

Pipette Tip *1000 ul*Catalog number: L55008T
Unit size: 500**Applications**

This pipette tip is best used with Medfuture Variable Volume Single Channel Pippettes of 100-1000 ul (L55008).