



343PR

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

A Geno Technology, Inc. (USA) brand name

Optimizer Buffer™ -IV

(Cat. # BKC-07)



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™-IV is designed to provide the optimal conditions necessary for carboxyl reactive reagents.

ITEM(S) SUPPLIED (Cat. # BKC-07)

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

STORAGE CONDITIONS

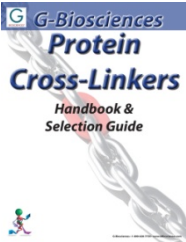
It is shipped at ambient temperature. Store refrigerated at 4°C. This product is stable for 1 year.

INSTRUCTIONS FOR USE:

1. Prior to experiment, the buffer is diluted to a 1X solution (i.e. 2ml Optimizer Buffer™ in 8ml distilled water)
2. Optimizer Buffer™-IV is designed to provide the optimal conditions for reactions with carboxyl 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: 1/24/2013 CMH

This page is intentionally left blank



www.GBiosciences.com