



296PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Nitrocellulose Membrane

(Cat. # 786-018NC, 786-056NC)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

G-Biosciences Nitrocellulose Membrane is made up high quality pure nitrocellulose, which ensures maximum binding capacity with low background. It does not contain any support fabric or detergents and is compatible with standard transfer conditions. It is ideal for Western, Southern, and Northern blotting, Dot/Slot blotting, Amino Acid Analysis etc. The nitrocellulose membranes are precut and Cat. # 786-018NC are supplied sandwiched with filter paper.

BINDING CAPACITY/INTERACTIONS

- ~200µg protein/cm²
- Hydrophobic and electrostatic binding interactions

ITEM(S) SUPPLIED

Description	Size: 8.5 x 7.5cm Cat. # 786-018NC	Size: 10 x 10cm Cat. # 786-056NC
Nitrocellulose Membrane	20	10
Pore size (µm)	0.2 µm	0.2 µm
Filter Paper Padding	Yes	No

STORAGE CONDITIONS

Nitrocellulose membranes must be stored properly to ensure maximum shelf life. Store at room temperature away. Keep away from chemical vapors, which may partially dissolve the nitrocellulose thereby disrupting the pore structure. Avoid direct sunlight, which may cause nitrocellulose membrane to dry out and become brittle.

WARNING: Nitrocellulose membranes are **highly flammable**. Keep away from heat and open flame. Flash point is approximately 200°C.

PROPER HANDLING

Gloves must always be worn when handling nitrocellulose membrane. Natural oil from fingers will prevent proper wetting of the nitrocellulose membrane. Also, proteins from the fingers/hands will bind instantly to the nitrocellulose and cause background spots when membrane is developed.

APPLICATIONS

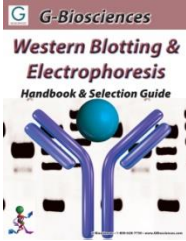
Nitrocellulose Membranes are suitable for Western, Southern, and Northern Blotting, Dot/Slot blotting, Amino Acid Analysis etc.

PREPARATION BEFORE USE

To ensure complete transfer with maximum binding and low background, the nitrocellulose must be thoroughly soaked in deionized water prior to transfer. Proteins and nucleic acids will not adsorb to dry areas.

RELATED PRODUCTS

Download our Western Blotting Handbook.



<http://info.gbiosciences.com/complete-western-blot-handbook--selection-guide>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 9/29/2015 CMH



www.GBiosciences.com