



243PR-02

**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™ KLH

Keyhole Limpet Hemocyanin  
Immunological Grade

(Cat. # 786-091)



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

## INTRODUCTION

KLH is a high purity grade keyhole limpet hemocyanin (KLH) prepared for immunotechnology research applications such as a carrier protein for antibody production. KLH is isolated from the mollusc *Megathura crenulata* and belongs to a family of giant respiratory proteins, the hemocyanins, which are found in molluscs and arthropods. The large molecular weight of KLH (400-3,000kD) makes it an excellent choice for an immunogen. The peptides are coupled to carrier proteins with the use of protein cross-linkers, normally through primary amine residues. KLH has a total of 300-600 lysine amines groups available for coupling, making KLH an outstanding carrier protein. KLH is supplied as a solution in convenient *OneQuant*<sup>™</sup> format of 2mg/vial.

## ITEM(S) SUPPLIED (Cat. # 786-091)

Description	Size
OneQuant <sup>™</sup> KLH (Immunological grade)	8 x 2mg/vial

## STORAGE CONDITIONS

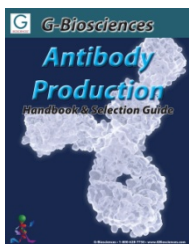
It is shipped at ambient temperature. Upon receiving, store it at 4°C.

## PROTOCOL

The protein concentration of the KLH solution is located on the product label.

## 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](http://www.GBiosciences.com) or contact us.

Last saved: 1/24/2014 CMH

*This page is intentionally left blank*



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