



BIOSCIENCES®

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

## OneQuant ActiveHOOK™ BSA

### 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.

### ITEM(S) SUPPLIED

Cat. #	Description	Size
786-093	OneQuant™ ActiveHOOK™ 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

- HOOK™ Peptide Coupling Kits:** Kits for coupling peptides to proteins, available as an amine or sulfhydryl reactive kit, supplied with or without carrier proteins.
- Carrier Proteins:** All available as OneQuant™ (8x2mg/vial) or as 10mg vials.
  - BSA (Cat.# 786-090/786-086)** (bovine serum albumin) An immunological grade carrier protein.
  - KLH (Cat.# 786-091/786-088)** (keyhole limpet hemocyanin) An immunological grade carrier protein.
  - HyperCarrier™ (Cat.# 786-091/786-096)** an activated, more immunogenic form of bovine serum albumin.
- ActiveHOOK™ Carrier Proteins:** Pre-activated carrier proteins. Ready for direct coupling to peptide sulfhydryl groups. All available as OneQuant™ (8x2mg/vial) or as 10mg vials.
  - ActiveHOOK™ KLH (Cat.# 786-094/786-089)** (keyhole limpet hemocyanin) Pre-activated immunological grade carrier protein.
  - ActiveHOOK™ HyperCarrier™ (Cat.# 786-095/786-097)** an activated, more immunogenic form of bovine serum albumin.

**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



think proteins! think G-Biosciences!

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