

Thermo Scientific Oxoid M.I.C.Evaluator Strips: An Effective Tool For Staphylococcal MIC Testing

S. Withey.

Thermo Fisher Scientific, Wade Road, Basingstoke, Hants, RG24 8PW, UK

Overview

Purpose: To compare CLSI agar/broth dilution with M.I.C.Evaluator™ (M.I.C.E.™) Strips for staphylococcal minimum inhibitory concentration (MIC) testing.

Methods: Organism MICs when following CLSI agar dilution and M.I.C.Evaluator strip methods were compared.

Results: All antibiotics tested showed >90% essential agreement for streptococci when compared with CLSI agar dilution.

Introduction

Staphylococcus aureus is one of the leading causes of nosocomial infections worldwide¹. Prevalence of *S. aureus* resistance shows regional variation, most notably with methicillin resistance of hospital-acquired *S. aureus* and coagulase-negative staphylococci (CNS). Although CNS are less pathogenic than *S. aureus*, they are important in line-associated bacteraemias and prosthetic device infections where they have been shown to be more resistant than *S. aureus* to teicoplanin².

This study looked at the use of 8 different M.I.C.Evaluator strips as an alternative to the CLSI agar dilution method for the MIC testing of staphylococci. The strips were: amikacin (MA0135), cefotaxime (MA0112), ceftriaxone (MA0122), clindamycin (MA0119), meropenem (MA0121), penicillin (MA0100),



FIGURE 1: M.I.C.E. strips on Mueller-Hinton Agar with 5% sheep blood.

teicoplanin (MA0118) and tigecycline (MA0124).

Methods

Using sterile 0.9% saline, 0.5 McFarland suspensions of methicillin-resistant *S. aureus* (MRSA), methicillin-sensitive *S. aureus* (MSSA), CNS and glycopeptide-intermediate *S. aureus* (GISA) isolates were prepared from cultures grown on Columbia Agar with Horse Blood (PB0122A). Using a multipoint inoculator, the McFarland suspensions were inoculated onto Mueller-Hinton Agar dilution plates. The same suspension was used to create a bacterial lawn on Mueller-Hinton Agar and using sterile forceps M.I.C.E. strips were applied. All plates were incubated at 36±1°C in ambient air following CLSI guidelines³. Results

were used to determine the percentage essential agreement (EA) between the CLSI reference method and M.I.C.E. strips.

Results

Amikacin, cefotaxime, daptomycin and tigecycline M.I.C.E. strips achieved 100% EA with the CLSI agar/broth dilution across all streptococci groups, with the other 3 compounds achieving an EA >90% for the group (as shown in table 1). *S. pneumoniae* were shown to have 100% EA for all 7 different M.I.C.E. strips.

Conclusion

Amikacin, cefotaxime, ceftriaxone, clindamycin, meropenem, penicillin, teicoplanin and tigecycline M.I.C.Evaluator strips offer an accurate, attractive alternative to the labour-intensive traditional agar dilution method for determining the MIC of staphylococci.

References

1. *Clin Infect Dis*. 2001 May 15;32 Suppl 2:S114-32. Survey of infections due to *Staphylococcus* species: frequency of occurrence and antimicrobial susceptibility of isolates collected in the United States, Canada, Latin America, Europe, and the Western
2. *Livermore DM*. Antibiotic resistance in staphylococci. *Int J Antimicrob Agents*. 2000 Nov;16 Suppl 1:S3-10.
3. CLSI Performance Standards for Antimicrobial Susceptibility Testing; Twenty-Second Informational Supplement M100-S22 January 2012

TABLE 1: Essential agreement (shown as a percentage) for each organism group. The number of organisms tested is shown in brackets.

M.I.C.E. Strip	Organism				
	All Staphylococci	MRSA	MSSA	CNS	GISA
Amikacin	100 (n=61)	100 (n=30)	100 (n=11)	100 (n=20)	
Cefotaxime	100 (n=30)	100 (n=20)	100 (n=5)	100 (n=5)	
Ceftriaxone	96.66 (n=30)	100 (n=10)	100 (n=9)	90.91 (n=11)	
Clindamycin	100 (n=101)	100 (n=33)	100 (n=33)	100 (n=35)	
Meropenem	93.33 (n=30)	100 (n=10)	100 (n=9)	81.82 (n=11)	
Penicillin	94.06 (n=101)	91.67 (n=36)	97.14 (n=35)	93.33 (n=30)	
Teicoplanin	93.38 (n=150)	96 (n=49)	100 (n=63)	78.95 (n=38)	
Tigecycline	97.01 (n=67)	100 (n=30)	100 (n=9)	95 (n=20)	87.5 (n=8)