


# Thermo Scientific Sensititre plate setup guide

## Non-fastidious Gram negative and Gram positive organisms

**1** Pick 3-5 colonies from agar plate and emulsify in 5 mL Thermo Scientific™ Demineralised Water



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

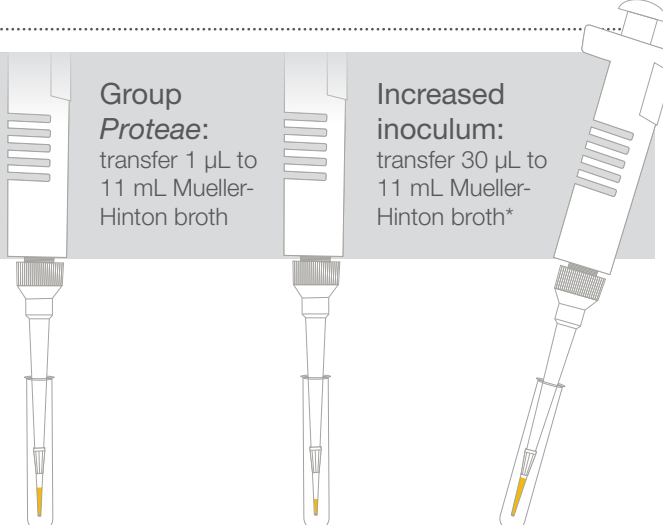


**3** Standard inoculum: transfer 10 µL to 11 mL Mueller-Hinton broth

Group *Proteae*: transfer 1 µL to 11 mL Mueller-Hinton broth


Increased inoculum: transfer 30 µL to 11 mL Mueller-Hinton broth\*

\*Depending on the strains tested an increase to 30 µL may aid in detecting resistance mechanisms



**4** Inoculate the panel with 50 µL per well using the Thermo Scientific™ Sensititre AIM™ Automated Inoculation Delivery System

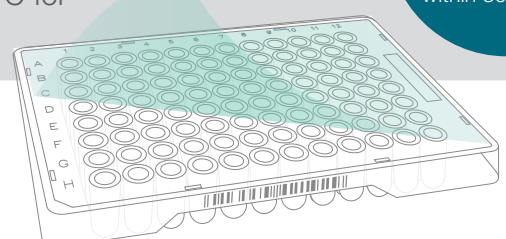
Use additional AIM Systems for increased throughput




**5** Cover all wells with the adhesive seal. Incubate aerobically at 34-36°C for 18-24 hours\*

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

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



**6** Read the results using Thermo Scientific™ Sensititre™ ARIS HiQ™ System, Sensititre™ Optiread™ system, Sensititre™ Vizion™ system or 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

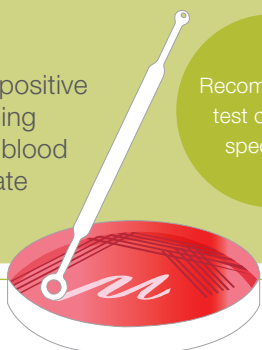
After inoculating the broth, transfer 100 µL from the remaining broth into a new 11 mL Mueller-Hinton broth. Inoculate 100 µL from this dilution onto a non-selective agar plate. Incubate overnight and count the colonies. Expected results should fall into the range of 50-500 colony forming units.

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 Demineralised Water	100 x 5 mL	T3339
Thermo Scientific Mueller-Hinton Broth	100 x 11 mL	T3462
Thermo Scientific Culti-Loops for quality control testing	Contact your Thermo Fisher Microbiology representative for options	



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](http://thermofisher.com/sensititre)

**ThermoFisher**  
SCIENTIFIC