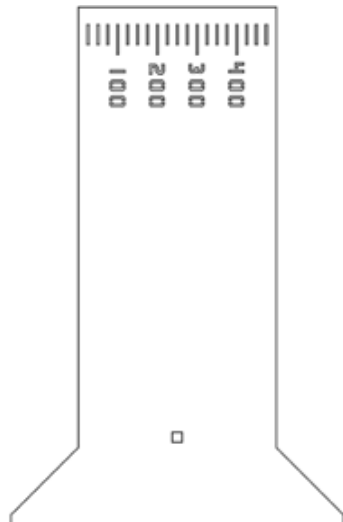


MicroTools™ for Measurements: Kit 3

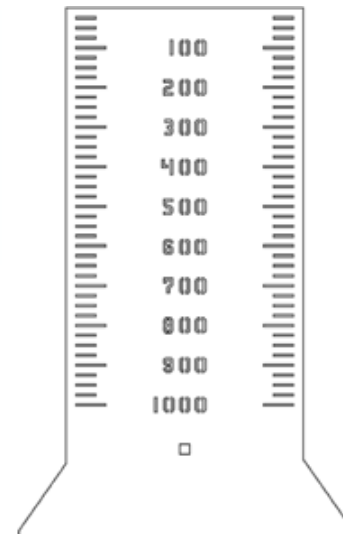
20 soft, flexible tools for linear, radial and angular measurements



1-5

Horizontal MicroRuler™

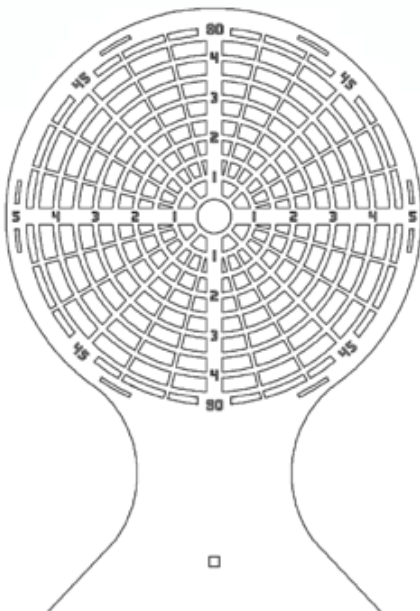
500 µm wide with 25 µm increments
For linear measurements



6-10

Vertical MicroRuler™

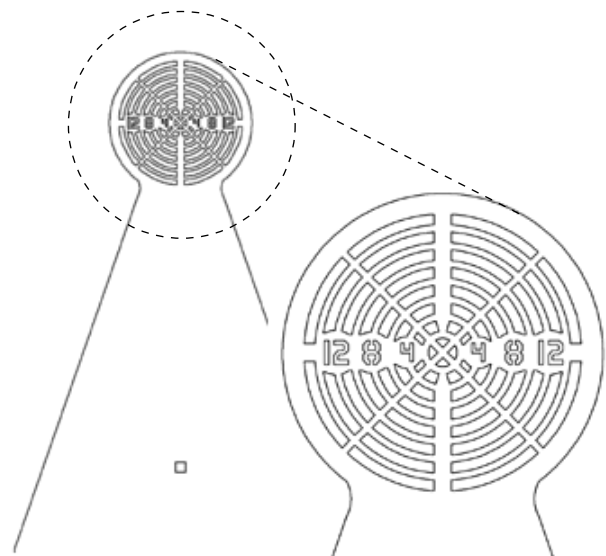
1000 µm length with 25 µm increments
For linear measurements



11-15

Large MicroProtractor™

1000 µm wide with 50 µm and 15° increments
For radial and angular measurements



16-20

Small MicroProtractor™

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