

## CERTIFICATE OF ANALYSIS

**Product Name:** Blot-Quick Blocker™

**Cat #** 786-011

**Lot #:** 201103

**Re-test date:** 03/2025

### TESTS

BlotQuickBlocker protocol (Cat# 786-011) was followed. Product tested with either Western blot/ Dot Blot analysis with both PVDF and /or Nitrocellulose (NC) membranes.

### METHODS

For Dot blot, *Genotech* Lab procedure

For SDS-PAGE & Western Blot, *Genotech* Lab procedure

For Detection method, *Genotech* Chemiluminescence kit (*femto-LUCENT* Kit, Cat.786-10)

Critical Agents: Mouse Liver tissue lysate, Probe- polyclonal Actin antibodies (primary) using, HRP-labeled secondary antibodies.

~~Dot Blot~~ / Western Blot (indicated method used)

<u>Method</u>	<u>Membrane</u>	<u>Background Clear</u>
Western Blot probed for Actin	PVDF / NC	Yes / <del>No</del>

**Storage:** Store at room temp.

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

**Date of print:** March 30, 2020

***G-Biosciences / Genotech***  
***9800 Page Avenue***  
St. Louis, MO 63132, USA  
[www.GBiosciences.com](http://www.GBiosciences.com)