

## Weights and Weight Sets



# Table of Contents

From Weights to Certified Testing Services	4-5
Your DAkkS Partner for Mass	6-7
Reference Line Weights	8-11
Reference Line Weight Sets	12-13
Proof Line Weights	14-17
Proof Line Weight Sets	18-19
Special and Service Weight Sets	20-21
Cylindrical Weights	22-23
Accessories	24-25
Weight Set Combinations	28





# From Weights to Certified Testing Services

## The Complete Line

Regular inspection and testing of weighing instruments are essential to ensure reliable weighing results. Sartorius offers highly accurate metrological weights and weight sets with nominal mass values from 1 mg to 1,000 kg, special and test weights, as well as the right accessories required for correct handling, storage and cleaning of weights.

For unique identification and classification, all Sartorius weights – except for wire and leaf weights – are laser-engraved with an individual three-character alphanumeric marking in accordance with the OIML R111-1 International Recommendation. Therefore, Sartorius weights are 100% traceable throughout their entire life cycles. On weights made of cast iron, this unique marking is stamped on the cover plate.



# The Perfect Test Weight for Your Weighing Instrument

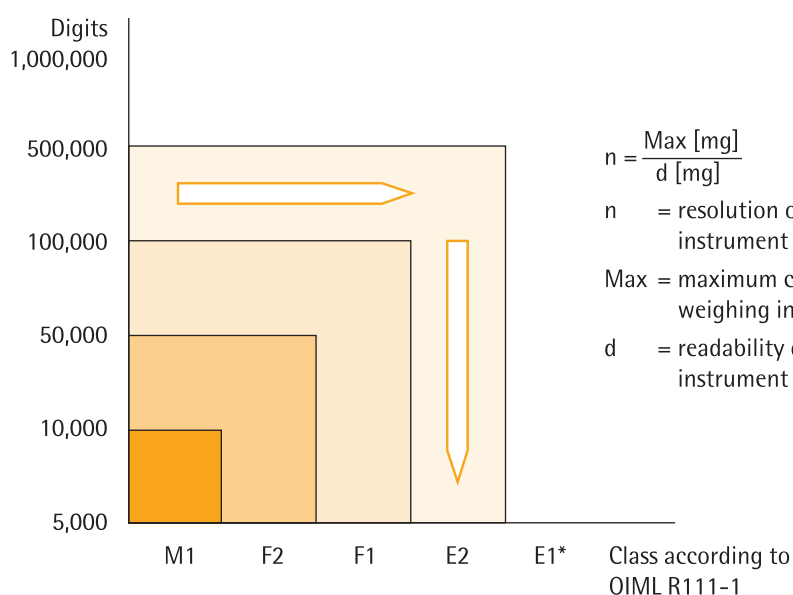
## Here's how to find the right weight

Just determine the number of digits specified for your weighing instrument's resolution, then check the graph below for the particular accuracy class that your weight must have.

The nominal mass value of your weight should be more than 80% of the maximum capacity of your weighing instrument.

Use the following chart to determine whether you need an individual weight or a weight set by comparing the nominal mass values.

Example: Suppose your weighing instrument has a maximum capacity of 2,200 g and a readability of 0.01 g. This yields 220,000 digits, which correspond to a class E2 weight. Select 2,000 g as the nominal mass value.



$$n = \frac{\text{Max [mg]}}{d \text{ [mg]}}$$

n = resolution of your weighing instrument (in digits)

Max = maximum capacity of your weighing instrument

d = readability of your weighing instrument

\* or E2, DAkkS-calibrated





# Your DAkkS Partner for Mass

Sartorius has calibration laboratories for weights and weighing instruments. These calibration laboratories are accredited by the DAkkS\* and meet the International Standard ISO IEC 17025, which is also recognized throughout Europe as a European Standard (EN ISO IEC 17025). Thus, weights calibrated by Sartorius are traceable to your national standard.

DAkkS assessments are made according to the applicable requirements and the corresponding rules of the European Co-operation for Accreditation (EA), the International Accreditation Forum (IAF) and the International Laboratory Accreditation Cooperation (ILAC).

\* German accreditation body for calibration, testing, inspection, certification of quality, products and services, etc.





**Recalibration of all weights,  
manufacturers and designs**

Depending on how frequently weights are used and which accuracy is required, they must be recalibrated at regular intervals so that they meet the requirements for reliable measuring, inspection and test equipment.

Sartorius offers recalibration service with DAkkS calibration certificates for all weights, regardless of class, design or brand.



Your local Sartorius Service Center will provide you with information on our recalibration service.

**Please send your weights to:**

Sartorius Lab Instruments GmbH & Co. KG  
DAkkS Calibration Laboratory  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

# Reference Line Weights

Sartorius Reference Line weights are the first choice for primary mass standards. They are supplied with a first-class finish featuring high gloss polished and corrosion-resistant surfaces. The quality of these weights exceeds the requirements of the OIML R111-1 International Recommendation. Sartorius Reference Line weights have only positive deviations from the nominal value\*.

OIML class E1 and E2 weights from 1 mg to 500 mg are available as wire weights; those with a nominal mass of 1 g and higher are designed as knob weights. They are manufactured from special stainless steel that has a density of  $8,000 \text{ kg/m}^3$  and comply with OIML R111-1 tolerances. The weights are non-magnetizable and are delivered in a high-quality wooden box\*\*.

