HEK293T host cell protein ELISA kit

Sensitive ELISA for the quantitative determination of host cell protein contamination in bulk products expressed in HEK293T expression systems.

The HEK293T host cell protein ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of HEK293T host cell protein contamination with results in 6 hours

Ordering Information

Order Online »

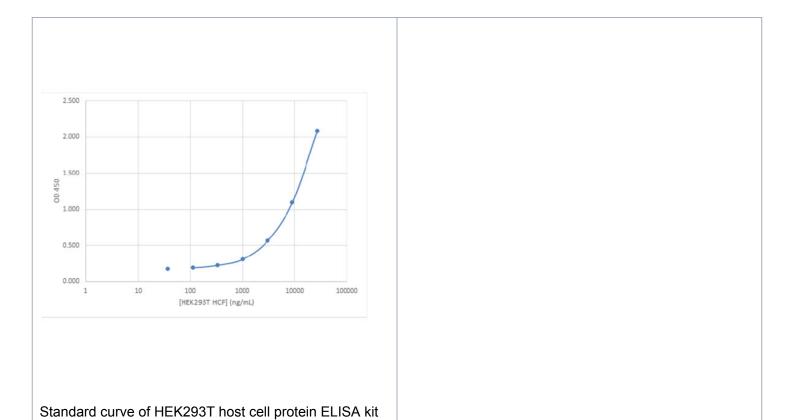
ENZ-KIT162-0001

96 wells

Manuals, SDS & CofA

View Online »

- Sensitive measurement of HEK293T host cell protein
- Low intra-assay variability allows for streamlined testing
- High throughput format with results in 6 hours
- Fully quantitative results that surpass semi-quantitative
 Western blot analysis



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HEK293T HCP ELISA kit

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of contamination with

HEK293T host cell protein.

Assay Time 6 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents 1 x 96 well plate, protein standards, 1x dilution buffer, 10x

PBS-T, reporting antibody, streptavidin-HRP conjugate,

TMB substrate, stop solution, plate sealer

Sensitivity 37 ng/ml (37-27000 ng/ml)

Species Reactivity Human

Last modified: May 29, 2024