

Turbidometric Growth Curve Measurements and Liquid Evaporation Studies with a Microplate Photometer

Reija-Riitta Harinen, Jorma Lampinen, Marika Raitio and Vuokko Kytöniemi, Thermo Fisher Scientific, Vantaa, Finland



Key words

- Thermo Scientific Multiskan FC
- Thermo Scientific SkanIt Software
- Photometer
- Turbidometry
- Antibiotic Susceptibility

Abstract

This paper describes how a filter-based microplate photometer, Thermo Scientific Multiskan FC, can be used reliably and easily to measure turbidity of bacterial samples in antibiotic susceptibility studies. The growth curve measurements were performed on a 96-well microplate. The performance of turbidometric measurements on a 96-well microplate and the liquid evaporation from 96- and 384-well microplates during long incubation were also studied.

Introduction

The turbidity of a liquid sample is traditionally measured either with a nephelometer or a photometer. The measure of turbidity (also called optical density, OD) refers to the scattering of light from the particles of the measured sample. When measuring turbidity of a bacterial culture, the amount of scattered light is directly proportional to the number of bacterial cells in the sample.

Bacterial growth curves are usually measured by using a photometer to detect the optical density of a sample. Photometers are designed to measure light absorbance by a sample, but they can also be utilized to study sample turbidity, if a nephelometer is not available. Light absorbance is a completely different kind of phenomenon

than light scattering from sample molecules. Therefore photometers are not optimal for turbidometric assays, but are still suitable for that purpose. Nephelometers are more sensitive in turbidometric measurements, but cannot be used for measuring absorbance.

Using microplates to cultivate bacteria facilitates the simultaneous measurement of a high number of samples. Therefore different growth conditions can be easily tested on a microplate format. The possibility to create a kinetic measurement protocol with specific interval times for kinetic readings, shaking the plate during the intervals and keeping the instrument in elevated temperature during a long kinetic measurement can all be achieved with a good microplate photometer.

Materials and Methods

The filter-based microplate photometer, Multiskan® FC, was used in three kinds of studies:

1. Bacterial growth curves

Salmonella Typhimurium was grown in the presence of varying concentrations of antibiotic (erythromycin) on a 96-well plate for 16 hours. The Multiskan FC was programmed by the Thermo Scientific SkanIt Software (PC software) to handle the entire kinetic protocol so the plate was automatically measured every 15 minutes and shaken during the kinetic interval times. The temperature of the instrument was kept at 37°C throughout the protocol execution.

2. Turbidometric performance

Salmonella Typhimurium was grown to late logarithmic phase (OD = 3.19). Then a dilution series of ½ dilutions was made from the

culture and 300 µl of each dilution was added to the wells of a 96-well microplate with 7 replicates. The OD of the dilution series was measured at 595 nm with Multiskan FC, and with a cuvette spectrophotometer for reference.

3. Evaporation studies

The idea was to study how much the liquid in the wells evaporates when the plate is incubated for hours inside the instrument in elevated temperature. The 384-well plate was filled with water (50 µl / well) and the plate was incubated at 37°C inside the Multiskan FC, both with and without a lid, for 24 hours. Evaporation from a 96-well plate (300 µl water / well) was followed by incubating the plate at 37 and 50°C without a lid. The liquid evaporation was followed by weighing the plate, and by measuring the changes in water absorbance at 975 nm at the time points of 0, 1, 3, 7 and 24 hours. The absorbance was measured with the Thermo Scientific Varioskan Flash, a multimode microplate reader.

The 96-well plates used in these measurements were Thermo Scientific 96 Well Immulon Microtiter Solid Microplates Medium/1B (# 3355), and the 384-well plates were Microtiter plates (# 95040000, round well, height 10 mm). The maximum height of a plate with a lid can be 15.3 mm with the Multiskan FC.