\* Valid at the time of delivery and without taking into account the measurement uncertainty

\*\* Unless otherwise specified







Nominal Value	Without DAkkS calibration certificate		With DAkkS calibration certificate	
	OIML Class E1	OIML Class E2	OIML Class E1	OIML Class E2
1 mg	YCW0111-00	YCW0121-00	YCW0111-02	YCW0121-02
2 mg	YCW0211-00	YCW0221-00	YCW0211-02	YCW0221-02
5 mg	YCW0511-00	YCW0521-00	YCW0511-02	YCW0521-02
10 mg	YCW1111-00	YCW1221-00	YCW1111-02	YCW1121-02
20 mg	YCW1211-00	YCW1121-00	YCW1211-02	YCW1221-02
50 mg	YCW1511-00	YCW1521-00	YCW1511-02	YCW1521-02
100 mg	YCW2111-00	YCW2121-00	YCW2111-02	YCW2121-02
200 mg	YCW2211-00	YCW2221-00	YCW2211-02	YCW2221-02
500 mg	YCW2511-00	YCW2521-00	YCW2511-02	YCW2521-02
1 g	YCW3111-00	YCW3121-00	YCW3111-02	YCW3121-02
2 g	YCW3211-00	YCW3221-00	YCW3211-02	YCW3221-02
5 g	YCW3511-00	YCW3521-00	YCW3511-02	YCW3521-02
10 g	YCW4111-00	YCW4121-00	YCW4111-02	YCW4121-02
20 g	YCW4211-00	YCW4221-00	YCW4211-02	YCW4221-02
50 g	YCW4511-00	YCW4521-00	YCW4511-02	YCW4521-02
100 g	YCW5111-00	YCW5121-00	YCW5111-02	YCW5121-02
200 g	YCW5211-00	YCW5221-00	YCW5211-02	YCW5221-02
500 g	YCW5511-00	YCW5521-00	YCW5511-02	YCW5521-02
1 kg	YCW6111-00	YCW6121-00	YCW6111-02	YCW6121-02
2 kg	YCW6211-00	YCW6221-00	YCW6211-02	YCW6221-02
5 kg	YCW6511-00	YCW6521-00	YCW6511-02	YCW6521-02
10 kg	YCW7111-00	YCW7121-00	YCW7111-02	YCW7121-02
20 kg	YCW7211-00	YCW7221-00	YCW7211-02	YCW7221-02
50 kg	YCW7511-00	YCW7521-00	YCW7511-02	YCW7521-02

# Reference Line Weights

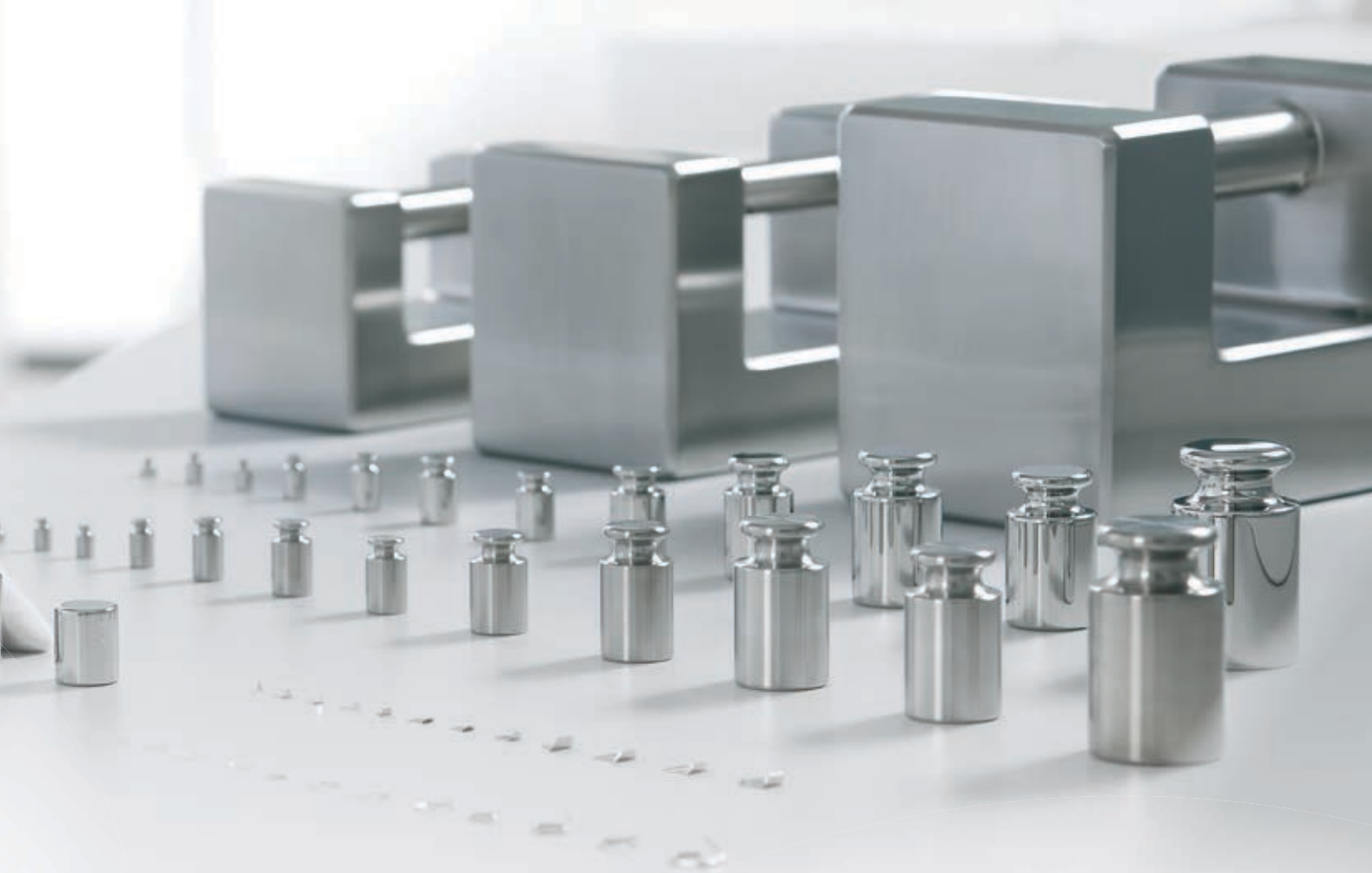
Sartorius Reference Line milligram weights of OIML class F1 are designed as wire weights. These weights are made of stainless steel that has a density of 7,950 kg/m<sup>3</sup>. The tolerances comply with OIML R111-1.

