

Use of Thermo Scientific FinnpiPETTE Novus Electronic Pipettes for Accurate and Efficient Preparation of Standard Curves

LHC Research Group, Thermo Fisher Scientific, Vantaa, Finland

Key Words

- Analyte concentration
- Bovine Serum Albumin (BSA)
- Thermo Scientific FinnpiPETTE Novus Electronic Pipette
- Thermo Scientific FinnpiPETTE Manual Pipette
- Thermo Scientific Varioskan Multimode Microplate Reader
- Standard Curve



Introduction

Accurate preparation and measurement of a standard curve is essential for the precise determination of analyte concentration. Because preparing standard curves is a daily routine in many biology and chemistry laboratories, pipette selection for this task is critical.

Examples of applications requiring standard curves

- Total protein determinations
- DNA determinations
- Glucose assays
- ELISA assays
- Enzymatic assays

This application note discusses the use of a Thermo Scientific FinnpiPETTE Novus Electronic Pipette for the preparation of standard curves. With 10 different functions, the versatile Novus Electronic pipette can perform most every standard curve required. The pipette's performance was compared to a manual Thermo Scientific FinnpiPETTE model to determine if there were appreciable differences in pipetting speed and accuracy during curve preparation.

Materials and Methods

1. Thermo Scientific FinnpiPETTE Novus Electronic Pipettes (30–300 μ l, 0.5–5 ml) (Cat. Nos. 46200500 and 4620070, respectively)
2. Thermo Scientific FinnpiPETTE Novus Manual Pipette (0.5–5 ml) (Cat. No. 4600110)
3. Thermo Scientific Finntip Pipette Tips (300 μ l, 5 ml) (Cat. Nos. 94060510 and 9402070, respectively)
4. Test Tubes
5. Black 96-well Thermo Scientific Microtiter Microplates (Cat. No. 9502867)
6. Bovine Serum Albumin (Sigma-Aldrich®, A7030)
7. Dulbecco's Phosphate Buffered Saline (Sigma-Aldrich, D1408)
8. Thermo Scientific Fluoraldehyde o-Phthalaldehyde Reagent Solution (OPA, 0026025)
9. Thermo Scientific Varioskan Multimode Microplate Reader (Cat. No. 5250000)

Testing was conducted using a 5 ml Novus Electronic Pipette and a 5 ml FinnpiPETTE manual pipette. BSA dilutions ranging from 12.5 to 100 μ g/ml were prepared with both pipettes.

Testing with the Novus pipette was conducted in two different ways. First, curves were prepared using the "Forward" (standard) pipetting function, then curves were prepared using the "Dilution & Mix" function with the pipetting speed set at "7" for both aspiration and dispensing.

The "Dilution & Mix" mode enables selected volumes of two liquids to be dispensed and mixed. After taking up the diluting buffer, an air buffer is aspirated before taking up the sample into the same pipette tip. The trigger button is then pressed and held, causing the pipette to "mix" (repeatedly aspirate and dispense) 70% of the selected volume.

The time required to prepare each dilution series was recorded (Fig. 3). The protein concentration was determined with the OPA-solution that allows for fast quantitation of protein or peptide in solution. This solution reacts with primary amino acids, resulting in highly fluorescent isoindole derivatives.

100 μ l of the BSA dilutions was pipetted as three replicates into black Microtiter® microplates, then 100 μ l of OPA solution was added. The microplates were centrifuged at 2,000 rpm for 2 minutes. Fluorescence was measured using the Varioskan® multimode microplate reader. Standard curves were prepared on three different days to verify repeatability.

Linear lines were used to connect the measurement points, then R²-values were calculated. An R²-value is a measure of the exactness of the linear regression; "1" represents a perfect fit between the data and the line connecting them, while "0" represents no statistical correlation between the data and the line.

Results

The BSA standard curves prepared with a Novus electronic pipette and those prepared with a Finnpiquette manual pipette are shown in Figures 1 and 2, respectively. The standard curves prepared with the electronic pipette using the forward pipetting function were similar to those prepared with the Dilute and Mix mode. While the R²-values for all standard curves were > 0.99, indicating a good correlation between the measurement points to a linear line, the curves prepared with the Novus electronic pipette were more convergent than those prepared with the manual pipette.

The time required to prepare the dilution series ranged from 4.5 to 6.5 minutes (Fig. 3). The Novus electronic pipette was 20–30% faster than the manual pipette, an advantage especially important when using sensitive analytes that can degrade or oxidize when dispensed slowly. The Novus electronic pipette's Dilute and Mix mode enabled faster preparation of standard curves and also resulted in a more efficient workflow. The need for a vortexer was eliminated, which reduces the risk of repetitive strain injuries (RSI).

Conclusion

The versatile Finnpiquette Novus Electronic Pipette was shown to be an excellent tool for the preparation of standard curves. The electronic pipette demonstrated the ability to perform more accurate curves in 20–30% less time than with a manual pipette. Curve accuracy was due in part to the electronic pipette's index finger operation, which helps to maintain optimal pipette positioning, and its motorized piston movement, which ensures constant pipetting speed.

