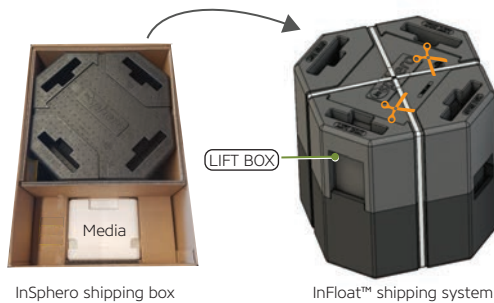


## Assay-ready solutions

# Unpacking 3D InSight™ Microtissues

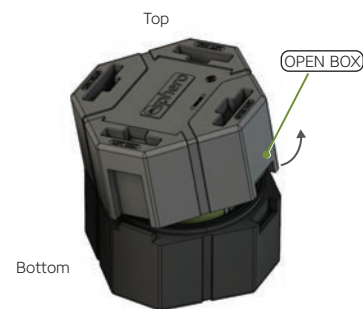
Assay-ready 3D InSight™ microtissues are delivered in Akura™ 96 plates carefully packaged for air transport in the InFloat™ shipping system, which is specially engineered to protect live microtissues from damage and keep them at a physiologically appropriate temperature during transit. Follow these 6 steps to open the black InFloat™ container and access your microtissues, then refer to the *Quick Start Guide* for instructions on handling microtissues before conducting assays.

- 1 Pick up the black InFloat™ container using the LIFT BOX handles.



Remove the documents from the InSphero shipping box, transfer the black InFloat™ container from the shipping box to a flat work surface and cut the straps at the locations indicated. Retrieve the media from the white box and store as indicated on the label.

- 2 Open the container using the OPEN BOX handles.



Carefully pull on the OPEN BOX handles, first on one side, then the other, until the container top is free. **Note:** The InFloat™ shipping system is black; the grey and green colors in these illustrations are used only to help distinguish different components.

- 3 Turn the container top upside down and place the floating sphere inside the top.



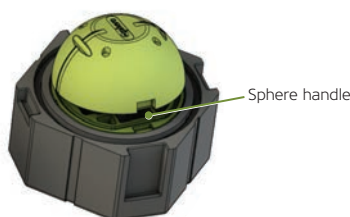
**!** Be careful not to tilt or drop the sphere when moving it.

- 4 Cut the strap on the sphere at the indentation and put the container bottom aside.

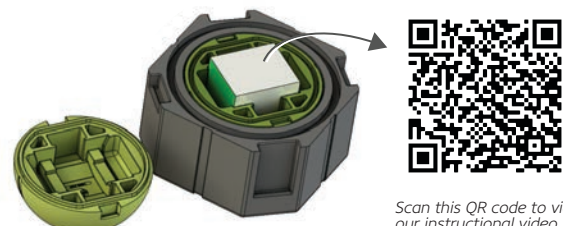


The liquid in the container bottom is water and may be discarded in a sink.

- 5 Open the sphere using the indented handles on the side of the sphere.



- 6 Remove the microtissue box and follow the instructions in the *Quick Start Guide*.



Scan this QR code to view our instructional video.