## E. coli host cell protein standards

Quantitative protein standards for Western blotting internal controls or ELISA standards.

*E. coli* host cell protein standards are native proteins that can be used as an internal control for the Western blot analysis of CHO Host Cell Proteins or as standards in an ELISA.

## **Ordering Information**

Order Online »

ENZ-PRT121-0025

25µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

Appearance Clear liquid.

Application Notes Protein standards for the E. coli host cell protein ELISA kit

or for Western blot analysis.

**Formulation** Liquid. In PBS containing 10 ppm Kathon.

**Source** Purified from *E. coli* host cells.

Last modified: May 29, 2024



eu@enzolifesciences.com