



## Optimizer Buffer™ -III

### 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:

- Optimizer Buffers (Cat. # BKC-04 – BKC-09):** Five different Optimizer Buffers™ are available, that are specific for other protein reactions.
- Tube-O-Reactor (Cat. # 786-027):** A complete kit for the exchange of buffers for Optimizer™ Buffer utilizing Tube-O-Dialyzer™. Ensures no sample loss.

**NOTE:** For other related products, visit our web site at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Rev 05.28.09/MM

