

# A Multi-Site Study Comparing a Commercially Prepared Dried MIC Susceptibility System to the CLSI/ISO Broth Microdilution Method for Ceftibuten-Clavulanic Acid using Gram-Negative Non-Fastidious Organisms

\*T.C. Lewis<sup>1</sup>, D.T. Staats<sup>1</sup>, N. M. Holliday<sup>1</sup>, C. C. Knapp<sup>1</sup>, S.B. Killian<sup>1</sup>, B.J. Olson<sup>2</sup>, C.L. Pike<sup>2</sup>, E.K. Donnerbauer<sup>2</sup>, N.P. Mehta<sup>2</sup>, T.R. Fritsche<sup>2</sup>, A. Gattis<sup>3</sup>, N. Waugh<sup>3</sup>, K.Doing<sup>3</sup>, J.Vlooswijk<sup>4</sup>, A. Cornelisse<sup>4</sup>, E. Scopes<sup>5</sup>, J. Williams<sup>5</sup>, A-M. Leonte<sup>5</sup>, K. Krause<sup>6</sup>

<sup>1</sup>Thermo Fisher Scientific, Cleveland, OH; <sup>2</sup> Marshfield Clinic, Marshfield, WI; <sup>3</sup>Centura Health, Colorado Springs, CO; <sup>4</sup>University Medical Center Utrecht, Utrecht, Netherlands; <sup>5</sup>Thermo Fisher Scientific, Basingstoke, UK; <sup>6</sup>Achaogen, Inc. South San Francisco, CA

## ABSTRACT

**Background:** Ceftibuten-clavulanic acid (T/CA) (Achaogen, Inc. South San Francisco, CA) is a combination of a 3<sup>rd</sup> generation cephalosporin (ceftibuten) and a beta-lactamase inhibitor (clavulanic acid) displaying activity against highly resistant gram-negative non-fastidious organisms, including Extended Spectrum Beta Lactamase (ESBL) producing strains of *Enterobacteriaceae*. A 4-site evaluation was designed to determine the accuracy and reproducibility of T/CA susceptibility testing against non-fastidious gram negative organisms using the Thermo Scientific™ Sensititre™ dried MIC susceptibility system compared with the CLSI (M07, M100)/ ISO 20776-1, ISO 20776-2 (CLSI/ISO) reference broth microdilution method (BMD). **Materials and Methods:** The Sensititre 18-24 Hour MIC or Breakpoint Susceptibility System with T/CA in the dilution range of 0.03/0.015-32/16µg/ml was tested against 455 recent clinical, challenge, and reproducibility *Enterobacteriaceae* isolates. The Sensititre dried MIC susceptibility system was inoculated per manufacturers' instructions. BMD was performed per CLSI/ISO guidelines. Recommended CLSI quality control (QC) organisms were tested daily and all results were within the published QC ranges. **Results:** Comparisons of the indicated gram-negative non-fastidious organisms MIC results on the Sensititre system to the CLSI/ISO BMD for automated and manual reads resulted in 96.6% and 97.5% essential agreement (EA; +/- 1 log<sub>2</sub> dilution) for T/CA, respectively. Overall agreement for the reproducibility (+/- 1 log<sub>2</sub> dilution of the modal MIC) using automated and manual reads was 99.7% and 99.0% respectively. **Conclusion:** The Sensititre susceptibility system demonstrates an equivalent level of performance compared to the CLSI/ISO BMD method when testing T/CA against gram negative non-fastidious organisms. The high level of agreement obtained by the Sensititre susceptibility system and the CLSI/ISO BMD method suggests that this is an acceptable method for susceptibility testing of T/CA.

## INTRODUCTION

Ceftibuten-clavulanic acid is combination of ceftibuten, an approved third generation cephalosporin, and clauanate, an approved beta-lactamase inhibitor. This agent, T/CA, has demonstrated *in vitro* activity and *in vivo* efficacy against complicated urinary tract infections (cUTI) caused by *Escherichia coli* and *Klebsiella pneumoniae* that produce extended spectrum beta lactamase (ESBL). This *in vitro* comparison study was done to validate the performance of ceftibuten-clavulanic acid on the commercially manufactured Thermo Scientific™ Sensititre™ 18 – 24 hour Dried Susceptibility Plate with the standard reference broth microdilution method recommended by the Clinical and Laboratory Standards Institute (CLSI M07/ M100) and ISO (20776-1). To establish equivalency for both auto and manual read methodologies, a series of studies were conducted at 4 trial sites including testing of clinical/challenge, reproducibility and quality control isolates.

