

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

*N. M. Holliday¹, T.C. Lewis¹, D.T. Staats¹, C. C. Knapp¹, S.B. Killian¹, B.J. Olson², C.L. Pike², E.K. Donnerbauer², N.P. Mehta², T.R. Fritsche², A. Gattis³, N. Waugh³, K.Doing³, J.Vlooswijk⁴, A. Cornelisse⁴, E. Scopes⁵, J. Williams⁵, A-M. Leonte⁵, N.E. Scangarella-Oman⁶
¹Thermo Fisher Scientific, Cleveland, OH; ² Marshfield Clinic, Marshfield, WI; ³Centura Health, Colorado Springs, CO; ⁴University Medical Center Utrecht, Utrecht, Netherlands; ⁵Thermo Fisher Scientific, Basingstoke, UK; ⁶GlaxoSmithKline, Collegeville, PA

ABSTRACT

Background: Gepotidacin (GEP) (Glaxo SmithKline, Collegeville, PA), a novel bacterial type II topoisomerase inhibitor, demonstrates activity against gram-positive and gram-negative pathogens associated with conventional and biothreat infections. A study was conducted at 4 different sites to determine the accuracy and reproducibility of GEP 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:** GEP, in a dilution range of 0.015-64µg/mL, was tested on the Sensititre system against 558 recent clinical and challenge isolates. 15 isolates were also included to assess MIC reproducibility. Microorganisms tested included 109 *Escherichia coli*, 143 *Klebsiella* spp., 107 *Enterobacter* spp., 61 *Proteus* spp., 26 *Serratia marcescens*, 101 *Pseudomonas aeruginosa*, and 1 *Acinetobacter baumannii*, and 25 *Moraxella catarrhalis*. 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 published ranges. **Results:** GEP MIC results for the non-fastidious gram-negative organisms tested on the Sensititre system compared with CLSI/ISO BMD for automated and manual reads resulted in 95.5% and 97% essential agreement (EA; +/- 1 log₂ dilution), respectively. Overall agreement for GEP MIC reproducibility (+/- 1 log₂ dilution of the modal MIC) using automated and manual reads was 98.7% and 99.4%, respectively. **Conclusion:** The Sensititre susceptibility system demonstrates an equivalent level of performance compared to the CLSI/ISO BMD method when testing GEP against non-fastidious gram-negative 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 GEP.

INTRODUCTION

Gepotidacin belongs to the triazaacenaphthylene class of antibiotics. This agent binds to DNA gyrase and topoisomerase IV in ways that distinguish it from other quinolone antibiotics. Gepotidacin has demonstrated *in vitro* activity and *in vivo* efficacy against gram-positive and gram-negative pathogens for skin and bio-threat infections including isolates resistant to existing classes of antimicrobials.

This *in vitro* comparison study was done to validate the performance of gepotidacin 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 and International Organization for Standardization (CLSI/ISO). 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 isolates, 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.

Gepotidacin was tested against: (Table 1.)

- 461 recent clinical isolates across the four sites
- 97 challenge isolates at a single testing site
- 15 reproducibility isolates at each site (tested in triplicate over a 3 day testing period)
- 1 Quality Control Strain (ATCC)



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.

● Essential agreement was calculated within plus or minus, one two-fold dilution of the Sensititre dried plates with the reference broth microdilution method.

● Each isolate was tested using a:

- Dried Sensititre 18–24 susceptibility plate containing Gepotidacin (0.015-64µg/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 plates were prepared and tested on each isolate according to the current CLSI/ISO standard method.

Table 1. Organisms Tested	Number Tested
Clinical Isolates (4 sites)	461
CDC Challenge Isolates (one site)	97
Reproducibility Isolates (4 sites) (3 x day for 3 days)	15 (540)
ATCC Quality Control Strains (20 replicates of each strain at 4 sites)	1 (80)
TOTAL	1178

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 2-8X10⁵ cfu/ml, Sensititre 5X10⁴-5X10⁵ cfu/ml

Table 2. Quality Control Strain	Gepotidacin CLSI QC Range (µg/ml)
<i>Escherichia coli</i> ATCC 25922	1-4

Results

Essential agreement for gepotidacin on the Sensititre susceptibility plate compared to the reference microdilution plate was calculated for each read method (Auto and Manual) using the +/- one log₂ dilution standard. Essential agreement rates are shown for non-fastidious gram-negative isolates in **Tables 3 and 4.**

Clinical Isolates and Challenge Organisms

The overall essential agreement for gepotidacin within ±1 log₂ dilution was **97.0%** for the manual method and **95.5%** for the auto read method.