OIML class F1 and F2 weights with nominal values of 1 g and higher are available as knob weights and made of special stainless steel that has a density of 7,950 kg/m<sup>3</sup>. Their tolerances likewise comply with OIML R111-1.

The weights are non-magnetized and, up to 20 g, come in a plastic case. Starting at 50 g, they are supplied in a high-quality wooden case\*.

\* Unless otherwise specified

Nominal Value	Without DAkkS calibration certificate		With DAkkS calibration certificate	
	OIML Class F1	OIML Class F2	OIML Class F1	OIML Class F2
1 mg	YCW0131-00	-	YCW0131-02	-
2 mg	YCW0231-00	-	YCW0231-02	-
5 mg	YCW0531-00	-	YCW0531-02	-
10 mg	YCW1131-00	-	YCW1131-02	-
20 mg	YCW1231-00	-	YCW1231-02	-
50 mg	YCW1531-00	-	YCW1531-02	-
100 mg	YCW2131-00	-	YCW2131-02	-
200 mg	YCW2231-00	-	YCW2231-02	-
500 mg	YCW2531-00	-	YCW2531-02	-
1 g	YCW3131-00	YCW314-00	YCW313-02	YCW314-02
2 g	YCW323-00	YCW324-00	YCW323-02	YCW324-02
5 g	YCW353-00	YCW354-00	YCW353-02	YCW354-02
10 g	YCW413-00	YCW414-00	YCW413-02	YCW414-02
20 g	YCW423-00	YCW424-00	YCW423-02	YCW424-02
50 g	YCW453-00	YCW454-00	YCW453-02	YCW454-02
100 g	YCW513-00	YCW514-00	YCW513-02	YCW514-02
200 g	YCW523-00	YCW524-00	YCW523-02	YCW524-02
500 g	YCW553-00	YCW554-00	YCW553-02	YCW554-02
1 kg	YCW613-00	YCW614-00	YCW613-02	YCW614-02
2 kg	YCW623-00	YCW624-00	YCW623-02	YCW624-02
5 kg	YCW653-00	YCW654-00	YCW653-02	YCW654-02
10 kg	YCW713-00	YCW714-00	YCW713-02	YCW714-02
20 kg	YCW723-00	YCW724-00	YCW723-02	YCW724-02
50 kg	YCW753-00	YCW754-00	YCW753-02	YCW754-02



OIML class F1 and F2 cylindrical weights with a lug are made of a special stainless steel that has a density of 7,950 kg/m<sup>3</sup> and their tolerances comply with OIML R111-1. The weights are delivered in a wooden case suitable for transportation by a pallet mover.

Nominal Value	Without DAkkS calibration certificate		With DAkkS calibration certificate	
	OIML Class F1	OIML Class F2	OIML Class F1	OIML Class F2
100 kg	YCW813-00	YCW814-00	YCW813-02	YCW814-02
200 kg	YCW823-00	YCW824-00	YCW823-02	YCW824-02
500 kg	YCW853-00	YCW854-00	YCW853-02	YCW854-02
1,000 kg	YCW913-00	YCW914-00	YCW913-02	YCW914-02



# Reference Line Weight Sets

The weights of the Sartorius Reference Line weight sets\* have the same features and properties as the individual weights of this series, along with the corresponding maximum permissible errors. These weight sets are supplied with gloves, tweezers and a brush in a high-quality wooden case.

OIML class E1 and E2 weights in the Reference Line sets are made of special stainless steel that has a density of  $8,000 \text{ kg/m}^3$ , and their tolerances comply with OIML R111-1. Nominal mass values up to 500 mg feature a wire weight shape.

OIML class F1 and F2 weights of Reference Line sets are made of special stainless steel that has a density of  $7,950 \text{ kg/m}^3$ . Their tolerances comply with OIML R111-1.

OIML class F1 Reference Line weight sets contain wire weights with nominal values of up to 500 mg; those of OIML class F2 have leaf weights instead and nominal values of up to 500 mg.

