

Pipette Tip *200 ul*

Catalog number: L55003T
Unit size: 1000

Applications

This pipette tip is best used with Medfuture Variable Volume Single Channel Pippettes of 5-200 ul (L55003 and L55007).