

Octet[®] HIS2 Biosensors

For Quantitation of
HIS-tagged Proteins

Key Features

- Rapid detection of HIS-tagged proteins
- High-specificity capture of HIS-tagged proteins for easy quantitation
- Allows rapid analysis of crude or purified samples



Overview

The polyhistidine tag, commonly known as HIS-tag, is fused to recombinant proteins as a means of facilitating detection and purification. The Octet[®] Anti-HIS (HIS2) biosensor consists of high affinity, high specificity Anti-HIS antibody from Maine Biotechnology Services (MBS) pre-immobilized on a Sartorius fiber optic biosensor. In conjunction with the Octet[®] systems, the Octet[®] HIS2 biosensor provides a rapid and label-free method for HIS-tagged protein quantitation and kinetic analysis. The high specificity of the antibody-based biosensor enables the direct quantitation of HIS-tagged proteins in crude lysates, column eluents, cell lysates and cell culture supernatants, serving as an alternative to traditional time-consuming analytical methods such as HPLC and ELISA.

Flexibility and Versatility

The Octet® HIS2 Biosensor is qualified specifically for quantitation applications. It enables scientists to quickly and easily detect HIS-tagged recombinant proteins for quantitation measurements. Together with the Octet® N1 system's ease of use or the Octet® platform's throughput, Anti-HIS biosensors greatly accelerate laboratory workflows

and reduce time to results. The Octet® N1 system further enables measurement of precious samples with sample volume requirements as low as 4 µL.

For technical information on Anti-HIS biosensors, see Technical Note 37, Octet® Anti-HIS (HIS2) Biosensor Quantitation Assays.

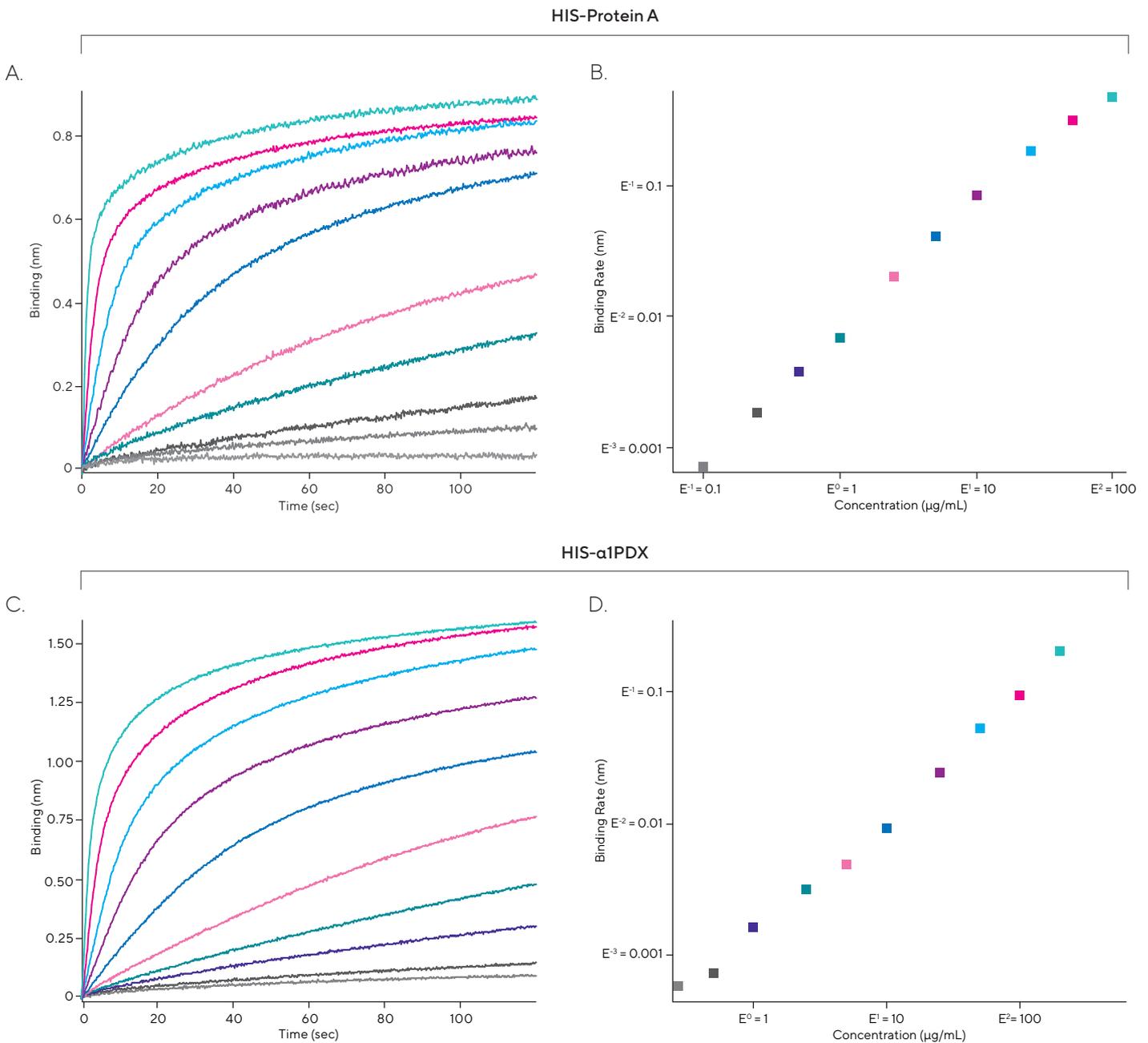


Figure 1: Detection of a HIS-Protein A (top) or HIS-α1PDX (bottom) standard using Anti-HIS biosensors on the Octet® RH16 system with assay parameters (1000 rpm, 2 minutes) for a standard dynamic range. (A) HIS-Protein A dose response. (B) HIS-Protein A calibration curve. (C) HIS-α1PDX dose response. (D) HIS-α1PDX calibration curve. Sample diluent was used as a matrix for all samples.

Range of Applications

The Anti-HIS biosensor offers researchers unparalleled ease of use and time to result in a wide range of laboratory applications such as:

- Rapid quantitation of any HIS-tagged proteins
- Easy protein expression monitoring
- Easy cell line development/optimization

Ordering Information

Part No.	UOM	Description
18-5114	Tray	One tray of 96 biosensors coated with the high-affinity, high specificity anti-HIS antibody for quantitative applications of all HIS-tagged proteins.
18-5115	Pack	Five trays of 96 biosensors coated with the high-affinity, high specificity anti-HIS antibody for quantitative applications of all HIS-tagged proteins.
18-5116	Case	Twenty trays of 96 biosensors coated with the high-affinity, high specificity anti-HIS antibody for quantitative applications of all HIS-tagged proteins.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 888 OCTET 75
Or +1 650 322 1360



For further contacts, visit
www.sartorius.com/octet-support