



347PR

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

A Geno Technology, Inc. (USA) brand name

Optimizer Buffer™ -VI

(Cat. #BKC-09)



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

INTRODUCTION 3

ITEM(s) SUPPLIED (Cat, # BKC-09) 3

STORAGE CONDITIONS 3

INSTRUCTIONS FOR USE 3

Related Products 3

INTRODUCTION

Optimizer Buffers™ provide the optimal conditions for specific protein reactions, including protein labeling, modification and cross-linking reactions. Optimizer Buffer™ -VI is an alkaline buffer specific for the formation of Schiff's base between amines and glyoxal groups.

ITEM(S) SUPPLIED (Cat, # BKC-09)

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

STORAGE CONDITIONS

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

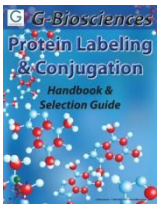
INSTRUCTIONS FOR USE

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

Optimizer Buffer™ -VI is designed to provide the optimal conditions for reactions requiring a strong alkaline buffer, particularly in the formation of Schiff's bases for coupling of amines to glyoxal 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 3/21/2014 SM



www.GBiosciences.com