## MATERIALS AND METHODS

●The Sensititre 18-24 hour MIC or breakpoint susceptibility system (Thermo Fisher Scientific, Oakwood Village, OH) is an *in vitro* diagnostic product for clinical susceptibility testing of both fastidious and non-fastidious organisms. ceftibuten-clavulanic acid was tested against: (Table 1.)

- 367 recent clinical gram-negative isolates across the four sites
- 77 challenge isolates at a single testing site
- 11 reproducibility isolates at each site (tested in triplicate over a 3 day testing period)
- 4 Quality Control Strains (ATCC) (Table 2.)



## MATERIALS AND METHODS Cont.

● Colony Counts and purity plates were performed on the inoculums of the Clinical, Challenge, Reproducibility and QC strains on each day of testing.

● Each isolate was tested using a:

- Dried Sensititre 18–24 susceptibility plate containing ceftibuten-clavulanic acid (0.03/0.015-32/16µg/ml). The dried plates were set up and tested by both automated and manual reading methodologies according to the manufacturer's instructions.

- Reference broth microdilution plate was prepared and tested on each isolate according to the current Clinical Laboratory Standards Institute and ISO standard method.

Table 1. Organisms Tested	Number Tested
<b>Clinical Isolates</b> (4 sites)	367
<b>Challenge Isolates</b> (one site)	77
<b>Reproducibility Isolates</b> (4 sites) (3 x day for 3 days)	11 (396)
<b>ATCC Quality Control Strains</b> (at least 20 replicates of each strain at 3 sites, 1 site yielded fewer results)	4 (323)
<b>TOTAL</b>	<b>1163</b>

### Quality Control

● Recommended CLSI quality control (QC) organisms were tested daily and were within the CLSI expected QC ranges.

● Colony counts were performed and fell within expected ranges  
Reference 2-8X10<sup>5</sup> CFU, Sensititre 5X10<sup>4</sup>-5X10<sup>5</sup> CFU

Table 2. Quality Control Strains	Expected CLSI QC Ranges (µg/ml)
<i>Escherichia coli</i> ATCC 25922	0.12/0.06-0.5/0.25
<i>Escherichia coli</i> NCTC 13353	0.25/0.12-1/0.5
<i>Pseudomonas aeruginosa</i> ATCC 27853	>32/16
<i>Klebsiella pneumoniae</i> ATCC 700603	0.06/0.03-0.25/0.12

## Results

Essential agreement for ceftibuten-clavulanic acid on the Sensititre susceptibility plate compared to the reference microdilution plate was calculated for each read method (Auto and Manual) using the ±1 log<sub>2</sub> dilution standard. Essential agreement rates are shown for gram-negative non-fastidious isolates in **Tables 3 and 4**.

### Clinical Isolates and Challenge Organisms

The overall essential agreement for ceftibuten-clavulanic acid within ±1 log<sub>2</sub> dilution was **96.6%** for the auto read method and **97.5%** for the manual method.

### Inter-laboratory Reproducibility

Reproducibility testing results for ceftibuten-clavulanic acid within ±1 log<sub>2</sub> dilution from the modal MIC was **99.7%** for the auto read method and **99.0%** for the manual method.

**Table 3. Summary Data and % Essential Agreement of Gram-Negative Non-Fastidious Clinical and Challenge Isolates Using the Manual Read Method**

The overall essential agreement for ceftibuten-clavulanic acid within +/- one log<sub>2</sub> dilution, was **97.5%** for the manual read method

