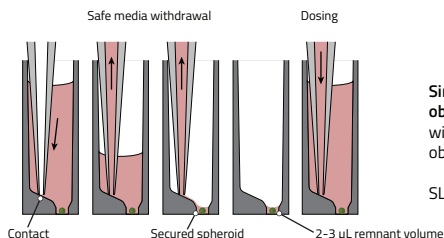
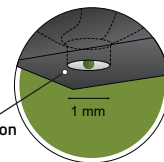


Akura™ 384 Microplate Solutions

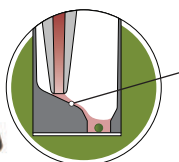
Unique Well Design



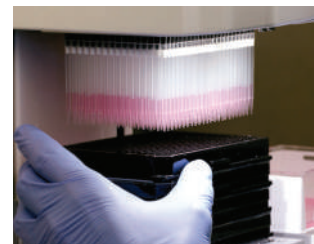
Simplify spheroid localization, observation, and ROI identification with standard 1 mm diameter observation chambers



SLAS/ANSI compliance assures automation compatibility

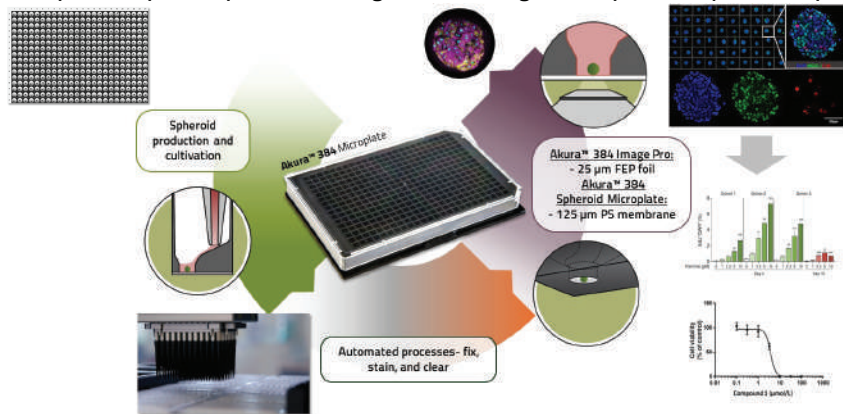


Prevent unintended spheroid aspiration with the unique SureXchange™ ledge design that shields the spheroid and retains consistent minimal remnant media volumes during medium exchange



End-to-end solution

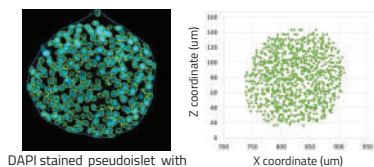
Incorporates spheroid production, long-term culturing and endpoint analysis in one plate



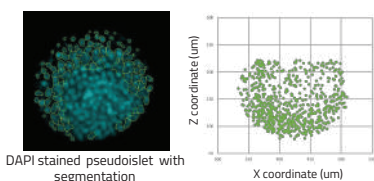
Akura™ 384 Spheroid Microplate: The Ultimate Spheroid Screening Plate

Image Segmentation

Akura™ 384 ImagePro
reduced light scatter, easy to segment



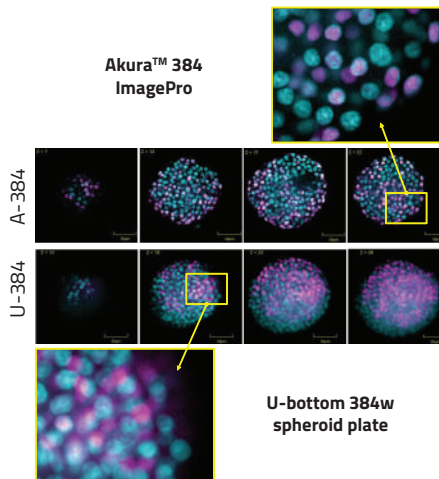
U-bottom 384w spheroid plate
excessive light scatter, difficult to segment



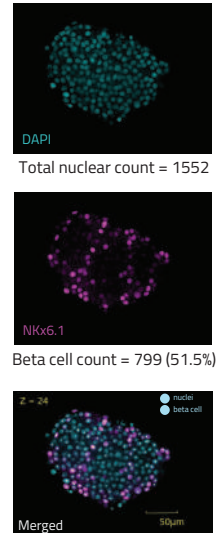
U-bottom plates produce significant light scatter and reduce practical imaging depth

Target Co-localization

Akura™ 384 ImagePro



Quantification of NKX6.1+ islet cells



Akura™ 384 ImagePro enables accurate nuclear segmentation and detection of co-localized nuclear signal throughout the entire spheroid volume resulting in the accurate quantification of the beta cell population within the pseudoislets

Akura™ 384 ImagePro:
Enables Superior High-Content Imaging Quality