


Thermo Scientific Sensititre plate setup guide

For *Streptococcus* spp. and *S. pneumoniae*

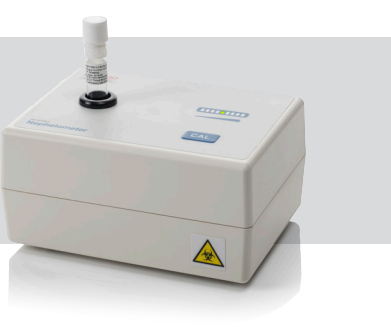
1

Pick 3-5 colonies from agar plate and emulsify in 5 mL Thermo Scientific™ Sensititre™ Mueller Hinton Broth



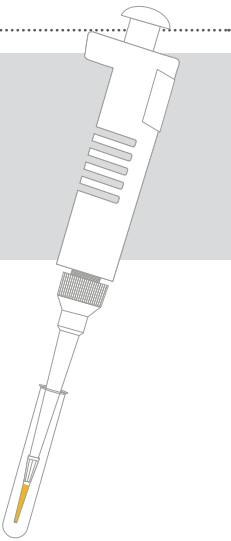
2

Adjust suspension to 0.5 McFarland equivalent using the Thermo Scientific™ Sensititre™ Nephelometer



3


Transfer 100 µL to 11 mL Thermo Scientific™ Mueller-Hinton Broth with Lysed Horseblood



4

Inoculate the panel with 100 µL per well using the Thermo Scientific™ Sensititre AIM™ Automated Inoculation Delivery System

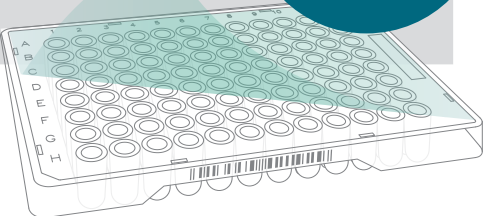
Use additional AIM Systems for increased throughput



5

Cover all wells with the perforated adhesive seal. Incubate in at 34-36 °C aerobically for 20-24 hours*


Complete panel set up (steps 1-5) and place into the incubator (step 6) within 30 minutes



* Sealed plates can be stacked up to 3 high in an incubator

6

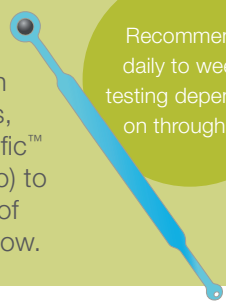
Read results using the Thermo Scientific™ Sensititre™ Vizion™ System or Thermo Scientific™ Sensititre™ Manual Viewbox



To meet your laboratory quality assurance guidelines

Quality control testing

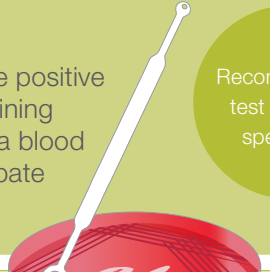
Analyze appropriate quality control organisms with known MIC values (e.g. ATCC strains, available from Thermo Scientific™ Culti-Loops™ product portfolio) to confirm reliable performance of your consumables and workflow.



Recommended daily to weekly testing depending on throughput

Colony count


After inoculating the broth, transfer 100 µL from the remaining broth into a new 11 mL Mueller-Hinton broth. Inoculate 100 µL from this dilution onto a non-selective agar plate. Incubate overnight and count the colonies. Expected results should fall into the range of 50-500 colony forming units.



Regular testing interval

Purity check

Streak 1 µL from the positive control well or remaining broth directly on to a blood agar plate and incubate overnight.



Recommended test on every specimen

Ordering information		
Product	Quantity	Cat. No.
Thermo Scientific 0.5 Polymer McFarland Standard	Each	E1041
Thermo Scientific Doseheads for Sensititre AIM	100/Pack	E3010
Thermo Scientific Mueller-Hinton Broth	100 x 5 mL	T3462-05
Thermo Scientific Mueller-Hinton Broth with Lysed Horseblood	10 x 11 mL	CP-112
Thermo Scientific Mueller-Hinton Broth with Lysed Horseblood for autoread	10 x 11 mL	CP114-10
Thermo Scientific Culti-Loops Quality Control Organisms	Contact your Thermo Fisher Microbiology representative for options	

ATCC Licensed Derivative

™ The ATCC Licensed Derivative® Emblem, the ATCC Licensed Derivative® word mark, and the ATCC catalog marks are trademarks of ATCC. Thermo Fisher Scientific Inc. is licensed to use these trademarks and to sell products derived from ATCC® cultures.

Find out more at thermofisher.com/sensititre

ThermoFisher
SCIENTIFIC

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Not all products are available in all countries. Please consult your local sales representative for details.
LT2553A 0820