Leaf weights from 1 mg to 5 mg are made of aluminum that has a density of  $2,650 \text{ kg/m}^3$ ; those with weights from 10 mg to 500 mg are made of nickel silver with a density of  $8,600 \text{ kg/m}^3$ . Their tolerances comply with OIML R111-1.





	Without DAkkS calibration certificate		With DAkkS calibration certificate	
Nominal Value	OIML Class E1	OIML Class E2	OIML Class E1	OIML Class E2
1 mg - 5 g	YCS011-351-00	YCS011-352-00	YCS011-351-02	YCS011-352-02
1 mg - 100 g	YCS011-511-00	YCS011-512-00	YCS011-511-02	YCS011-512-02
1 mg - 200 g	YCS011-521-00	YCS011-522-00	YCS011-521-02	YCS011-522-02
1 mg - 1 kg	YCS011-611-00	YCS011-612-00	YCS011-611-02	YCS011-612-02
1 mg - 5 kg	YCS011-651-00	YCS011-652-00	YCS011-651-02	YCS011-652-02
1 g - 1 kg	YCS31-611-00	YCS31-612-00	YCS31-611-02	YCS31-612-02
1 g - 5 kg	YCS31-651-00	YCS31-652-00	YCS31-651-02	YCS31-652-02
1 g - 10 kg	YCS31-711-00	YCS31-712-00	YCS31-711-02	YCS31-712-02

	Without DAkkS calibration certificate		With DAkkS calibration certificate	
Nominal Value	OIML Class F1	OIML Class F2	OIML Class F1	OIML Class F2
1 mg - 100 g	YCS01-513-00	YCS01-514-00	YCS01-513-02	YCS01-514-02
1 mg - 200 g	YCS01-523-00	YCS01-524-00	YCS01-523-02	YCS01-524-02
1 mg - 1 kg	YCS01-613-00	YCS01-614-00	YCS01-613-02	YCS01-614-02
1 mg - 5 kg	YCS01-653-00	YCS01-654-00	YCS01-653-02	YCS01-654-02
1 g - 1 kg	YCS31-613-00	YCS31-614-00	YCS31-613-02	YCS31-614-02
1 g - 5 kg	YCS31-653-00	YCS31-654-00	YCS31-653-02	YCS31-654-02
1 g - 10 kg	YCS31-713-00	YCS31-714-00	YCS31-713-02	YCS31-714-02

\* For a list of individual weight combinations in each set, see page 28.



# Proof Line Weights

Sartorius Proof Line weights are high-quality working standards for everyday use. These weights are made of high quality material with polished or precision-lathed surfaces and comply all respects with the requirements on shape, material and markings specified by the OIML R111-1 International Recommendation.

OIML class E2 milligram weights are designed as leaf weights. With a nominal mass ranging from 1 mg to 5 mg, these weights are made of aluminum that has a density of 2,650 kg/m<sup>3</sup>; those with nominal values ranging from 10 mg to 500 mg are manufactured from nickel silver with a density of 8,600 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

Weights with nominal values of 1 g and higher are available as knob weights and made of a special stainless steel that has a density of 7,950 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

Proof Line weights are non-magnetizable and supplied in a high-quality plastic case\*. These plastic cases are suitable for use in cleanrooms.

	Without DAkkS calibration certificate	With DAkkS calibration certificate
Nominal Value	OIML Class E2	
1 mg	YCW0121-AC-00	YCW0121-AC-02
2 mg	YCW0221-AC-00	YCW0221-AC-02
5 mg	YCW0521-AC-00	YCW0521-AC-02
10 mg	YCW1121-AC-00	YCW1121-AC-02
20 mg	YCW1221-AC-00	YCW1221-AC-02
50 mg	YCW1521-AC-00	YCW1521-AC-02
100 mg	YCW2121-AC-00	YCW2221-AC-02
200 mg	YCW2221-AC-00	YCW2121-AC-02
500 mg	YCW2521-AC-00	YCW2521-AC-02
1 g	YCW3121-AC-00	YCW3121-AC-02
2 g	YCW3221-AC-00	YCW3221-AC-02
5 g	YCW3521-AC-00	YCW3521-AC-02
10 g	YCW4121-AC-00	YCW4121-AC-02
20 g	YCW4221-AC-00	YCW4221-AC-02
50 g	YCW4521-AC-00	YCW4521-AC-02
100 g	YCW5121-AC-00	YCW5121-AC-02
200 g	YCW5221-AC-00	YCW5221-AC-02
500 g	YCW5521-AC-00	YCW5521-AC-02
1 kg	YCW6121-AC-00	YCW6121-AC-02
2 kg	YCW6221-AC-00	YCW6221-AC-02
5 kg	YCW6521-AC-00	YCW6521-AC-02
10 kg	YCW7121-AC-00	YCW7121-AC-02

\* Unless otherwise specified



Sartorius Proof Line knob weights of OIML classes F1 and F2 with nominal values of 1 g and higher are made of stainless steel that has a density of 7,950 kg/m<sup>3</sup>. The tolerances of these knob weights comply with OIML R111-1 and they are supplied in high-quality plastic screw-top boxes. Knob weights from 10 g and up feature an adjustment chamber.



