



996PR-03

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

A Geno Technology, Inc. (USA) brand name

2D-Xtract™

Chaotropic Extraction Buffer

(Cat. # 786-501)



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

INTRODUCTION

2D-Xtract™ is a dry, urea-based and ready-to-use protein solubilization buffer. 2D-Xtract™ is a proprietary modification of urea based chaotropic extraction buffers. Normal urea based buffers solubilize proteins effectively however can modify the native charge of the proteins, due to carbamylation. A process that modifies amino groups, inducing changes in the isoelectric point of proteins leading to artifactual 2D results.

2D-Xtract™ is a dry, pre-mixed formulation of urea, thiourea, CHAPS and non detergent sulfobetaine (ND SB 201) for maximum solubilizing strength. The dry format allows researchers to freshly rehydrate as much or as little as is required and therefore prevent urea induced carbamylation.

ITEM(S) SUPPLIED (Cat. # 786-501)

Description	Size
2D-Xtract™	25gm
DILUENT-III	30ml

STORAGE CONDITION

The kit is shipped at ambient temperature. Upon arrival, store it as specified on the bottle labels.

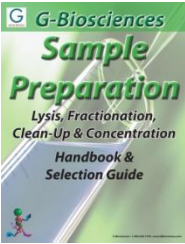
INSTRUCTION FOR USE

For optimal performance of 2D-Xtract™, vigorously shake the powder before use and prepare fresh immediately prior to solubilization.

1. Vigorously shake the 2D-Xtract™ bottle for 15 seconds.
2. For 100µl final 2D-Xtract™ buffer, weigh 50mg dry 2D-Xtract™ in to a suitable tube.
3. Add 55µl Diluent-III for every 50mg of 2D-Xtract™ and alternatively vortex and incubate at room temperature until a clear solution is achieved.
4. If additional reagents are required, including reducing agents, inhibitors, carrier ampholytes and dyes, these are added at this point. Vortex to mix.
5. Use the prepared 2D-Xtract solution immediately and discard any unused buffer.

RELATED PRODUCTS

Download our Sample Preparation Handbook.



<http://info.gbiosciences.com/complete-protein-sample-preparation-handbook>

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

Last saved: 7/3/2014 CMH



www.GBiosciences.com