



341PR

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

A Geno Technology, Inc. (USA) brand name

Optimizer Buffer™ -III

(Cat. # BKC-06)



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™-III is designed for sulfhydryl reactive groups, maleimides and pyridyl disulfides, and has reagents that prevent oxidation of free sulfhydryls.

ITEM(S) SUPPLIED (Cat. # BKC-06)

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

STORAGE CONDITIONS

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

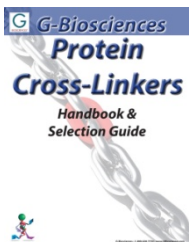
INSTRUCTIONS FOR USE:

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

Optimizer Buffer™ -III is designed to provide the optimal conditions for sulfhydryl reactions involving maleimide and pyridyl disulfide groups.

RELATED PRODUCTS

Download our Protein Cross-linkers Handbook.



<http://info.gbiosciences.com/complete-protein-cross-linkers-handbook/>

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

Last saved: 12/14/2012 CMH

This page is intentionally left blank



www.GBiosciences.com