Result

1. Bacterial growth curves

The bacterial growth curves of Salmonella Typhimurium with serial dilutions of erythromycin are presented in figure 1. The differences in the growth speed caused by antibiotic inhibition

could be clearly and easily detected by the OD measurement. The SkanIt® Software controlling the Multiskan FC, automatically created the kinetic curves during the measurement. The software's built-in effective dose (ED) calculation can also be utilized to set the software up to calculate the antibiotic concentration causing the 50% percent (or other user-defined percent) growth inhibition. The ED calculation tool is practical especially in cytotoxicity assays.

linearity of the Multiskan FC goes up to 3.0 Abs. The results of turbidometric measurement do not directly correlate to linearity performance of absorbance due to different nature of absorbance and light scattering. The OD of 1/10 dilution of the culture was measured with a cuvette spectrophotometer to be 0.319. Based on this measurement, the theoretical OD values for the whole dilution series were calculated. The theoretical values

different from the 1 cm pathlength of a standard cuvette. Therefore the microplate measurement values were corrected to correlate to 1 cm pathlength. The calculated differences between theoretical and measured values show that OD values measured with Multiskan FC correlate better to the theoretical values in wider part of the dilution series than the cuvette spectrophotometer.

3. Evaporation studies

The water weight loss from a 96- and 384-well plate due to evaporation during incubation is presented in table 2. The results are presented as water weight loss in grams and as a percentage of the initial weight. A long incubation time clearly causes the liquid to evaporate when the plate is not covered with a lid. But the amount of evaporation without a lid is still not significant for example at a 7 hour time point either with 96- or 384-well plates. The lid obviously inhibits evaporation efficiently, therefore 24 hours at 37°C causes only < 9 % of the plate liquid to evaporate on a 384 well plate.

The changes in water absorbance due to evaporation on a 96-well plate are presented in figure 2. It is clearly seen that evaporation increases towards the edges of the plate. Figure 3 presents the changes in water absorbance on a 384-well plate. The lid expectedly slows down evaporation very efficiently. Incubation at high temperature may cause water to condensate

on the lid and disturb the measurement through the lid. The condensation would have been seen as random increase of absorbance, which clearly was not seen in these results. Results show the incubator in the Multiskan FC performs well.

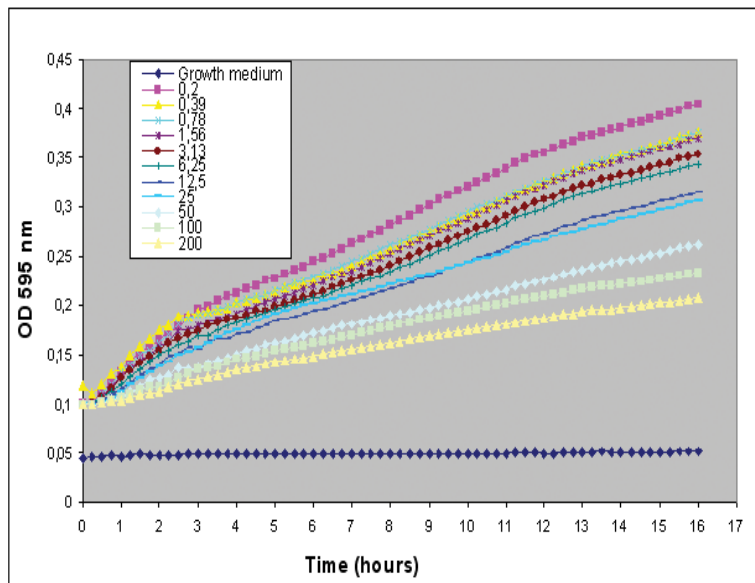


Figure 1. The growth curves of Salmonella Typhimurium at varying concentrations (0.2 – 200 µg/ml) of erythromycin.