Proof Line OIML class F1 leaf weights ranging from 1 mg to 5 mg are made out of aluminum that has a density of 2,650 kg/m<sup>3</sup>; those with nominal values from 10 mg to 500 mg are manufactured from nickel silver that has a density of 8,600 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

Milligram weights of the Proof Line range are delivered in a high-quality plastic case.

These plastic cases are suitable for use in cleanrooms.

Nominal Value	Without DAkkS calibration certificate		With DAkkS calibration certificate	
	OIML Class F1	OIML Class F2	OIML Class F1	OIML Class F2
1 mg	YCW013-00	-	YCW013-02	-
2 mg	YCW023-00	-	YCW023-02	-
5 mg	YCW053-00	-	YCW053-02	-
10 mg	YCW113-00	-	YCW113-02	-
20 mg	YCW123-00	-	YCW123-02	-
50 mg	YCW153-00	-	YCW153-02	-
100 mg	YCW213-00	-	YCW213-02	-
200 mg	YCW223-00	-	YCW223-02	-
500 mg	YCW253-00	-	YCW253-02	-
1 g	YCW313-AC-00	YCW314-AC-00	YCW313-AC-02	YCW314-AC-02
2 g	YCW323-AC-00	YCW324-AC-00	YCW323-AC-02	YCW324-AC-02
5 g	YCW353-AC-00	YCW354-AC-00	YCW353-AC-02	YCW354-AC-02
10 g	YCW413-AC-00	YCW414-AC-00	YCW413-AC-02	YCW414-AC-02
20 g	YCW423-AC-00	YCW424-AC-00	YCW423-AC-02	YCW424-AC-02
50 g	YCW453-AC-00	YCW454-AC-00	YCW453-AC-02	YCW454-AC-02
100 g	YCW513-AC-00	YCW514-AC-00	YCW513-AC-02	YCW514-AC-02
200 g	YCW523-AC-00	YCW524-AC-00	YCW523-AC-02	YCW524-AC-02
500 g	YCW553-AC-00	YCW554-AC-00	YCW553-AC-02	YCW554-AC-02
1 kg	YCW613-AC-00	YCW614-AC-00	YCW613-AC-02	YCW614-AC-02
2 kg	YCW623-AC-00	YCW624-AC-00	YCW623-AC-02	YCW624-AC-02
5 kg	YCW653-AC-00	YCW654-AC-00	YCW653-AC-02	YCW654-AC-02
10 kg	YCW713-AC-00	YCW714-AC-00	YCW713-AC-02	YCW714-AC-02

# Proof Line Weights

Sartorius Proof Line knob weights of OIML class M1 are made of stainless steel, precision-lathed and have a density of 7,850 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

OIML class M2 knob weights of the Sartorius Proof Line series are manufactured from brass, precision-lathed and have a density of 8,400 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1

All Proof Line weights are supplied in high-quality plastic screw-top boxes that are





suitable for use in cleanrooms. Knob weights from 10 g and up feature an adjustment chamber.

OIML M1 and M2 block weights and ton weights of the Proof Line series are made of cast iron that is black lacquered and has a density of 7,200 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

Proof Line OIML class M1 block weights are also available in a choice of stainless steel that has a density of 7,850 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

Nominal Value	Without DAkkS calibration certificate		With DAkkS calibration certificate	
	OIML Class M1	OIML Class M2	OIML Class M1	OIML Class M2
1 g	YCW315-AC-00	YCW316-AC-00	YCW315-AC-02	YCW316-AC-02
2 g	YCW325-AC-00	YCW326-AC-00	YCW325-AC-02	YCW326-AC-02
5 g	YCW355-AC-00	YCW356-AC-00	YCW355-AC-02	YCW356-AC-02
10 g	YCW415-AC-00	YCW416-AC-00	YCW415-AC-02	YCW416-AC-02
20 g	YCW425-AC-00	YCW426-AC-00	YCW425-AC-02	YCW426-AC-02
50 g	YCW455-AC-00	YCW456-AC-00	YCW455-AC-02	YCW456-AC-02
100 g	YCW515-AC-00	YCW516-AC-00	YCW515-AC-02	YCW516-AC-02
200 g	YCW525-AC-00	YCW526-AC-00	YCW525-AC-02	YCW526-AC-02
500 g	YCW555-AC-00	YCW556-AC-00	YCW555-AC-02	YCW556-AC-02
1 kg	YCW615-AC-00	YCW616-AC-00	YCW615-AC-02	YCW616-AC-02
2 kg	YCW625-AC-00	YCW626-AC-00	YCW625-AC-02	YCW626-AC-02
5 kg	YCW655-AC-00	YCW656-AC-00	YCW655-AC-02	YCW656-AC-02
10 kg	YCW715-AC-00	YCW716-AC-00	YCW715-AC-02	YCW716-AC-02



