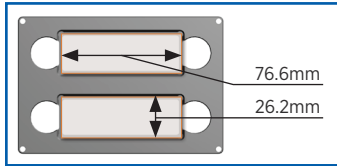


EVOS® VESSEL HOLDERS & STAGE PLATES

All models

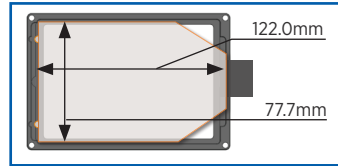
AMEPVH001

holds two 25mm x 75mm standard microscope slides, chamber slides, etc.



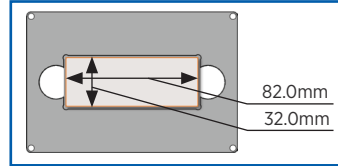
AMEPVH008

holds one Greiner T-75 flask; 75cm²



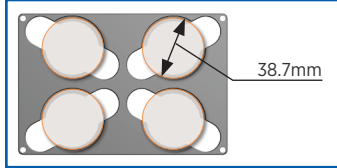
AMEPVH017

holds one KOVA Glasstic Slide 10



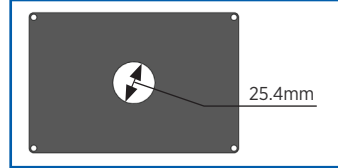
AMEPVH002

holds four 35mm Petri dishes



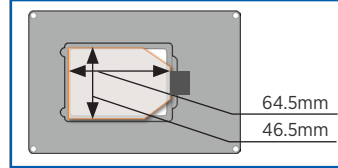
AMEPVH009

Universal stage insert



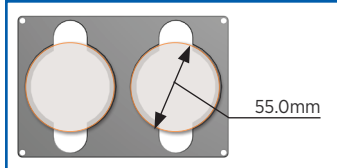
AMEPVH018

holds one Nunc T-25 flask; 25cm²



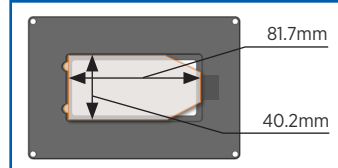
AMEPVH003

holds two 60mm Petri dishes



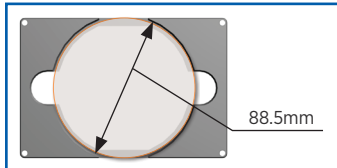
AMEPVH010

holds one BD/Greiner T-25 flask; 25cm²



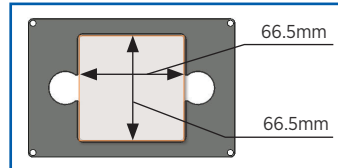
AMEPVH004

holds one 100mm Petri dish



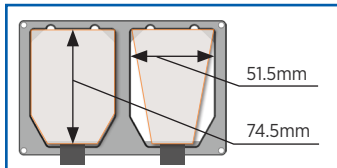
AMEPVH011

holds one NUNC/SPL IVF 4-well dish



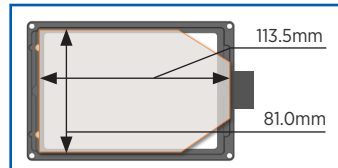
AMEPVH005

holds two 25cm² flasks; rectangular or triangular



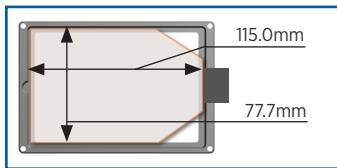
AMEPVH012

holds one SPL T-75 flask; 75cm²



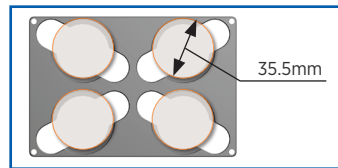
AMEPVH006

holds one Nunc T-75 flask; 75cm²



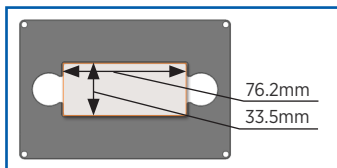
AMEPVH013

holds four Ibidi 35mm Petri dishes



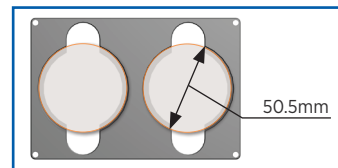
AMEPVH007

holds one Hemocytometer



AMEPVH014

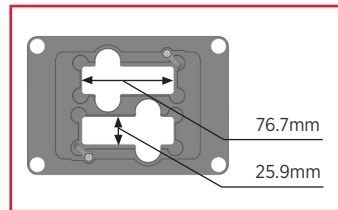
holds two Ibidi 50mm Petri dishes



FL Auto

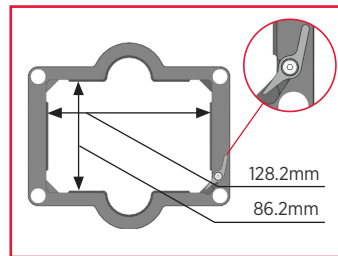
AMEPVH021

Securely holds two 25mm x 75mm standard microscope slides, chamber slides, etc.



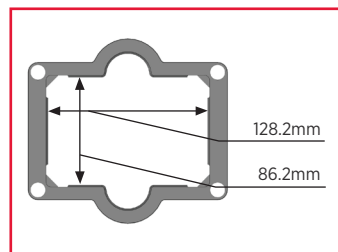
AMEPVH022

Intermediate Plate for Automated Stage. Securely holds multiwell vessels with convenient lever Adaptor for AMEPVH001 and AMEPVH009



AMEPVH023

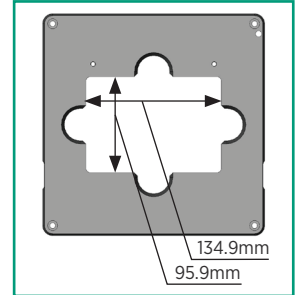
holds multiwell vessels Adaptor for AMEPVH001 and AMEPVH009



FL, FL Color & XL

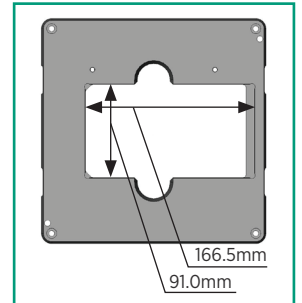
AMEP4684

Stage Plate for heating tray, Tokai Hit #MATS-UAXKW-D



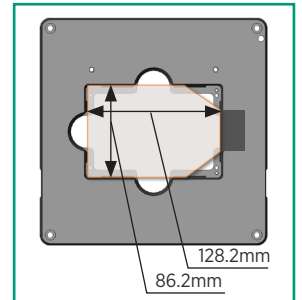
AMEP4685

Stage Plate for heating stage, BioFlux by Fluxion



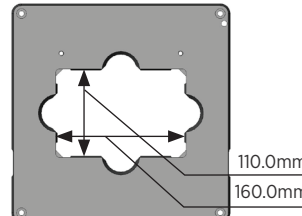
AMEP4686

Stage Plate for multiwell vessels; also holds one Corning T-75 flask



AMEP4691

Stage Plate with 110mm x 160mm opening (Use with AMEP4692 for standard sizes)



AMEP4692

Stage Plate Adaptor 110mm x 160mm opening for standard size (sold separately)