2. Turbidometric performance

Table 1 presents the results of OD measurements of a bacterial dilution series. Multiskan FC was compared to a cuvette spectrophotometer of another manufacturer. The photometric

are here compared to the actual measurement values given by the cuvette spectrophotometer (in yellow) and a microplate photometer, Multiskan FC (in green). The sample pathlength on microplate wells is naturally

Dilution	Theoretical OD	OD (cuvette spectrophotometer)	Difference (%) to theoretical OD	OD (Multiskan FC)	OD changed to correlate to 1 cm pathlength	Difference (%) to theoretical OD
1	3.190	1.941	-64.3	1.287	2.253	-41.6
2	1.595	1.291	-23.5	0.761	1.331	-19.8
3	0.798	0.894	10.8	0.421	0.737	-8.2
4	0.399	0.514	22.4	0.229	0.400	0.4
5	0.199	0.273	27.0	0.113	0.199	-0.4
6	0.0997	0.138	27.8	0.059	0.103	3.0
7	0.0498	0.07	28.8	0.029	0.051	1.7
8	0.0249	0.035	28.8	0.015	0.026	2.4
9	0.0125	0.021	40.7	0.007	0.013	2.6
10	0.0062	0.007	11.0	0.004	0.008	17.6
11	0.0031	-	-	0.003	0.005	41.4
12	0.0016	0.003	48.1	0.002	0.003	53.3

Table 1. The optical density (OD) measurements of the bacterial dilution series.

Time (hours)	Water (g)	Water weight loss (g) at 37°C	Water weight loss (%) at 37°C
0	29.1	-	-
1	28.9	0.3	1.0
3	28.4	0.8	2.7
7	27.5	1.6	5.6
24	24.3	4.8	16.6
Time (hours)	Water (g)	Water weight loss (g) at 50°C	Water weight loss (%) at 50°C
0	29.2	-	-
1	28.6	0.5	1.8
3	27.5	1.6	5.6
7	25.6	3.5	12.0
24	18.9	10.3	35.3

Water (g)	Water weight loss (g) without a lid	Water weight loss (%) without a lid
19.4	-	-
18.9	0.5	2.5
18.1	1.4	7.1
16.6	2.9	15.9
10.8	8.6	51.9
Water (g)	Water weight loss (g) with a lid	Water weight loss (%) with a lid
19.4	-	-
19.3	0.1	0.5
19.1	0.3	1.6
18.8	0.6	3.2
17.8	1.6	8.6

Table 2. Water evaporation during 24 hour incubation from a 96-well plate (in yellow) at 37° and 50° C without a lid, and from a 384-well plate (in blue) at 37° C with and without a lid.

