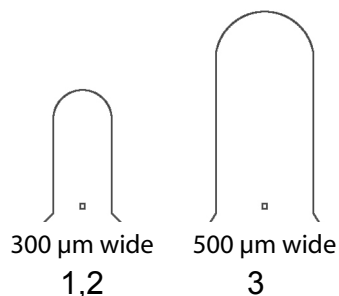


MicroTools Kit 1:

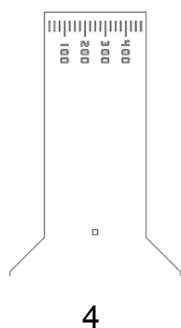
MicroTools™ for Manipulation and Measurement

20 soft, flexible and durable tools to simplify your sample handling tasks



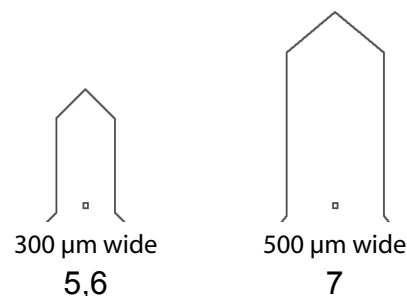
MicroSpoons™

Push, detach from surfaces, separate from a cluster, lift, transfer
 Push down to flatten tip and slide under sample.



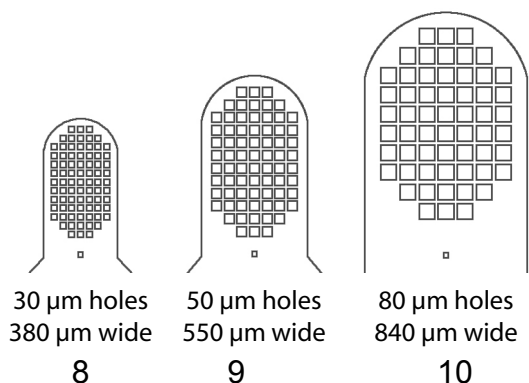
Horizontal MicroRuler™

500 µm wide with 25 µm increments
 For linear measurements



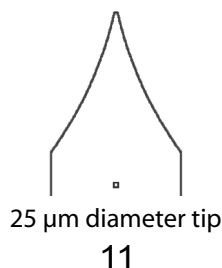
MicroChisels™

Push, detach from surfaces, separate from a cluster, lift, transfer
 Push down to flatten tip and slide under sample.



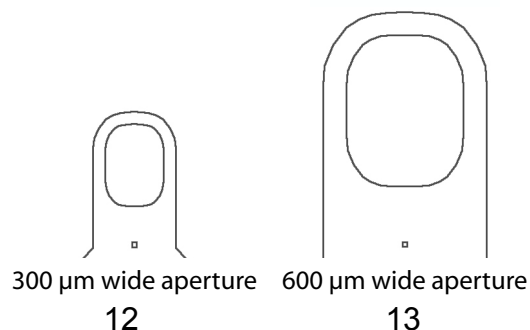
MicroSieves™

Sieve, lift, transfer and soak



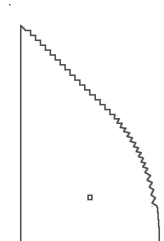
MicroTips™

Poke, separate, lift and reposition on a MicroMount



MicroLassos™

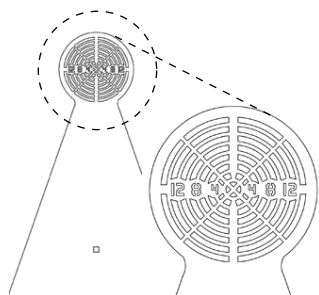
Lasso and pull to detach from surfaces



25 µm teeth
 840 µm wide

MicroSaws™

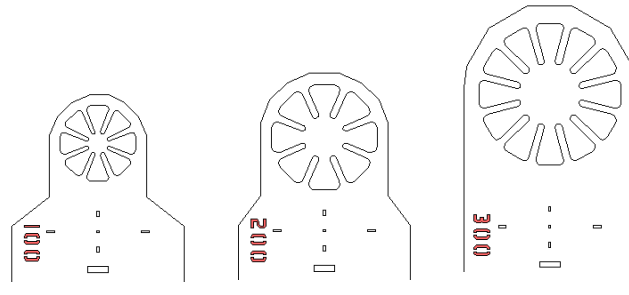
Cut protein skins and dense lipid phases



16-17

Small MicroProtractor™

320 µm wide with 20 µm and 45° increments
 For radial and angular measurements



18

19

20

MicroGrippers™

Push down onto your sample to firmly capture it for room-temperature crystallography..