



007PR

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 ActiveHOOK™ BSA*

(Cat. #786-093)



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

INTRODUCTION ..... 3

ITEMS SUPPLIED ..... 3

STORAGE CONDITIONS ..... 3

PROTOCOL ..... 3

RELATED PRODUCTS ..... 3

## INTRODUCTION

*OneQuant ActiveHOOK™* BSA is a pre-activated form of immunological grade BSA (MW 67kD) that has been activated by sulfoSMCC and is ready for direct coupling with peptides through the cysteine residues. Immunological grade BSA is a high purity grade bovine serum albumin (BSA) prepared for immunotechnology research applications such as a carrier protein for antibody production. *ActiveHOOK™* BSA is supplied lyophilized in *OneQuant™* format; 2mg/vial.

## ITEMS SUPPLIED

Cat. #	Description	Size
786-093	<i>OneQuant™ ActiveHOOK™</i> BSA	8 x 2mg/vial

## STORAGE CONDITIONS

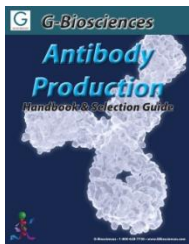
It is shipped at ambient temperature. Upon receiving, store at 4°C.

## PROTOCOL

Reconstitute with pure water or a buffer of choice to desired concentration. See related products for additional protocol information.

## RELATED PRODUCTS

Download our Antibody Production Handbook.



<http://info2.gbiosciences.com/complete-antibody-production-handbook>

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

Last saved: 3/4/20142014 TNN



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