

# His-Tag Protein ELISA Kit

Our His-Tag Protein ELISA Kit allows you to detect and quantify His-tagged protein samples simply and reliably by comparing your unknown samples to a known recombinant standard. Sensitivity range of the kit is 4 µg/mL to 1 ng/mL, or 950 nM to 50 pM of 6xHis-tag residues.

## Ordering Information

[Order Online »](#)

|             |          |
|-------------|----------|
| CBI-AKR-130 | 96 wells |
|-------------|----------|

## Manuals, SDS & CofA

[View Online »](#)

- Quantify proteins with His-tag at either N-terminus or C-terminus
- Detect as low as 1 ng/mL of protein or 50 pM 6xHis-tag residues
- Suitable for cell or tissue lysates

## Handling & Storage

Long Term Storage      +4°C

Shipping                      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application                      Colorimetric detection, ELISA

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)