



009PR

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

A Geno Technology, Inc. (USA) brand name

ActiveHOOK™ HyperCarrier™

(Cat. # 786-097)



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

INTRODUCTION

ActiveHOOK™ HyperCarrier™ is a pre-activated form of HyperCarrier™ that has been activated by sulfoSMCC and is ready for direct coupling with peptides through the cysteine residues. HyperCarrier™ is a cationized form of bovine serum albumin. 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). ActiveHOOK™ HyperCarrier™ is supplied lyophilized in 10mg/vial size.

ITEM(S) SUPPLIED (Cat. # 786-097)

Description	Size
ActiveHOOK™ HyperCarrier™	1 x 10mg

STORAGE CONDITION

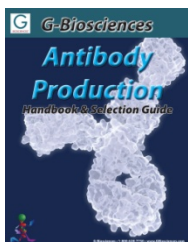
Shipped at ambient temperature. Upon receiving, store it at 4°C.

PROTOCOL

Reconstitute with pure water or a buffer of choice to desired concentration.

RELATED PRODUCTS

Download our Antibody Production Handbook



<http://info.gbiosciences.com/complete-Antibody-Production-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 7/26/2012 CMH

This page is intentionally left blank



www.GBiosciences.com