



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

A Geno Technology, Inc. (USA) brand name

Optimizer Buffer™-VIII [5X]

(Cat. # BKC-11)



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™-VIII [5X] is designed for use in citraconic anhydride modification. Citraconic Anhydride reacts with a primary amine to create an amide linkage and a terminal carboxylate. The linkage is stable at neutral to alkaline pH and is hydrolyzed in acidic conditions (pH3-4).

ITEM(S) SUPPLIED (Cat. # BKC-11)

Description	Size
Optimizer Buffer™-VIII [5X]	2 x 25ml

STORAGE CONDITIONS

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

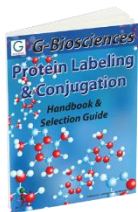
INSTRUCTIONS FOR USE

Prior to experiment, the buffer is diluted to a 1X solution (i.e. 2ml Optimizer Buffer™-VIII [5X] in 8ml distilled water).

Optimizer Buffer™-VIII is designed for use in citraconic anhydride modification. Dissolve the amine containing sample or protein in 1X Optimizer Buffer™-VIII buffer. Add 5-10 fold molar excess of citraconic anhydride over the amount of amines. Incubate at room temperature for 1-2 hrs or overnight at 4°C.

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.

This page is intentionally left blank.



www.GBiosciences.com