



626PR-01

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

A Geno Technology, Inc. (USA) brand name

Bolton-Hunter Reagent

Sulfo-SHPP

(Cat. # BC92)



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

INTRODUCTION

G-Biosciences water-soluble Bolton-Hunter Reagent (Sulfo-SHPP) conjugates tyrosine-like groups to end-terminal α -amino groups or ϵ -amino groups of lysine to increase the number of tyrosyl groups that can be iodinated by iodine-125 labeling procedures.

Radioactive iodine (^{125}I) is routinely used by researchers to label proteins. The iodination of proteins can be performed enzymatically or chemically. The Bolton-Hunter reagent is designed to aid the labeling of proteins with radioactive iodine.

ITEM(S) SUPPLIED

Cat. #	Description	Size
BC92	Sulfo SHPP (Water soluble Bolton-Hunter reagent)	100mg

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at -20°C .

SPECIFICATIONS

- Synonyms: Sulfosuccinimidyl-3-(4-hydroxyphenyl) propionate; Sulfo SHPP; Water Soluble Bolton-Hunter Reagent
- CAS Number: 106827-57-0
- Molecular Weight: 365.3
- Ideal for proteins with masked or no tyrosine residues
- Optimal reaction at pH 8.5
- Water soluble

PROTOCOL FOR IODINATION OF A PURIFIED PROTEIN

The following protocol is a guide to modifying a protein before iodination.

Material required

- 200mM borate buffer, pH9.0 (Modification buffer)
- Protein
- Phosphate Buffered Saline [1X] (Cat. # 786-377)

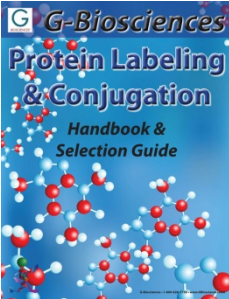
1. Dissolve 190mg of protein in 2ml 200mM borate buffer, pH9.0.
2. Immediately before use dissolve 5mg of Water-Soluble Bolton-Hunter Reagent in 1ml 200mM borate buffer, pH9.0.

NOTE: Alternatively, add 0.5mg directly to the protein solution and then go to step 4.

3. Add 100µl sulfoSHPP solution to the protein solution.
4. Incubate on ice for 3 hours with periodic mixing.
5. Remove the non-reacted sulfo-SHPP by dialyzing against PBS or other suitable buffer. We recommend using our Tube-O-DIALYZER™ to ensure 100% sample recovery
6. Radiolabel with ¹²⁵I using your established procedures.

RELATED PRODUCTS

Download our Protein Conjugation and Labeling 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: 5/19/2015 CMH



www.GBiosciences.com