

E.coli HCP ELISA Detection Kit

Overview

Cat.No. HG-HCP002

This kit is designed for the quantitative detection of HCP (host cell protein) content in biopharmaceuticals expressed on E.coli by using a double-antibody sandwich method.

This kit can be used to detect all components of HCP (host cell protein) in E.coli.

Specification

Assay range: 10-810ng/mL

Limit of quantitation: 10ng/mL

Limit of detection: 3.3ng/mL

Precision: CV% \leq 10%, RE% \leq ±15%

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

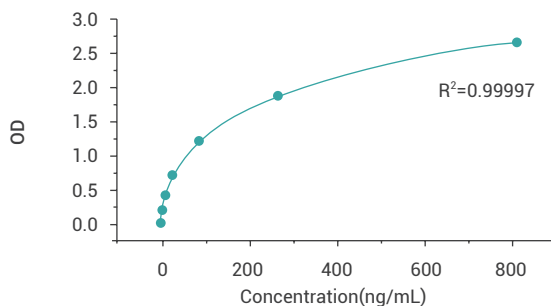
20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com info@eaglebio.com



Standard curve



Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
810	2.659	2.658	2.659
270	1.902	1.908	1.905
90	1.213	1.194	1.204
30	0.720	0.723	0.722
10	0.408	0.41	0.409
3.3	0.228	0.231	0.230
0	0.047	0.050	0.049