

# 384-Well Tilted-Bottom Microplate

*Reduced sample volumes and improved sensitivity for Octet assays*

## Key features

- Assay with volumes as low as 40  $\mu\text{L}$  per well
- Reduces variation in baseline signals, and improves assay sensitivity
- Use with Octet® HTX, RED384 and QK384 systems
- 384-well, black, polypropylene microplate
- Standard SBS format

The 384TW microplate is a black, polypropylene, 384-well tilted-bottom plate designed for ForteBio's Octet family of label-free biomolecular interaction analysis instruments. It can be used on the Octet QK384, RED384 and HTX instruments, all of which accommodate the 384-well microplate format. The 384TW microplate enables use of samples as small as 40  $\mu\text{L}$  in Octet assays and reduces variability in the background signal, thus improving assay sensitivity, particularly beneficial for peptide, small molecule and fragment analysis.

Test concentrations	Measured avg. concentrations ( $\mu\text{g}/\text{mL}$ )		Percent CV	
	96-well	384TW	96-well	384TW
2000	2000.00	1999.98	12%	7%
1500	1508.78	1503.58	7%	3%
1000	1015.90	1007.08	7%	4%
700	700.20	700.70	2%	3%
500	503.03	502.53	3%	4%
300	300.50	300.88	2%	3%
30	30.10	30.05	3%	3%
10	10.01	10.00	3%	3%
3	3.00	3.00	3%	3%
1	1.00	1.00	1%	6%
0.5	0.50	0.51	4%	9%

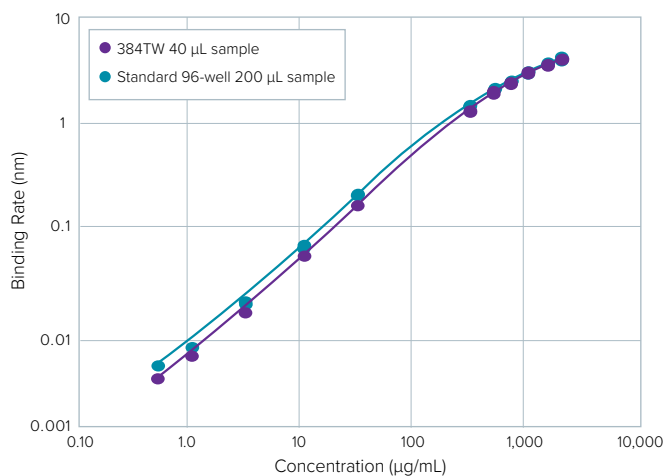
**Table 1:** Measured concentrations and CVs show good dynamic range, recovery and precision using the 384TW vs. 96-well standard microplate.

## Assays using limited sample volume

When assaying precious samples, the ability to limit the volume of sample per experiment is critical. The 384TW microplate provides the ability to perform assays using as little as 40  $\mu\text{L}$  per well. This is a 3-fold decrease over the volume needed in a standard 384-well plate and a 5-fold decrease over the sample volume needed in a 96-well plate. The assay performance (dynamic range, precision, accuracy) for kinetic characterization and quantitation in the 384TW microplate are equivalent to that in a black, polypropylene 384-well standard flat-bottom microplate.

Figure 1 shows results from a Protein A biosensor-based quantitation assay to detect human IgG (hIgG) in solution. The assay was run on Octet RED384 system using the standard Basic Quantitation settings with Regeneration module, available in the Octet Data Acquisition software.

Table 1 compares data obtained in a 96-well flat-bottom microplate (Figure 1) using 200  $\mu\text{L}$  of sample to that obtained in a 384TW microplate using only 40  $\mu\text{L}$  of sample shows that assay performance is equivalent in both microplates.



**Figure 1:** Calibration curve across the dynamic range of 0.5–2000  $\mu\text{g}/\text{mL}$  hIgG; results from 384TW microplates compares very well to those from standard flat-bottom 96-well microplates.

	384-well flat-bottom plate	384TW plate
Baseline signal average (PBS)	-4.8 pm	-2.0 pm
Baseline signal standard deviation	10.4 pm	4.3 pm
Lower limit of signal detection	26.4 pm	10.9 pm

**Table 2:** 384TW data for Super Streptavidin biosensors on the Octet RED384 system. The data shows a more than 2-fold improvement in detection limit for the 384TW microplate.

## Improved sensitivity in small molecule and peptide analyses

In high-sensitivity screening and kinetic applications, the limit of detection is often defined as 3 times the standard deviation of the baseline signal. The 384TW microplate decreases system artifacts when used on the Octet QK384, RED384 and HTX systems, decreasing the variation (and thus the standard deviation) of the baseline signal, allowing smaller signals to be resolved from baseline noise.

Table 2 contains data obtained on the Octet RED384 system using Super Streptavidin (SSA) biosensors. It includes the average signal and standard deviation of the baseline of these biosensors. The averages and standard deviations for each data point are calculated from 1920 replicates. The lower limit of detection shown is defined as 3 times the standard deviation over the average baseline. For many screening campaigns, this value defines the smallest signal contribution by a positive binding sample. In this experiment the 384TW microplate was found to be advantageous, having a 2-fold lower limit of detection than a standard 384-well flat-bottom plate.

## Instructions for use

The 384TW microplate is used on the Octet HTX, RED384 and QK384 systems in the same way as a flat-bottom 384-well microplate with the following exceptions:

- Minimum sample volume is 40 µL.
- Maximum sample volume is 100 µL.
- When pipetting into the wells, avoid trapping bubbles of air in them. Bubbles in the well can result in spikes in the assay data. An effective method for removing any bubbles that occur is to centrifuge the filled plate in a swinging plate rotor at 1000 rpm for 2 minutes prior to the assay.
- Assays should be set up with the biosensor offset at 3 mm.

- Set the “Delay experiment start” setting available in the Octet Data Acquisition software to a minimum of 10 minutes (600 seconds) to allow the plate to equilibrate inside the instrument prior to the first data point.
- Check the “Shake sample plate while waiting” option in the Octet Data Acquisition software. Shaking the plate during the equilibration time helps to eliminate trapped air.

## 384TW Microplate specifications\*

Specification	Details
Applications	Low volume and high sensitivity applications
Instrument compatibility	Octet HTX, RED384 and QK384 systems
Material	Polypropylene
Color	Black
Dimensions	SBS standard
Run time on Octet RED384, QK384 and HTX systems	Screening applications: <10 hours at 40 µL (biosensor offset 3 mm) Kinetics & quantitation applications: <4 hours at 40 µL (biosensor offset 3 mm )
Evaporation rate	Non-linear rate of 30% loss at 6 hours for initial volume of 40 µL of PBS
Maximum volume	100 µL
Minimum volume	40 µL
Sterilization	No

## Ordering information

Part No.	UOM	Description
18-5080	Pack	384-well microplate, black, tilted-bottom, polypropylene, 10 per pack. Enables low volume samples, lowers background for peptide, small molecule and fragment analysis applications.
18-5076	Case	384-well microplate, black, tilted-bottom, polypropylene, 100 per case. Enables low volume samples, lowers background for peptide, small molecule and fragment analysis applications.

\* All specifications are subject to change without notice.



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

**ForteBio 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