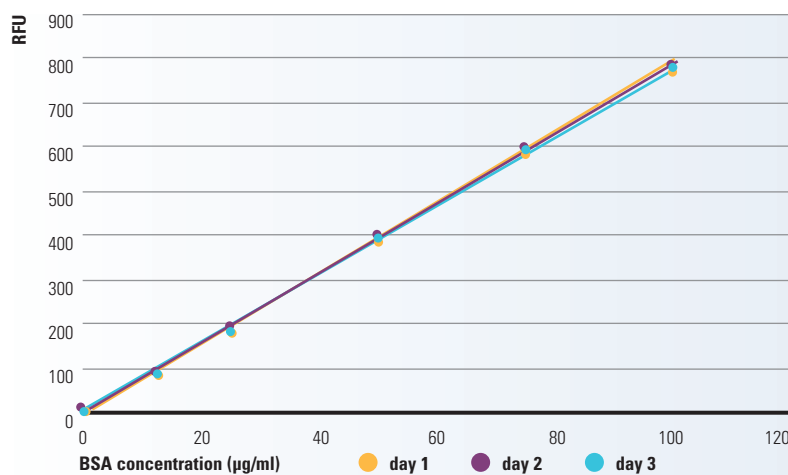


Figure 1. BSA standard curves prepared on three different days with a Finnpiquette Novus Electronic Pipette using the Dilute and Mix mode

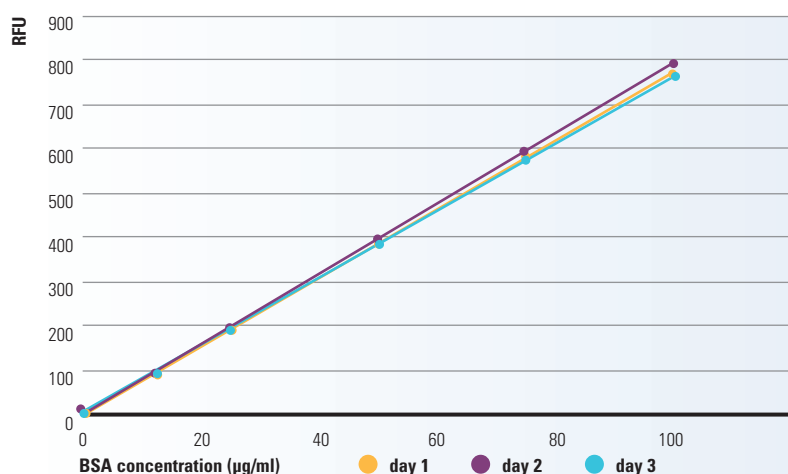


Figure 2. BSA standard curves prepared on three different days with a manual Finnpiquette using the forward pipetting technique

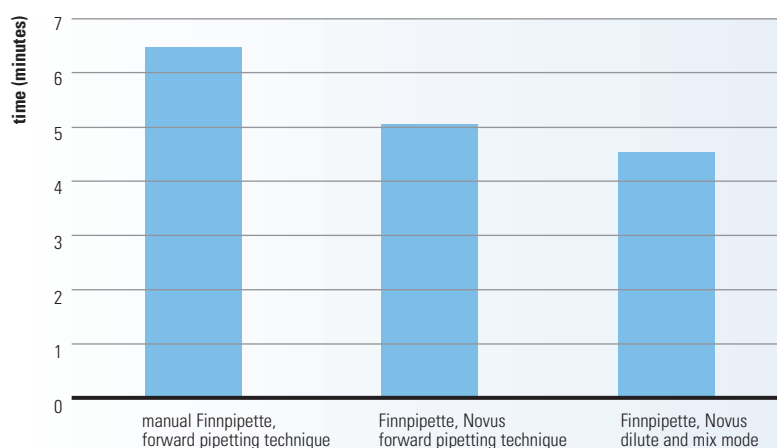


Figure 3. Time required to prepare a dilution series with a manual Finnpiquette and a Novus electronic pipette using different modes

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

North America:

USA / Canada
+1 800 522 7763

Europe:
Austria
+43 1 801 40 0

Belgium
+32 2 482 30 30

France
+33 2 2803 2180

Germany national toll free
08001-536 376

Germany international
+49 6184 90 6940

Italy
+39 02 02 95059 448

Netherlands
+31 76 571 4440

Nordic countries
+358 9 329 100

Russia/CIS
+7 (495) 739 76 41

Spain/Portugal
+34 93 223 09 18

Switzerland
+41 44 454 12 12

UK/Ireland
+44 870 609 9203

Asia:
China

+86 21 6885 4588 or
+86 10 8419 3588

India toll free
1800 22 8374

India
+91 22 6716 2200

Japan
+81 45 453 9220

Other Asian countries
+852 2885 4613

Countries not listed:
+49 6184 90 6940 or
+33 2 2803 2180

www.thermoscientific.com/finnpiquette

© 2010 Thermo Fisher Scientific Inc. All rights reserved. Sigma-Aldrich is a registered trademark of Sigma-Aldrich. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Lit. no. 151 7390