



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

A Geno Technology, Inc. (USA) brand name

Efficient™ Western Transfer Buffer

High Efficiency Transfer Buffer for Western Blotting

(Cat. # 786-019, 786-019G)



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

INTRODUCTION

Efficient™ Western Transfer Buffer is a high efficiency transfer buffer for Western blotting. The pre-made and ready-to-use [20X] transfer buffer is prepared for optimal conductivity and efficient protein transfer without generating excessive heat or transfer distortion.

ITEM(S) SUPPLIED (Cat. # 786-019)

Description	Size
Efficient™ Western Transfer Buffer [20X]	1L
Efficient™ Western Transfer Buffer [20X]	1 gallon

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

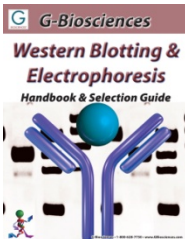
- Deionized water
- Methanol
- SDS (optional)

PROTOCOL

1. Take 50ml 20X Efficient™ Western Transfer Buffer and add to 750ml deionized or ultrapure water.
2. Add 200ml methanol to make a final volume of 1 liter. Methanol is 20% (v/v).
OPTIONAL: Add SDS to a final concentration of 0.01% (w/v) to facilitate transfer.
3. For best results, chill the 1X Efficient™ Western Transfer Buffer to 4°C before use.
4. Use the 1X Efficient™ 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.

This page is intentionally left blank



www.GBiosciences.com