



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

A Geno Technology, Inc. (USA) brand name

High Molecular Weight Western Transfer Buffer

For the Improved Transfer of >70kD Proteins

(Cat. # 786-423)



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

INTRODUCTION

Based on Efficient™ Western Transfer Buffer, the High Molecular Weight Transfer buffer is designed to facilitate the transfer of notoriously difficult high molecular weight proteins (>70kDa) during Western blotting.

ITEM(S) SUPPLIED (Cat. # 786-423)

Description	Size
High Molecular Weight Western Transfer Buffer [5X]	1L

STORAGE CONDITIONS

It is shipped at ambient temperature. Store it at room temperature upon arrival. Stable for one year when stored and used properly.

ADDITIONAL ITEM(S) REQUIRED

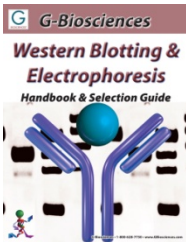
- Deionized water
- Methanol

PROTOCOL

1. Take 200ml 5X High Molecular Weight Western Transfer Buffer and add to 600ml deionized or ultrapure water.
2. Add 200ml methanol to make a final volume of 1 liter. Methanol is 20% (v/v).
3. For best results, chill the 1X High Molecular Weight Western Transfer Buffer to 4°C before use.
4. Use the 1X High Molecular Weight Western Transfer Buffer as per the Western blotting equipment instructions. Suitable for wet and semi-dry transfer.

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: 2/22/2013 CMH

This page is intentionally left blank



www.GBiosciences.com