

### Get the most out of your electronic pipette

Work faster, smarter & improve reproducibility in cellular and molecular biology



#### 96-well or 384-well plate assays

- PCR
- Enzyme assays
- Cell culture – drug or reagent additions



#### Use "Plate Tracker"

For keeping track of one's position when pipetting on micro-well plates. Never accidentally pipette into the wrong well again.



#### Cell culture: Mixing and seeding cells

Use Pipetting with Mixing mode: Adjustable pipetting speed for sensitive cell types. Number of mixings can be varied to ensure even distribution of cells.



#### PCR, cell assays, immunoassays, fluorescence assays, ELISA

Use Multi-Dispensing mode: Used for long pipetting series. Ensures fast and reproducible pipetting. Can be combined with Plate tracker.



#### Avoiding bubbles

Use Reverse Pipetting mode: Adjust the excess volume to save on expensive reagents. Used for fluorescence-, absorbance-, or cell-based assays, proteins, foamy or viscous liquids.



#### Checking protein yield, sample or reagent volumes

Use manual mode: Manually controlled aspiration for accurate and rapid checking of volume.



#### Standard curves and calibration curves

Use "Sequential Pipetting" mode: Ensures faster preparation of curves in spectrometric or colorimetric assays and immunoassays.



#### Sample pooling

Use Multi-Aspiration mode: Efficient and fast sample pooling. Used for NGS, PCR, protein assays and changing media on cell cultures in 96-well plates.