



## HyperCarrier™

### INTRODUCTION

HyperCarrier™ is a cationized form of bovine serum albumin (MW: 69kD). Cationized proteins are taken into antigen presenting cells (APC) quicker and more efficiently than normal proteins. The positive charges on the cationized protein give it greater affinity for the negatively charged cell surface of the APCs. The cationized protein is internalized via activated receptor mediated endocytosis as opposed to pinocytosis (*Border et al. (1982) J. Clin. Invest. 69:451*). The peptides are coupled to carrier proteins with the use of protein cross-linkers, normally through primary amine residues. HyperCarrier™ is supplied lyophilized in a 10mg/vial size format.

### ITEMS INCLUDED

Cat. #	Description	Size
786-096	HyperCarrier™	1 x 10mg

### STORAGE CONDITION

Shipped at ambient temperature. Upon receiving, store at 4°C.

### PROTOCOL

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

### RELATED PRODUCTS

- 1. HOOK™ Peptide Coupling Kits:** Kits for coupling peptides to proteins, available as an amine or sulfhydryl reactive kit, supplied with or without carrier proteins.
- 2. Carrier Proteins:** All available as OneQuant™ (8x2mg/vial) or as 10mg vials.
- 3. BSA (Cat# 786-090/786-086)** (bovine serum albumin) An immunological grade carrier protein.
- 4. KLH (Cat# 786-091/786-088)** (keyhole limpet hemocyanin) An immunological grade carrier protein.
- 5. 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.
- 6. ActiveHOOK™ BSA (Cat# 786-093/786-087)** (bovine serum albumin) Pre-activated immunological grade carrier protein.
- 7. ActiveHOOK™ KLH (Cat# 786-094/786-089)** (keyhole limpet hemocyanin) Pre-activated immunological grade carrier protein.
- 8. ActiveHOOK™ HyperCarrier™ (Cat# 786-095/786-097)** Pre-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.

LU 11.05.08-SA/MM/IA

