



Fortebio Analytics (Shanghai) Co., Ltd.
No.88 Shang Ke Road Zhangjiang Hi-tech Park
Shanghai, 201210, P.R. China
fortebiosupport@moldev.com , www.fortebio.com/support

Certificate of Analysis

Streptavidin (SA) Biosensors

Sales Numbers: 18-5019 to 18-5021

Part Number: 18-0009

Lot Number: 1908082

Date of Manufacturing: Aug-2019

Test Date: Aug-2019

Test Instrument: Octet RED384

Test Temperature: 30°C

Expiration Date: Aug-2020

TEST

SPECIFICATION

RESULT

Baseline Drift (nm) over 1 hour, 1000rpm

-0.025 – 0.060 (Lot Mean)
≤ 0.045 (Lot Std Dev)

0.054 (n=46)
0.008 (n=46)

Biotin-Protein A Binding (nm)

≥ 1.35 (Lot Mean)
≤ 12% (Lot %CV)

1.88 (n=182)
3.6% (n=182)

Verified By: 

Date: 8/15 2019

Approved By: 

Date: 8/15 2019