



227PR

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

A Geno Technology, Inc. (USA) brand name

OneQuant™ HyperCarrier™

(Cat. # 786-092)



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

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. OneQuant HyperCarrier™ is supplied lyophilized in 2mg/vial size format.

ITEM(S) SUPPLIED (Cat. # 786-092)

Description	Size
OneQuant™ HyperCarrier™	8 x 2mg/vial

STORAGE CONDITIONS

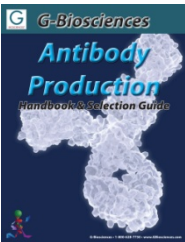
It is shipped at ambient temperature. Upon receiving store at 4°C.

PROTOCOL

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

RELATED PRODUCTS

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: 9/5/2012 CMH

This page is intentionally left blank



www.GBiosciences.com