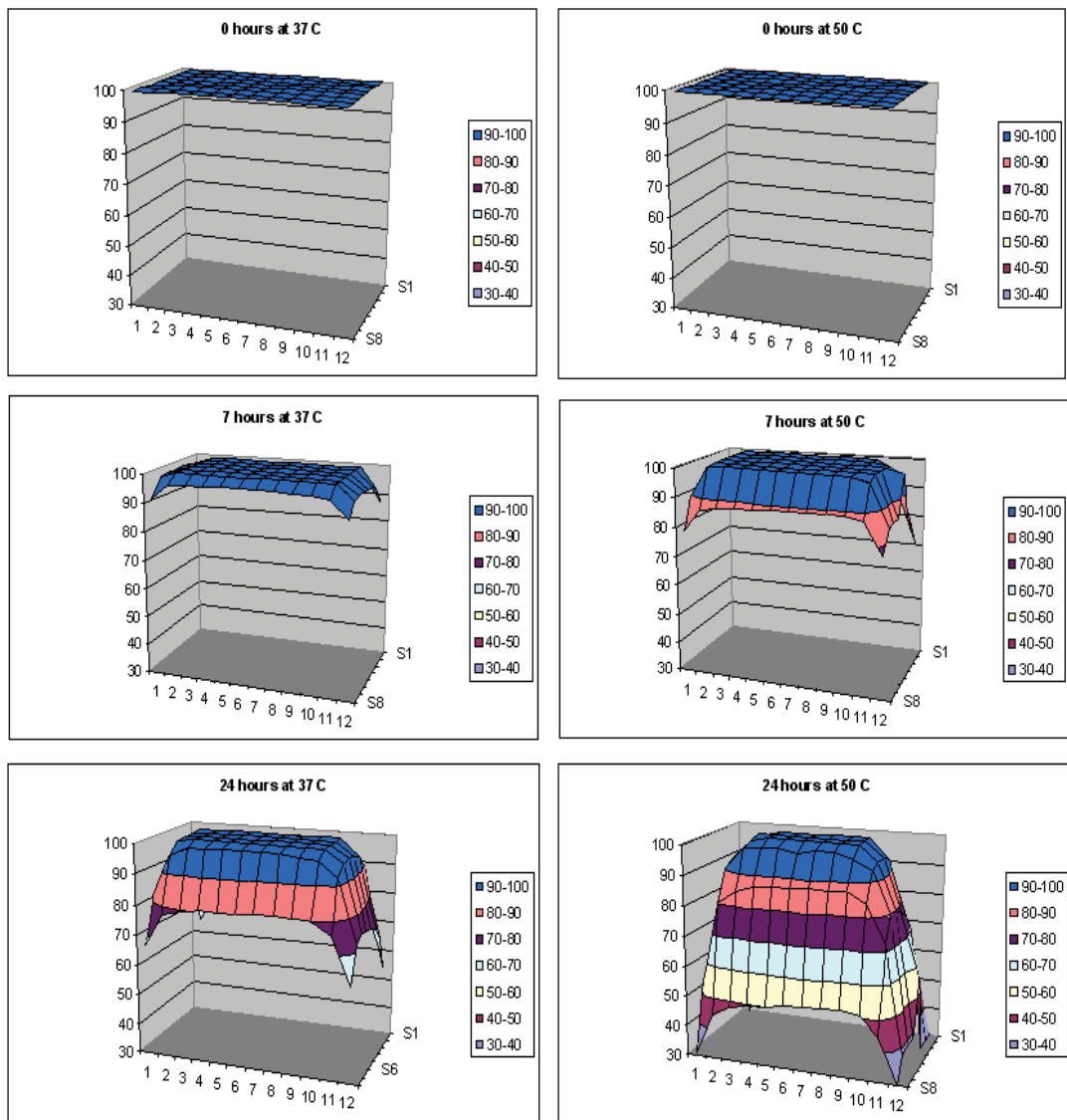


Figure 2. Evaporation on different parts of a 96-well plate during incubation at 37° and 50° C without a lid. The effects of evaporation on the wells are presented as a plate shaped graph (y-axis = % of absorbance at 975 nm, x-axis = columns of the plate, z-axis = rows of the plate) to indicate on which parts of the plate the evaporation is the fastest (the absorbance decreases the most).

