



339PR

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

A Geno Technology, Inc. (USA) brand name

Optimizer Buffer™ - II

(Cat. #BKC-05)



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

INTRODUCTION

Optimizer Buffers™ provide the optimal conditions for specific protein reactions, including protein labeling, modification and cross-linking reactions.

Optimizer Buffer™ -II has a high pH and is specifically designed for the sulfhydryl reactive haloacetyls, such as iodoacetyl.

ITEM(S) SUPPLIED (Cat. # BKC-05)

Description	Size
5X Optimizer Buffer™ -II	2 x 25ml

STORAGE CONDITIONS

It is shipped at ambient temperature. Store refrigerated at 4°C. This product is stable for 1 year.

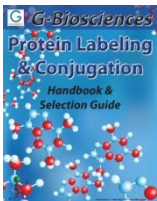
INSTRUCTIONS FOR USE

Prior to experiment, the buffer is diluted to a 1X solution (i.e. 2ml Optimizer Buffer™ -II in 8ml distilled water).

Optimizer Buffer™ -II is designed to provide the optimal conditions for sulfhydryl reactions involving iodoacetyl groups.

RELATED PRODUCTS

Download our Protein Labeling & Conjugation Handbook.



<http://info2.gbiosciences.com/complete-protein-labeling-conjugation-handbook>

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

Last saved: 06/12/2014 ADB

This page is intentionally left blank.



www.GBiosciences.com