

# SMALL YET MIGHTY



**Provectus™ 6500**  
Ultra Microbalance

  
**PerkinElmer®**  
*For the Better*

# THE WEIGHT OF ACCURACY

Weighing the tiniest samples with minimal error is all about balance. When it comes to accurately weighing the smallest amounts of your valuable samples, you need an efficient, error-free process that saves you time, money, and materials. The **Provectus 6500 ultra microbalance** ensures your sample preparation will be precise from the start, saving you from time-consuming errors down the road.

It is not the readability that determines the accuracy of a weighing instrument, but it's the repeatability or minimum weight capability. We know samples are the core of every analysis, and the Provectus 6500 ultra microbalance makes it possible to weigh extremely small quantities of your precious samples with ease and high accuracy.



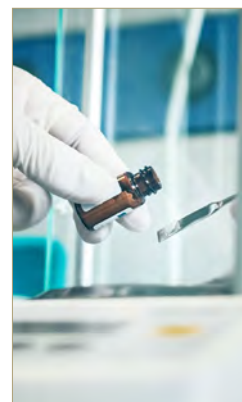
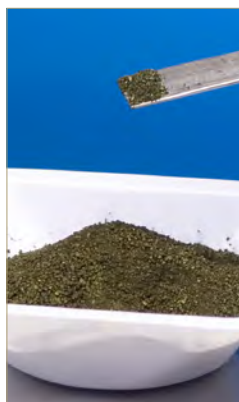
**Provectus 6500**  
Ultra Microbalance

INTRODUCTION

# ANALYZE RIGHT FROM THE START

Whether you're weighing insect wings or preparing samples and standards for major analytical techniques, our Provectus 6500 ultra microbalance is perfect for a wide range of applications. Applications include, but are not limited to, use in production areas for calibrating precision micropipettes, weighing precious metals and food, dispensing modern drugs, and preparing costly instrument standards in the environmental and biotechnology areas.

**Click on each application for more information.**



**Provectus 6500**  
Ultra Microbalance

APPLICATIONS

ORGANIC ELEMENTAL AND NITROGEN ANALYSIS

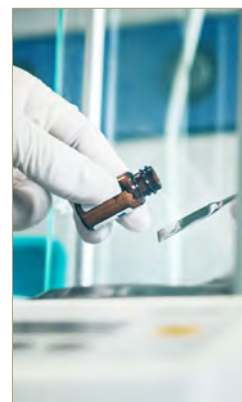
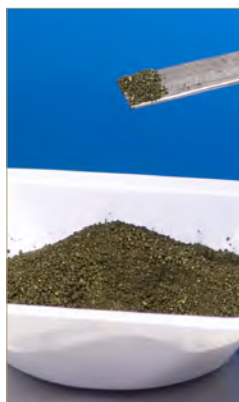
## Organic Elemental and Nitrogen Analysis

Simplify your application and minimize transcription errors when you connect the Provectus 6500 to our 2400 Series II CHNS/O Elemental Analyzer for an automatic transfer of sample weights.

# ANALYZE RIGHT FROM THE START

Whether you're weighing insect wings or preparing samples and standards for major analytical techniques, our Provectus 6500 ultra microbalance is perfect for a wide range of applications. Applications include, but are not limited to, use in production areas for calibrating precision micropipettes, weighing precious metals and food, dispensing modern drugs, and preparing costly instrument standards in the environmental and biotechnology areas.

**Click on each application for more information.**



## THERMAL ANALYSIS

### Thermal Analysis

In differential scanning calorimetry (DSC) and thermogravimetric analysis (TGA), where samples ranging in weight from less than a milligram to several hundred milligrams are common, weighing accuracy and precision are of paramount importance. The Provectus 6500 ultra microbalance allows automatic and accurate weight measurements beyond this weight range.

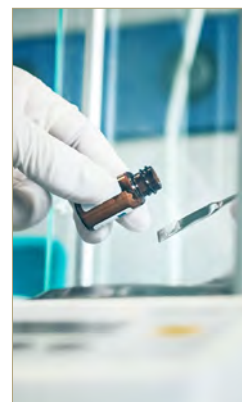
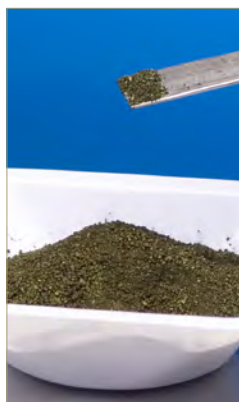
**Provectus 6500**  
Ultra Microbalance

## APPLICATIONS

# ANALYZE RIGHT FROM THE START

Whether you're weighing insect wings or preparing samples and standards for major analytical techniques, our Provectus 6500 ultra microbalance is perfect for a wide range of applications. Applications include, but are not limited to, use in production areas for calibrating precision micropipettes, weighing precious metals and food, dispensing modern drugs, and preparing costly instrument standards in the environmental and biotechnology areas.

**Click on each application for more information.**



BIOTECHNOLOGY/PHARMACEUTICALS

## Biotechnology/Pharmaceuticals

Accurate weighing of small particles reduces the cost of expensive or quantity-limited samples for analysis, and the Provectus 6500 ultra microbalance delivers.

