



Certificate of Analysis / Data Sheet

Product: KLH (Keyhole Limpet Hemocyanin)
Catalog No: 786-088, 786-091, 786-069, & 786-073
Release Date: 12/11/19
Retest Date: 12/11/23
Size: 2mg & 10mg
Storage: 4°C
Lot No: 191112

Specifications

CAS Number: 9013-73-3

Appearance: Faint blue to blue liquid

NOTE: Solution in phosphate buffered saline, pH 7.2, containing 1mM CaCl₂ and 0.5mM MgCl₂ with 15mM sodium azide as preservative

Test	Specification	Result
Appearance (Turbidity)	Clear to slightly hazy	Slightly Hazy
Copper ((w/w) <i>Cu as a percentage of protein</i>)	0.20 – 0.30%	0.21%
mg protein/ml (UV)	3.0 – 7.0	5.2

Based in E1% at 280nm = 15.7

This document has been produced electronically and is valid without signature

Date of Print: December 24, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

