

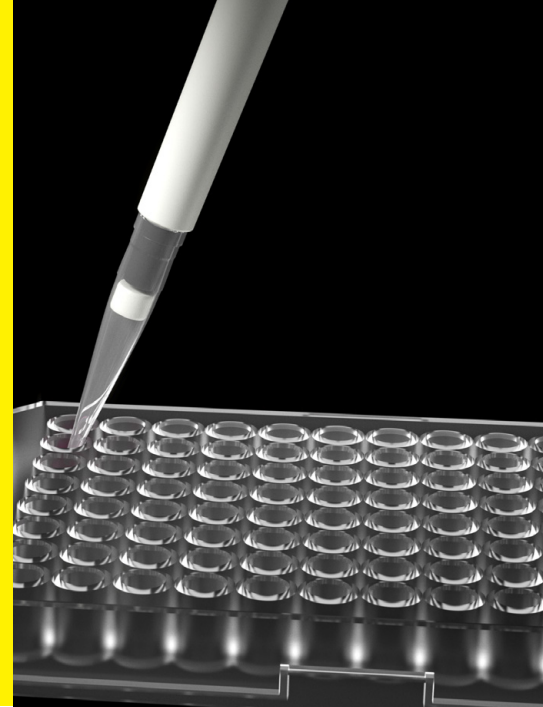
## Pipetting Software: Enhancing Efficiency and Accuracy

Pipetting is a daily task in laboratory operations, requiring frequent adjustments to the settings and other routine functions. Traditional pipetting methods often involve manual adjustments and repetitive tasks, which can lead to human error, inconsistencies, and inefficiencies. There is a clear need for software tools designed for streamlining pipette and protocol management to simplify these processes, ensure accuracy, and enhance productivity in the lab.

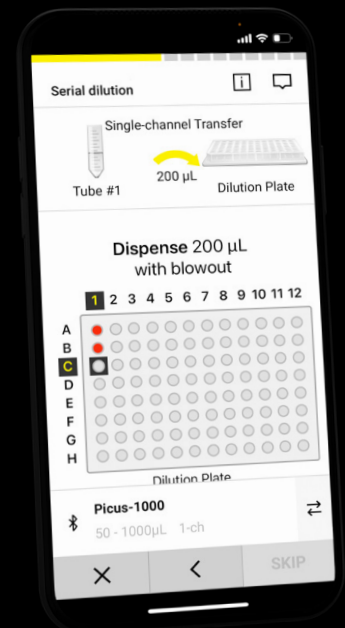
The Sartorius Pipetting software is a companion tool designed to work with the Sartorius Picus® 2 pipette, enhancing workflow management and other critical tasks. This intuitive and user-friendly software provides step-by-step guidance for pipetting workflows, ensuring accuracy and reproducibility for both experienced scientists and novices

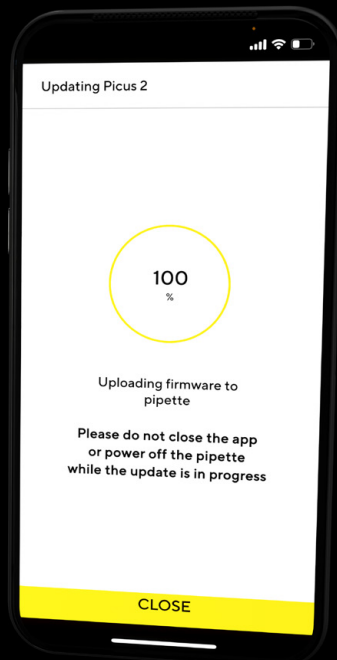
### Efficiency Benefits

- **Automated Workflow Management:** Define and execute complex protocols without manual adjustments, reducing repetitive tasks and minimizing human error.
- **Real-Time Guidance:** Get step-by-step instructions during pipetting procedures, ensuring correct sequence and appropriate volumes.
- **Bluetooth Connectivity:** Easily connect to the Picus® 2 pipette via Bluetooth, facilitating direct communication and automatic adjustment of pipette settings based on the selected protocol.
- **Comprehensive Protocol Library:** Access to a growing library of predefined protocols, simplifying sample preparation with clear instructions for common tasks.
- **Traceability and Record Keeping:** Automatically log all executed workflows, creating a reliable record of pipetting activities to support your compliance requirements.
- **User-Friendly Interface:** Quickly learn how to operate the Picus® 2 effectively, even with minimal experience.



Download Now  
at No Cost:



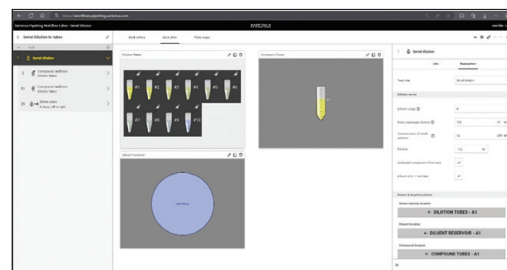


## Did You Know?

With the Sartorius Pipetting mobile application, you can update your Sartorius Picus® 2 pipettes, allowing you to enjoy the latest device improvements and new features

## How Does it Work?

- **Plan or Edit Workflows:** Use the Sartorius Pipetting workflow editor, accessible via a web browser on a PC or laptop, to plan a new pipetting workflow or edit existing example workflows. You can also share these workflows with colleagues for collaborative efficiency.
- **Execute Workflows via Mobile App:** Your workflows are visible in the Sartorius Pipetting mobile application, where you can initiate runs. Some workflows feature adjustable parameters, such as the number of samples or pipetted volume, making them highly versatile for everyday lab work.
- **Automated Pipette Settings:** The system automatically transfers pipetting settings to the most suitable Picus® 2 pipette in your set, eliminating the need for manual adjustments and ensuring consistent execution of steps. It also generates a run report for each workflow execution, which you can download as a PDF from the Sartorius Pipetting workflow editor.



## Example Workflows

New workflow templates are published regularly, and you can always create custom workflows to meet your specific needs. Available workflows for common laboratory tasks include the following:

- Serial dilution in plates and tubes
- Various reformatting workflows
- Media exchange and compound addition workflows

## Ideal for Repetitive or Complex Tasks

The Sartorius Pipetting software is perfect for repetitive pipetting workflows, such as those used in compound screening and quality control (QC). It is also ideal for long and tedious workflows where the potential for errors is high. Using the software will help your lab eliminate user-to-user variance, ensuring reliable and reproducible results.

## Download Now at No Cost

Experience the incredible Sartorius Pipetting software at no cost and streamline your laboratory operations today! Download the Sartorius Pipetting mobile app from the App Store or Play Store and access the Sartorius Pipetting workflow editor at <https://workflows.pipetting.sartorius.com>.