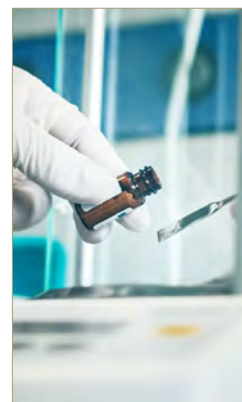
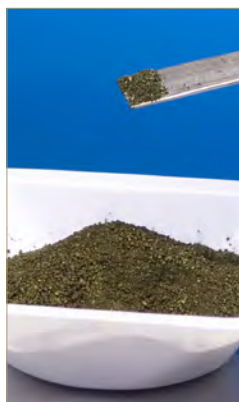
**Provectus 6500**  
Ultra Microbalance

APPLICATIONS

# ANALYZE RIGHT FROM THE START

Whether you're weighing insect wings or preparing samples and standards for major analytical techniques, our Provectus 6500 ultra microbalance is perfect for a wide range of applications. Applications include, but are not limited to, use in production areas for calibrating precision micropipettes, weighing precious metals and food, dispensing modern drugs, and preparing costly instrument standards in the environmental and biotechnology areas.

**Click on each application for more information.**



**Provectus 6500**  
Ultra Microbalance

APPLICATIONS

ENVIRONMENTAL

## Environmental

When weighing airborne or waterborne contaminants, it's common to collect samples onto a filter, weighing the filter before and after collection. The Provectus 6500 ultra microbalance has an optional filter stirrup to simplify routine weighing of filters.

# EQUAL BALANCE BETWEEN FORM AND FUNCTION



For industrial companies, universities, and pharmaceutical labs, accurate and precise measurements are vital to overall results. The Provectus 6500 is a compact, high-precision microprocessor-controlled ultra microbalance capable of weighing samples in the range of 0 mg to 5000 mg to an ultimate resolution of 0.1  $\mu\text{g}$ . This allows you to measure even the smallest weight changes.

With the Provectus 6500, you can weigh your materials directly in the sampling pans, avoiding sample transfer and possible loss of materials caused by using weighing papers. Often sold with our CHNS/O elemental analyzer, the ultra microbalance can be conveniently connected, automatically transferring weight measurements to the analyzer.

**Provectus 6500**  
Ultra Microbalance

TECHNOLOGY

# PRECISE ACCURATE PROVECTUS

Ultra microbalances are crucial to your research. You need precise measurements for the most accurate results, and the Provectus 6500 is up to the task. With benefits such as automation, data transfer, and overall convenience, you can be confident in your results from the first step to the last.

**Provectus 6500**  
Ultra Microbalance

TECHNOLOGY

## Automation

The easy-to-use, single-point automated calibration routine, automated single-keystroke tare, and range selection procedures simplify your process and improve performance.

## Data Transfer

Connect to the CHNS/O elemental analyzer for direct weight readings, or link to computers and other devices for easy record keeping and data sharing. Replace time-consuming, error-prone manual data capture and transfer process with automatic recording of your weight data for use in subsequent calculations or steps later in the process.

## Convenience

The weighing chamber is separate from the controller, enabling you to place the chamber in a comfortable, optimal position. Remote weighing is also possible, as the balance chamber and control units are separate. The controller can be conveniently placed outside your safety cabinet.

# MINIATURE SIZE, MAXIMUM PERFORMANCE

From its diminutive stature to its pinpoint accuracy to its user-friendly design, the Provectus 6500 is a must for your lab. See what makes this ultra microbalance ultra smart.

## ACCURATE MEASURING

The glass door minimizes air currents on the sample.

## SMALL FOOTPRINT

Stackable design saves valuable bench space in your lab.

## SLEEK STYLE

It features a modern, compact design.



## VERSATILE

Separate balance and control panel allow remote weighing inside the fume hood or dry box.

## CONVENIENT

One of the easiest microbalances to operate, it includes:

- Self-leveling tripod base
- Easy-glide door with cushioned glass door stoppers
- Solid back panel for improved depth perception
- Easy stirrup handling with newly designed push button
- Easy left- or right-hand measurements

**Provectus 6500**  
Ultra Microbalance

SPECIFICATIONS

# SMALL DETAILS, BIG DIFFERENCES

With a range of accessories – from filler paper stirrups to microspatulas – the Provectus 6500 ultra microbalance is as versatile as your lab. Start with the accessories you need now and add more as you need them.

## INCLUDED ACCESSORIES

Standard stirrup kit
Tare weight set (1 mg to 1 g)
Calibration weight
Nonmagnetic forceps
Instruction manual
CHNS/O Interface Cable
Foil pans (20 per pack)

## ADDITIONAL ACCESSORIES

Filter paper stirrups
Lightweight stirrups
Antistatic forceps kits
Thick-handled microspatula
Specifically shaped tweezers
Flat-bottomed weighing dishes

**Provectus 6500**  
Ultra Microbalance

ACCESSORIES

Visit us online at [www.perkinelmer.com](http://www.perkinelmer.com) to learn more about our solutions.

**PerkinElmer, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.