### Combined Total Isolates

Ceftibuten-clavulanic acid	Number of Isolates		Essential Agreement		% Essential Agreement	
	All	Evaluable	Total	Evaluable	Total	Evaluable
<b>Organism Group</b>						
<i>Escherichia coli</i>	102	86	97	81	95.1%	94.2%
<i>Klebsiella pneumoniae</i>	94	59	90	56	95.7%	94.9%
<i>Klebsiella oxytoca</i>	43	9	43	9	100.0%	100.0%
<i>Enterobacter cloacae</i>	62	35	61	34	98.4%	97.1%
<i>Klebsiella aerogenes</i>	39	29	39	29	100.0%	100.0%
<i>Citrobacter freundii</i>	26	22	25	21	96.2%	95.5%
<i>Serratia marcescens</i>	26	24	26	24	100.0%	100.0%
<i>Proteus vulgaris</i>	23	1	23	1	100.0%	100.0%
<i>Proteus mirabilis</i>	29	3	29	3	100.0%	100.0%
<b>Total</b>	<b>444</b>	<b>268</b>	<b>433</b>	<b>258</b>	<b>97.5%</b>	<b>96.3%</b>

**Table 4. Summary Data and % Essential Agreement of Gram-Negative Non-Fastidious Clinical and Challenge Isolates Using the Auto Read Method**

The overall essential agreement for ceftibuten-clavulanic acid within +/- one log<sub>2</sub> dilution, was **96.6%** for the auto read method

### Combined Total Isolates

Ceftibuten-clavulanic acid	Number of Isolates		Essential Agreement		% Essential Agreement	
	All	Evaluable	Total	Evaluable	Total	Evaluable
<b>Organism Group</b>						
<i>Escherichia coli</i>	102	86	97	81	95.1%	94.2%
<i>Klebsiella pneumoniae</i>	94	59	89	55	94.7%	93.2%
<i>Klebsiella oxytoca</i>	43	8	43	8	100.0%	100.0%
<i>Enterobacter cloacae</i>	62	32	60	30	96.8%	93.8%
<i>Klebsiella aerogenes</i>	39	26	39	26	100.0%	100.0%
<i>Citrobacter freundii</i>	26	21	25	20	96.2%	95.2%
<i>Serratia marcescens</i>	26	23	25	23	96.2%	100.0%
<i>Proteus vulgaris</i>	23	1	22	1	95.7%	100.0%
<i>Proteus mirabilis</i>	29	3	29	3	100.0%	100.0%
<b>Total</b>	<b>444</b>	<b>259</b>	<b>429</b>	<b>247</b>	<b>96.6%</b>	<b>95.4%</b>



## RESULTS Cont.

**Table 5. Inter-laboratory Reproducibility % Essential Agreement ±1 log<sub>2</sub> dilution from the Modal Value**

Ceftibuten-clavulanic acid	Auto Read	Manual Read
Between-site total isolates tested	396	396
Between-site isolates within +/- 1 well from mode	395	392
Between-site reproducibility ratio	395	392
Between-site reproducibility %	<b>99.7%</b>	<b>99.0%</b>
Total essential agreement	395/396	392/396
<b>Essential agreement %</b>	<b>99.7%</b>	<b>99.0%</b>

## CONCLUSIONS

This study validates that the Sensititre 18–24 hour susceptibility system (both auto read and manual read) demonstrated an equivalent level of performance compared to the CLSI M07/M100 and ISO 20776-1 reference broth microdilution plate when testing ceftibuten-clavulanic acid against gram-negative non-fastidious clinical and challenge isolates. This study suggests that this is an acceptable method for susceptibility testing of ceftibuten-clavulanic acid.

## REFERENCES

Clinical and Laboratory Standards Institute. 2015. *Methods for dilution antimicrobial susceptibility tests for bacteria that grow aerobically; approved standard-tenth edition*. Approved document M07-A10. Wayne, PA: CLSI.

Clinical and Laboratory Standards Institute. 2018. *Performance Standards for Antimicrobial Susceptibility Testing; Twenty-eighth Informational Supplement M100-S28*. Wayne, PA: CLSI.

FDA Guidance for Industry and FDA Class II Special Controls Guidance Document: Antimicrobial Susceptibility Test (AST) Systems, August 28, 2009.

*Clinical laboratory testing and in vitro diagnostic test systems - Susceptibility testing of infectious agents and evaluation of performance of antimicrobial susceptibility test devices - Part 1: Reference method for testing the in vitro activity of antimicrobial agents against rapidly growing aerobic bacteria involved in infectious diseases* (ISO 20776-1 :2006).

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

**ThermoFisher**  
SCIENTIFIC