Inter-laboratory Reproducibility

Reproducibility testing results for gepotidacin within ±1 log₂ dilution from the modal MIC was **98.7%** for the auto read method and **99.4%** for the manual read method.

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

The overall essential agreement for gepotidacin within +/- one log₂ dilution, was **97.0%** for the manual read method

Combined Total Isolates

Gepotidacin	Number of Isolates		Essential Agreement		% Essential Agreement	
	All	Evaluable	Total	Evaluable	Total	Evaluable
Organism Group						
<i>Escherichia coli</i>	104	104	102	101	98.1%	97.1%
<i>Klebsiella pneumoniae</i>	94	93	89	88	94.7%	94.6%
<i>Klebsiella oxytoca</i>	43	43	41	41	95.3%	95.3%
<i>Enterobacter cloacae</i>	67	66	66	65	98.5%	98.5%
<i>Klebsiella (Enterobacter) aerogenes</i>	39	39	39	39	100.0%	100.0%
<i>Serratia marcescens</i>	26	26	26	26	100.0%	100.0%
<i>Proteus vulgaris</i>	24	24	24	24	100.0%	100.0%
<i>Proteus mirabilis</i>	36	36	36	36	100.0%	100.0%
<i>Moraxella catarrhalis</i>	25	24	23	22	92.0%	91.7%
<i>Pseudomonas aeruginosa</i>	100	99	95	94	95.0%	94.9%
Total	558	554	541	536	97.0%	96.8%

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

The overall essential agreement for gepotidacin within +/- one log₂ dilution, was **95.5%** for the auto read method

Combined Total Isolates

Gepotidacin	Number of Isolates		Essential Agreement		% Essential Agreement	
	All	Evaluable	Total	Evaluable	Total	Evaluable
Organism Group						
<i>Escherichia coli</i>	104	104	101	100	97.1%	96.2%
<i>Klebsiella pneumoniae</i>	94	93	88	87	93.6%	93.5%
<i>Klebsiella oxytoca</i>	43	43	41	41	95.3%	95.3%
<i>Enterobacter cloacae</i>	67	66	66	65	98.5%	98.5%
<i>Klebsiella (Enterobacter) aerogenes</i>	39	39	37	37	94.9%	94.9%
<i>Serratia marcescens</i>	26	26	24	24	92.3%	92.3%
<i>Proteus vulgaris</i>	24	24	22	22	91.7%	91.7%
<i>Proteus mirabilis</i>	36	36	36	36	100.0%	100.0%
<i>Moraxella catarrhalis</i>	25	23	23	22	92.0%	95.7%
<i>Pseudomonas aeruginosa</i>	100	99	95	94	95.0%	94.9%
Total	558	553	533	528	95.5%	95.5%



RESULTS Cont.

Table 5. Inter-laboratory Reproducibility, % Essential Agreement and ±1 log₂ dilution from the Modal Value

Gepotidacin	Auto Read	Manual Read
Between-site total isolates tested (N)	540	540
Between-site isolates within +/- 1 well from mode (n)	533	537
Between-site reproducibility %	98.7%	99.4%
Total essential agreement (n)	533	537
Essential agreement %	98.7%	99.4%

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/ISO) reference broth microdilution plate when testing gepotidacin against non-fastidious gram-negative clinical and challenge isolates. This study suggests that this is an acceptable method for susceptibility testing of gepotidacin.

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-seventh 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 rapid growing aerobic bacteria involved in infectious diseases (ISO 20776-1 :2006).

These studies were supported by GSK and funded through OTA HHSO100201300011C with HHS/BARDA

© 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