		Without DAkkS calibration certificate		With DAkkS calibration certificate	
Nominal Value		OIML Class M1	OIML Class M2	OIML Class M1	OIML Class M2
<b>Block weights, cast iron</b>					
	5 kg	YCW6559-AC-00	YCW6569-AC-00	YCW6559-AC-02	YCW6569-AC-02
	10 kg	YCW7159-AC-00	YCW7169-AC-00	YCW7159-AC-02	YCW7169-AC-02
	20 kg	YCW7259-AC-00	YCW7269-AC-00	YCW7259-AC-02	YCW7269-AC-02
	50 kg	YCW7559-AC-00	YCW7569-AC-00	YCW7559-AC-02	YCW7569-AC-02
<b>Block weights, stainless steel</b>					
	5 kg	YCW6554-AC-00	-	YCW6554-AC-02	-
	10 kg	YCW7154-AC-00	-	YCW7154-AC-02	-
	20 kg	YCW7254-AC-00	-	YCW7254-AC-02	-
	50 kg	YCW7554-AC-00	-	YCW7554-AC-02	-
<b>Block weights, cast iron</b>					
	100 kg	YCW8159-AC-00	-	YCW8159-AC-02	-
	200 kg	YCW8259-AC-00	-	YCW8259-AC-02	-
	500 kg	YCW8559-AC-00	-	YCW8559-AC-02	-
	1,000 kg	YCW9159-AC-00	-	YCW9159-AC-02	-
<b>Ton weights, cast iron</b>					
	100 kg	YCW8157-AC-00	-	YCW8157-AC-02	-
	200 kg	YCW8257-AC-00	-	YCW8257-AC-02	-
	500 kg	YCW8557-AC-00	-	YCW8557-AC-02	-
	1,000 kg	YCW9157-AC-00	-	YCW9157-AC-02	-





# Proof Line Weight Sets

The weights of the Sartorius Proof Line weight sets\* have the same features and properties as the individual weights of this series, along with the corresponding maximum permissible errors.

These weight sets are supplied with gloves, tweezers and a brush in a high-quality plastic case.

Proof Line milligram weights are provided as leaf weights. With a nominal mass ranging from 1 mg to 5 mg, they are made of aluminum with a density of  $2,650 \text{ kg/m}^3$ ; those with nominal values of 10 mg to 500 mg are made of nickel silver that has a density of  $8,600 \text{ kg/m}^3$ . Their tolerances comply with OIML R111-1.

\* For a list of individual weight combinations in each set, see page 28.





Without DAkkS calibration certificate

Nominal Value	OIML Class E2	OIML Class F1	OIML Class F2	OIML Class M1
1 mg - 500 mg	YCS011-252-AC-00	YCS01-253-AC-00	YCS01-254-AC-00	YCS01-255-AC-00
1 mg - 100 g	-	-	-	YCS01-515-AC-00
1 mg - 200 g	-	-	-	YCS01-525-AC-00
1 mg - 1 kg	-	-	-	YCS01-615-AC-00
1 mg - 5 kg	-	-	-	YCS01-655-AC-00
1 g - 100 g	YCS31-512-AC-00	YCS31-513-AC-00	YCS31-514-AC-00	-
1 g - 200 g	YCS31-522-AC-00	-	-	-
1 g - 1 kg	YCS31-612-AC-00	YCS31-613-AC-00	YCS31-614-AC-00	YCS31-615-AC-00
1 g - 5 kg	YCS31-652-AC-00	YCS31-653-AC-00	YCS31-654-AC-00	YCS31-655-AC-00
1 g - 10 kg	-	-	-	YCS31-715-AC-00

With DAkkS calibration certificate

Nominal Value	OIML Class E2	OIML Class F1	OIML Class F2	OIML Class M1
1 mg - 500 mg	YCS011-252-AC-02	YCS01-253-AC-02	YCS01-254-AC-02	YCS01-255-AC-02
1 mg - 100 g	-	-	-	YCS01-515-AC-02
1 mg - 200 g	-	-	-	YCS01-525-AC-02
1 mg - 1 kg	-	-	-	YCS01-615-AC-02
1 mg - 5 kg	-	-	-	YCS01-655-AC-02
1 g - 100 g	YCS31-512-AC-02	YCS31-513-AC-02	YCS31-514-AC-02	-
1 g - 200 g	YCS31-522-AC-02	-	-	-
1 g - 1 kg	YCS31-612-AC-02	YCS31-613-AC-02	YCS31-614-AC-02	YCS31-615-AC-02
1 g - 5 kg	YCS31-652-AC-02	YCS31-653-AC-02	YCS31-654-AC-02	YCS31-655-AC-02
1 g - 10 kg	-	-	-	YCS31-715-AC-02



# Special Weight Sets and Service Weight Sets



All weights of the Sartorius service weight sets\* have the same features and properties as the individual weights of the Sartorius Reference Line series, along with the corresponding maximum permissible errors. These service weight sets are supplied with gloves, tweezers and a brush in a service case.

\* For a list of individual weight combinations in each set, see page 28.

## Service Weight Sets

OIML Class	Nominal Value	Without DAkkS calibration certificate	With DAkkS calibration certificate
		Order No.	Order No.
E2	1 mg - 5 kg	YSS0121-353-00	YSS0121-353-02
E2	100 g - 5 kg	YSS5128-6528-00	YSS5128-6528-02
F1	1 g - 5 kg	YSS3138-6538-00	YSS3138-6538-02



The Sartorius density reference set\* consists of stackable disc weights. All weights are made of special stainless steel that has a density of  $8,000 \text{ kg/m}^3$ . Their tolerances comply with OIML R111-1. Density reference weights are non-magnetizable and are supplied in a high-quality wooden case with a calibration certificate, including density determination data.

\* For weight set compositions, see page 28



### Density Reference Set

OIML Class	Nominal Value	Order No.
E1	1 g - 1 kg	YCS31-612-09

# Cylindrical Weights

Sartorius cylindrical weights of OIML classes E2, F1 and F2 are made of special stainless steel that has a density of 7,950 kg/m<sup>3</sup>. Their tolerances comply with OIML R111-1.

