



319PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# OneQuant™ HOOK™ Sulfo-NHS-LC-Biotin

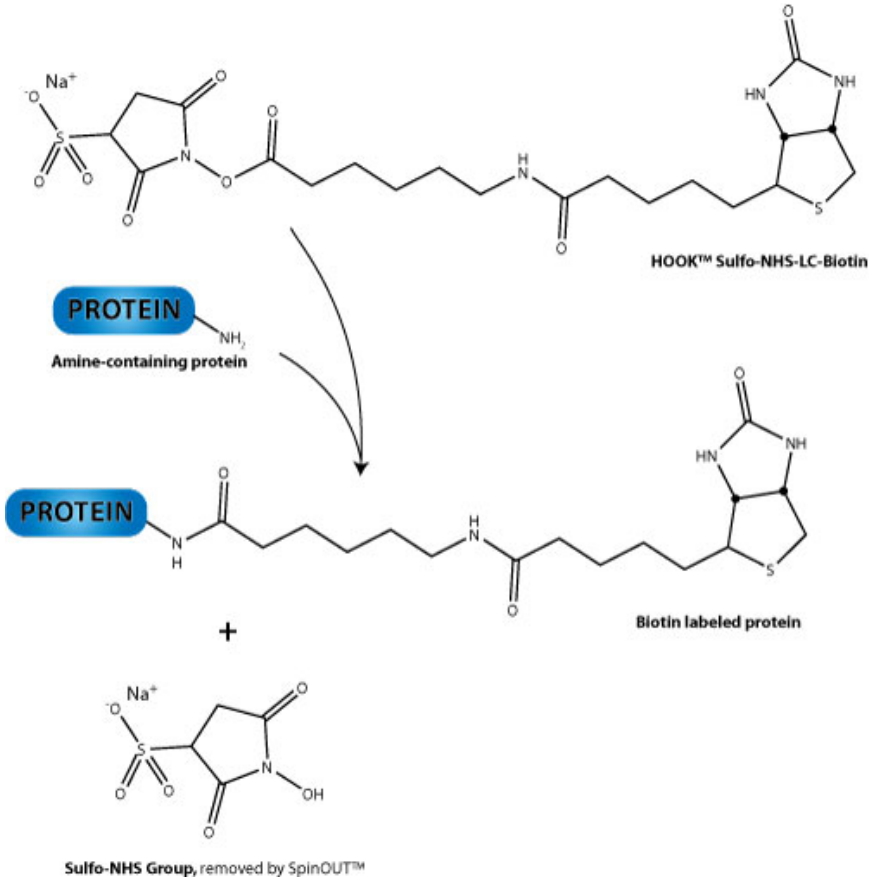
(Cat. # 786-084)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Sulfo-NHS-LC-Biotin is a commonly used amine reactive biotin-labeling reagent. OneQuant™ Sulfo-NHS-LC-Biotin are single aliquots of Sulfo-NHS-LC-Biotin reagent in convenient, premeasured microtubes. Each microtube contains 1mg of the biotin reagent. Simply dissolve one vial with appropriate solvent and add to the reaction.



## SPECIFICATIONS

Chemical name: Sulfosuccinimidyl-6-(biotinamido)hexanoate

Formula:  $\text{C}_{20}\text{H}_{29}\text{N}_4\text{O}_9\text{S}_2\text{Na}$

Molecular Weight: 556.59

CAS No: 127062-22-0

Spacer Arm Length: 22.4Å

Reaction pH: 7-9

Water soluble

## ITEM(S) SUPPLIED (Cat. #: 786-084)

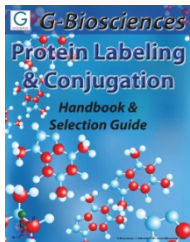
Description	Size
OneQuant™ Sulfo-NHS-LC-Biotin	1mg/vial, 1 strip of 8 vials

## STORAGE CONDITIONS

It is shipped at ambient temperature. Upon receipt store product at 4°C protected from moisture. Stable for one year.

## RELATED PRODUCTS

Download our Protein Conjugation and Labeling Handbook.



<http://info.gbiosciences.com/complete-protein-labeling-conjugation-handbook>

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

Last saved: 10/8/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)