

Pipette Tip *10 ml*Catalog number: L55010T
Unit size: 300**Applications**

This pipette tip is best used with Medfuture Variable Volume Single Channel Pippettes of 1-10 ml (L55010 and L55011).