These weights are non-magnetizable and have a polished surface. They are supplied in high-quality plastic cases that are suitable for use in cleanrooms.

Without DAkkS calibration certificate

Nominal Value	OIML Class E2	OIML Class F1	OIML Class F2
1 g	YCW3128-00	YCW3138-00	-
2 g	YCW3228-00	YCW3238-00	-
5 g	YCW3528-00	YCW3538-00	-
10 g	YCW4128-00	YCW4138-00	-
20 g	YCW4228-00	YCW4238-00	-
50 g	YCW4528-00	YCW4538-00	-
100 g	YCW5128-00	YCW5138-00	YCW5148-00
200 g	YCW5228-00	YCW5238-00	YCW5248-00
500 g	YCW5528-00	YCW5538-00	YCW5548-00
1 kg	YCW6128-00	YCW6138-00	YCW6148-00
2 kg	YCW6228-00	YCW6238-00	YCW6248-00
5 kg	YCW6528-00	YCW6538-00	YCW6548-00
10 kg	-	YCW7138-00	YCW7148-00







With DAkkS calibration certificate

Nominal Value	OIML Class E2	OIML Class F1	OIML Class F2
1 g	YCW3128-02	YCW3138-02	-
2 g	YCW3228-02	YCW3238-02	-
5 g	YCW3528-02	YCW3538-02	-
10 g	YCW4128-02	YCW4138-02	-
20 g	YCW4228-02	YCW4238-02	-
50 g	YCW4528-02	YCW4538-02	-
100 g	YCW5128-02	YCW5138-02	YCW5148-02
200 g	YCW5228-02	YCW5238-02	YCW5248-02
500 g	YCW5528-02	YCW5538-02	YCW5548-02
1 kg	YCW6128-02	YCW6138-02	YCW6148-02
2 kg	YCW6228-02	YCW6238-02	YCW6248-02
5 kg	YCW6528-02	YCW6538-02	YCW6548-02
10 kg	-	YCW7138-02	YCW7148-02



# Accessories for Cleaning and Handling of Weights

According to the OIML R111-1 International Recommendation, special care must be taken in handling weights. Sartorius supplies a wide range of work-saving accessories to enable you to meet such special requirements.

To ensure correct handling and cleaning of weights, Sartorius offers brushes, gloves

and handling tools in various designs and sizes.

For proper storage of weights, Sartorius provides glass bell jars and plastic screw-top boxes in different sizes. Wood and plastic cases are available in a choice of designs on request.

Accessories		Order No.
Brush	Small, 100 mm	YAW11
	Medium, 115 mm	YAW12
	Large, 150 mm	YAW13
	Extra large, 250 mm	YAW14
Gloves	Cotton	YAW21
	Leather	YAW22
Forceps with silicone-coated tips	115 mm for 1 mg – 5 g	YAW31
	160 mm for 1 g – 200 g	YAW32
	230 mm for 1 g – 1 kg	YAW33
Tweezers with carbon tips	105 mm for 1 mg – 50 g	YAW35
	115 mm for 1 mg – 5 g	YAW36
	160 mm for 1 g – 200 g	YAW37
	230 mm for 1 g – 1 kg	YAW38
Weight forks	For 500 g	YAW41
	For 1 kg	YAW42
	For 2 kg	YAW43
Weight fork Fo(u)rk		
Fork, complete incl. adapters	For 500 g – 5 kg	YAW70
Fork	For 1 kg and 5 kg	YAW71
Adapter	For 500 g	YAW72
	For 2 kg	YAW73



