

Anti-HIS (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 Dip and Read™ Anti-HIS (HIS2) Biosensor consists of high affinity, high specificity Anti-HIS antibody from Maine Biotechnology Services (MBS) pre-immobilized on a ForteBio fiber optic biosensor. In conjunction with the Octet® and BLItz® systems, the Anti-HIS (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 Anti-HIS (HIS2) Dip and Read 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 BLItz system's ease of use or the Octet platform's throughput, Anti-HIS biosensors greatly accelerate laboratory workflows and reduce time to results. The BLItz 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 (Anti-HIS Biosensor Quantitation Assays).

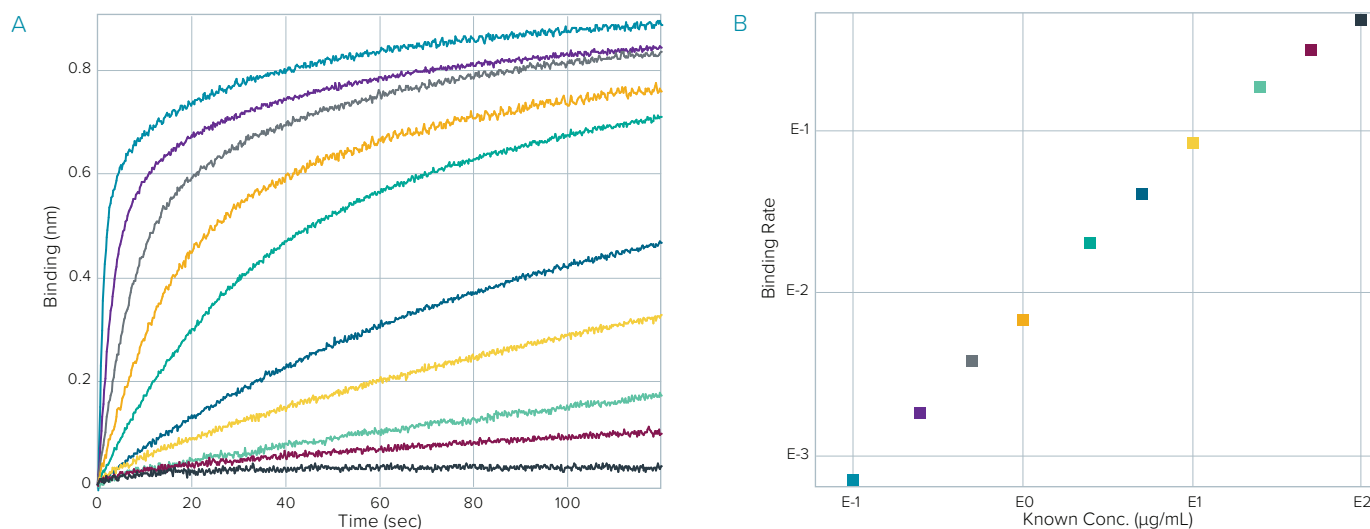
Range of applications

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



HIS-Protein A



HIS-α1PDX

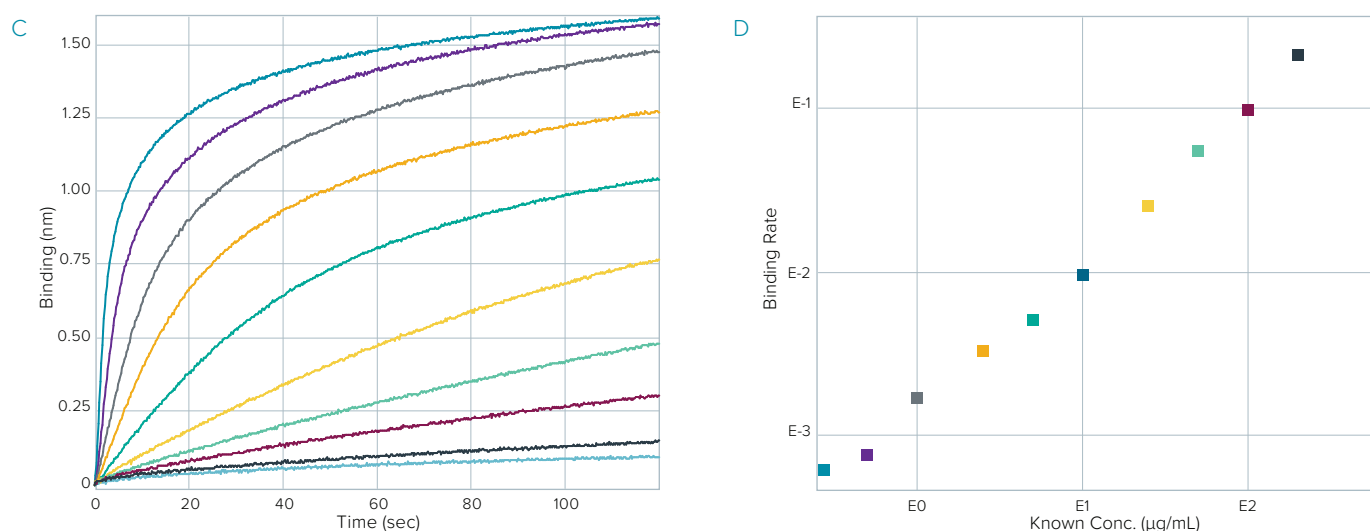


Figure 1: Detection of a HIS-Protein A (top) or HIS-α1PDX (bottom) standard using Anti-HIS biosensors on the Octet RED384 system with assay parameters (1000 rpm, 2 minutes) for a standard dynamic range. A) HIS-Protein A dose response. B) HIS-Protein A calibration curve. C) HIS-α1PDX dose response. D) HIS-α1PDX calibration curve. Sample diluent was used as a matrix for all samples.

Ordering information

Part No.	UOM	Description
18-5114	Tray	One tray of 96 Anti-HIS (HIS2) Biosensors
18-5115	Pack	Five trays of 96 Anti-HIS (HIS2) Biosensors
18-5116	Case	Twenty trays of 96 Anti-HIS (HIS2) Biosensors



FortéBio
 47661 Fremont Boulevard
 Fremont, CA 94538
 888.OCTET-75 or 650.322.1360
 fortebio.info@moldev.com

FortéBio Analytics (Shanghai) Co., Ltd.
 No. 88 Shang Ke Road
 Zhangjiang Hi-tech Park
 Shanghai, China 201210
 salesops.china@moldev.com

Molecular Devices (UK) Ltd.
 660-665 Eskdale
 Winnersh Triangle
 Wokingham, Berkshire
 RG41 5TS, United Kingdom
 +44 118 944 8000
 uk@moldev.com

Molecular Devices (Germany) GmbH
 Bismarckring 39
 88400 Biberach an der Riss
 Germany
 +00800 665 32860