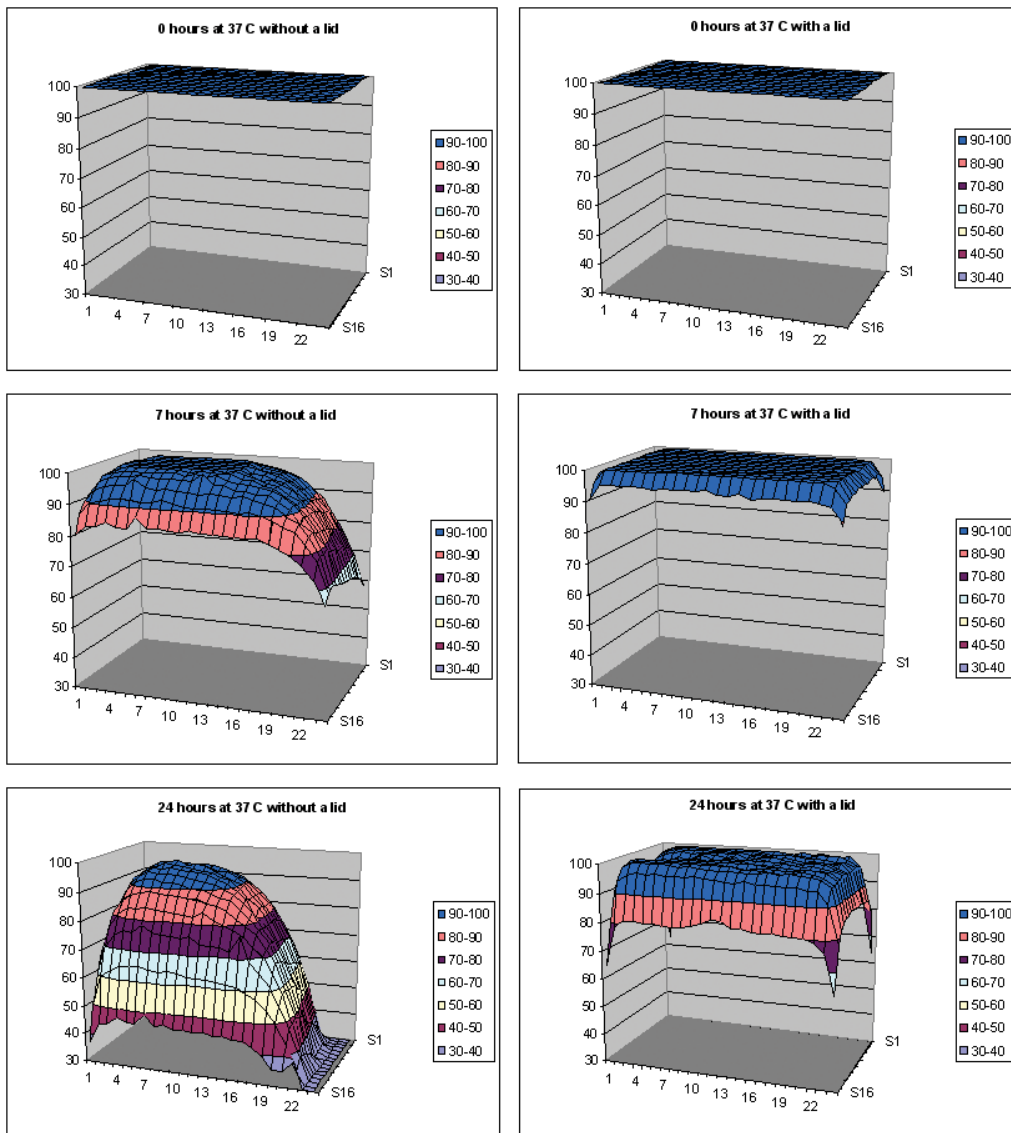


Figure 3. Evaporation on different parts of a 384-well plate during incubation at 37° C with and without a lid. (y-axis = % of absorbance at 975 nm, x-axis = columns of the plate, z-axis = rows of the plate).

Conclusion

The Multiskan FC is well suited for turbidometric bacterial measurements. The OD of the samples in kinetic measurements can be reliably monitored at a wide OD range. The liquid evaporation from microplate wells in elevated temperature must be taken into account, when the incubation time

is long. The effects of evaporation can be minimized e.g. by adding the samples to the plate by leaving the edge wells empty. Filling the empty edge wells with water also decreases evaporation from the actual sample wells, as the evaporation occurs primarily from the edge wells.

Further information

For further information about the Multiskan FC microplate photometer, and SkanIt Software please refer to the following web page:
www.thermoscientific.com/readingroom
www.thermoscientific.com/mpi

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

North America:
USA / Canada
 1 800 345 0206

Europe:
Austria
 +43 1 801 40 0

Belgium
 +32 53 73 42 41

France
 +33 2 2803 2180

Germany national toll free
 08001-536 376

Germany international
 +49 6184 90 6940

Italy
 +39 02 02 95059 448

Netherlands
 +31 76 571 4440

Nordic countries
 +358 9 329 100

Russia/CIS
 +7 (495) 739 76 41

Spain/Portugal
 +34 93 223 09 18

Switzerland
 +41 44 454 12 12

UK/Ireland
 +44 870 609 9203

Asia:
China

+86 21 6865 4588 or
 +86 10 8419 3588

India toll free
 1800 22 8374

India
 +91 22 6716 2200

Japan
 +81 45 453 9220

Other Asian countries
 +852 2885 4613

Countries not listed:
 +49 6184 90 6940 or
 +33 2 2803 2180

www.thermoscientific.com/readingroom
www.thermoscientific.com/mpi

© 2010 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.