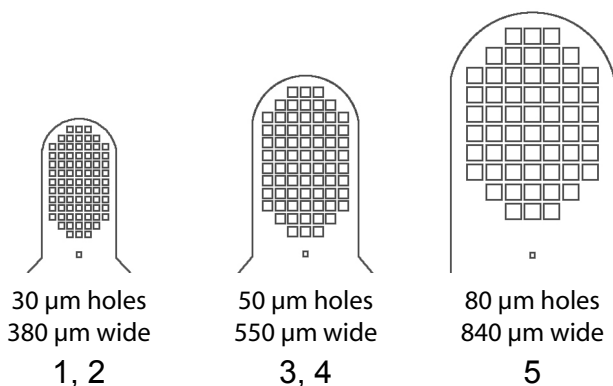


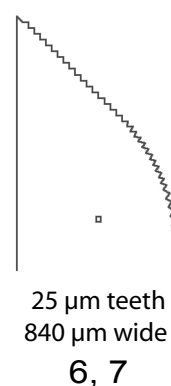
MicroTools™ for Manipulation: Kit 2

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



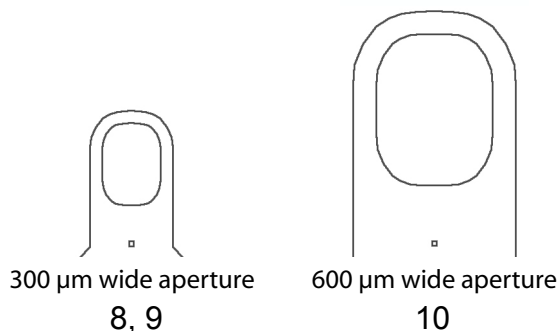
MicroSieves™

Sieve, lift, transfer and soak



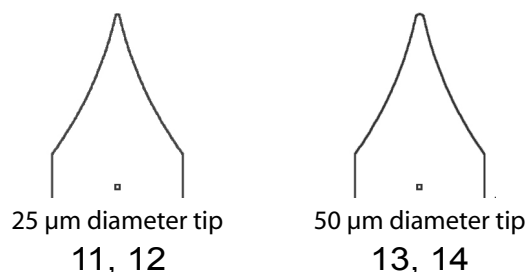
MicroSaws™

Cut protein skins and dense lipid phases



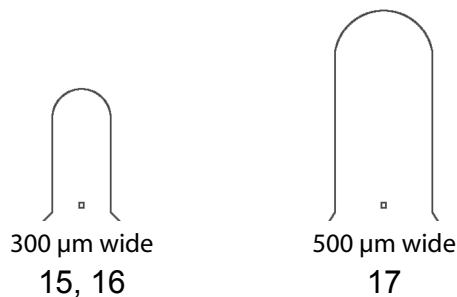
MicroLassos™

Lasso and pull to detach from surfaces



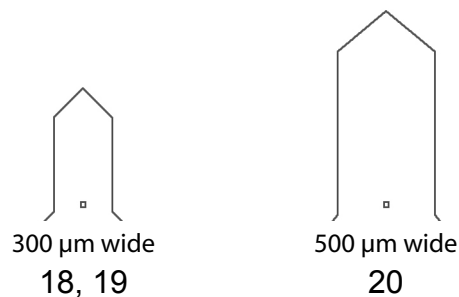
MicroTips™

Poke, separate, lift and reposition on a MicroMount



MicroSpoons™

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



MicroChisels™

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