



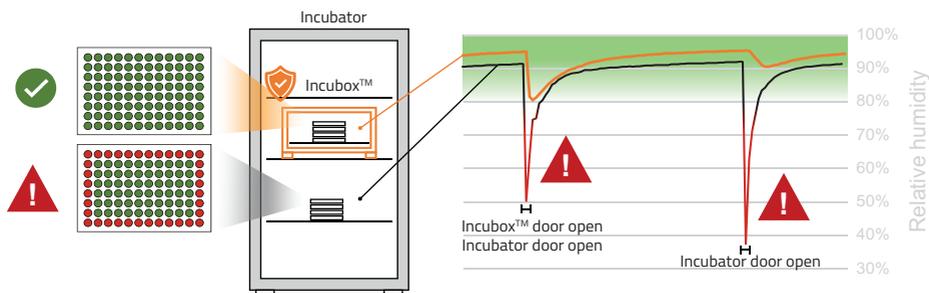
InSphero Incubox™

Quick Start Guide

Thank you for choosing InSphero's Incubox™ to maintain stable culture conditions of your cell culture. This Quick Start Guide contains important information to get you started immediately. For detailed instructions please refer to the Product Manual and additional resources on shop.insphero.com.

Introduction

Edge effects due to media evaporation are a common problem in long-term cell cultures, especially in busy labs with many projects in which incubator doors are frequently opened and closed. The Incubox™ creates a high-humidity microclimate inside conventional cell culture incubators and protects your sensitive plates from large humidity drops and fluctuations – the major cause for evaporation effects.



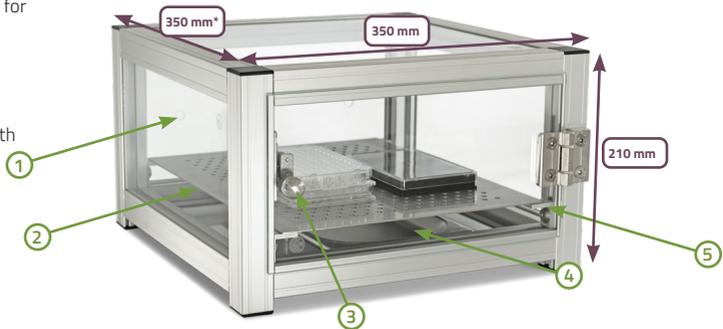
- **Maintain a stable medium volume and accurate compound concentration** due to minimal evaporation over time.
- **Reduce well-to-well variation** by minimizing evaporation-induced reagent gradients across the plate.
- **Use all wells of your plate** by eliminating edge effects in your culture plate.
- **Save on plasticware** by using up to 38% more wells of your plate.

Key Features

1. Fits in a wide range of cell culture incubators.
2. Holds up to 20 microplates.
3. Temperature-stable and compatible with decontamination procedures.
4. Inert materials are compatible with a wide range of solvents.
5. Separate water bath to quickly re-establish high humidity.
6. Gas-equilibration holes ensure stable O₂ and CO₂ conditions.
7. Easy one-handed door opening and closing.

Components of the Incubox™

1. Three holes on backside for gas exchange
2. Stainless steel shelf
3. Magnetic doorknob
4. Stainless steel water bath
5. Shelf ledge



*Depth incl. doorknob 375 mm

Figure 1: Photo of Incubox™ filled with Akura™ Plates.

Step-by-Step Protocol

Please follow the steps below to ensure proper function of the Incubox™ and optimal culture conditions:

1. Carefully clean all components of the Incubox™ with Ethanol to remove dust.
2. Place the Incubox™ directly on a shelf inside your incubator. Make sure it stands stable and at least 5 cm away from the backside of the incubator. Don't block the holes on the backside. We recommend placing the Incubox™ on one of the center shelves for easy access of your plates and to avoid blocking the ventilator on the top of some incubator models.
3. **Optional:** Use the high-temperature sterilization cycle of your incubator to ensure complete elimination of microbial contaminants.
4. Remove the shelf from the Incubox™ and fill water bath with approximately 500 ml of sterile water including a water bath stabilizer, i.e. Aqua resist. (To re-fill water bath, remove plates, lift the shelf to one side and re-fill bottom container).
5. Place the shelf back onto the ledge.
6. Close Incubox™ and incubator. Wait at least one hour to ensure stable temperature and high relative humidity.
7. Add your plates and/or culture dishes and start your experiment.

For more detailed information, please refer to the Incubox™ Product Manual.



InSphero AG
Schlieren, Switzerland

InSphero Inc.
Brunswick, ME, USA

If you have more questions please refer to the FAQs section at shop.insphero.com

Sign up for the latest news and updates at insphero.com/newsletter

Follow us on LinkedIn and Twitter

InSphero is ISO 9001:2015 certified

All rights reserved. © 2021 InSphero AG. 3D InSight, Akura, InFloat and Incubox are trademarks of InSphero AG. For life science research only. Not for use in diagnostic procedures.