tested with either Western blot/ Dot Blot analysis.
- Protein Free Blocking Buffer protocol (Cat. # 786-665) was followed.
- Product tested with both PVDF and Nitrocellulose (NC) membranes.

#### METHODS

G-Bioscience Lab procedures

- Detection method – GBiosciences Protein Estimation kit (*NI Protein Assay Kit, Cat# 786-005*)
- SDS-PAGE of total tissue protein & Western Blot or Dot Blot Method
- Detection of protein Actin on the western blot or dot blot. Method - *G-Bioscience Chemiluminescence kit (femto-LUCENT kit, Cat. # 786-10)*

#### Critical Agents/Tests:

- Multiple tissue lysates (total protein) for Western Blot / Dot Blot
- Probe- Actin antibodies (*primary*) polyclonal, using HRP-labeled antibodies (*secondary*)

#### Type of membrane tested

PVDF  
Nitrocellulose (NC)

#### Background Clear

Yes / ~~No~~  
Yes / ~~No~~

**Storage:** Store at 4°C.

**Shelf Life:** One year, if used and stored as recommended.

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

Date of print: December 5, 2018

**Geno Technology, Inc. /G-Biosciences**

9800 Page Avenue, St. Louis, MO 63132, USA

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



think proteins! think G-Biosciences!

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