# Accessories for Storage of Weights

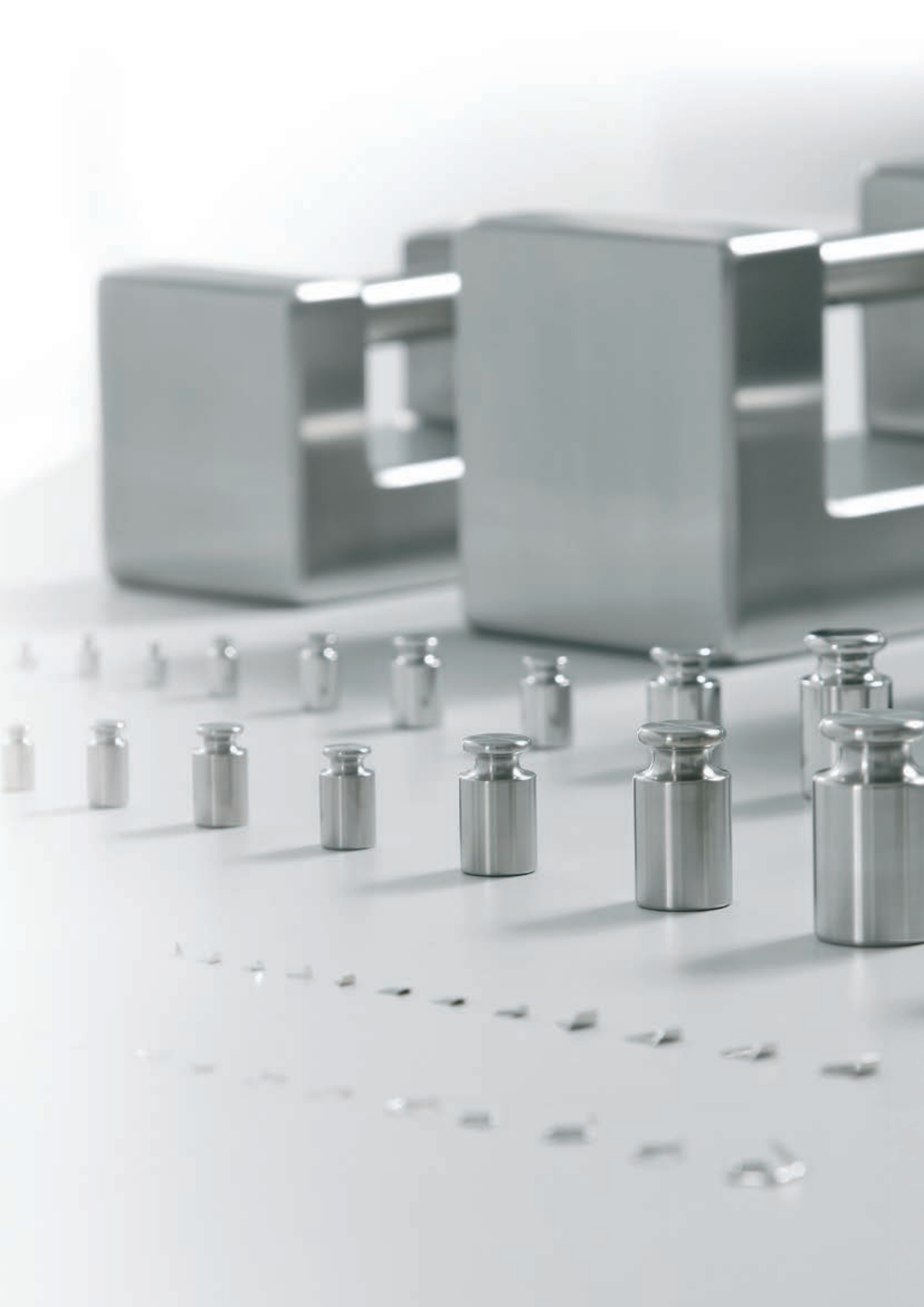
Accessories		Order No.
Handles for lifting weights	For 5 kg	YAW50
	For 10 kg	YAW51
	For 20 kg	YAW52
	For 50 kg	YAW53
Glass bell jar with support plate	For 1 mg - 5 g	YAW00
	For 1 mg - 50 g (100 g or 200 g)	YAW01
	For 100 g - 1 kg (2 kg)	YAW02
	For 2 kg - 5 kg	YAW03
	For 10 kg	YAW04
	For 20 kg	YAW05
	For 50 kg	YAW06
Plastic screw-top boxes with closed foam inlay are suitable for use in cleanrooms	For 1 mg - 100 mg weights	YAW01GL
	For 200 mg - 500 mg weights	YAW05GL
	For 1 g - 20 g weights*	YAW20GL
	For 50 g weights*	YAW50GL
	For 100 g weights*	YAW100GL
	For 200 g weights*	YAW200GL
	For 500 g weights*	YAW500GL
	For 1 kg weights*	YAW1000GL
	For 2 kg weights*	YAW2000GL
	For 5 kg weights*	YAW5000GL
	For 10 kg weights*	YAW10000GL

\* For knob weights only; for information on boxes for cylindrical weights, please contact Sartorius (equal price)



# Maximum permissible errors for weights according to OIML R111-1:2004

+/- in mg	Class								
Nominal Values	E1	E2	F1	F2	M1	M1-2	M2	M2-3	M3
1 mg	0.003	0.006	0.02	0.06	0.2				
2 mg	0.003	0.006	0.02	0.06	0.2				
5 mg	0.003	0.006	0.02	0.06	0.2				
10 mg	0.003	0.008	0.025	0.08	0.25				
20 mg	0.003	0.010	0.03	0.10	0.3				
50 mg	0.004	0.012	0.04	0.12	0.4				
100 mg	0.005	0.016	0.05	0.16	0.5		1.6		
200 mg	0.006	0.020	0.06	0.20	0.6		2.0		
500 mg	0.008	0.025	0.08	0.25	0.8		2.5		
1 g	0.010	0.030	0.10	0.3	1.0		3		10
2 g	0.012	0.040	0.12	0.4	1.2		4		12
5 g	0.016	0.050	0.16	0.5	1.6		5		16
10 g	0.020	0.060	0.20	0.6	2.0		6		20
20 g	0.025	0.080	0.25	0.8	2.5		8		25
50 g	0.030	0.10	0.30	1.0	3.0		10		30
100 g	0.05	0.16	0.5	1.6	5		16		50
200 g	0.10	0.30	1.0	3.0	10		30		100
500 g	0.25	0.8	2.5	8.0	25		80		250
1 kg	0.5	1.6	5	16	50		160		500
2 kg	1.0	3.0	10	30	100		300		1,000
5 kg	2.5	8.0	25	80	250		800		2,500
10 kg	5	16	50	160	500		1,600		5,000
20 kg	10	30	100	300	1,000		3,000		10,000
50 kg	25	80	250	800	2,500	5,000	8,000	16,000	25,000
100 kg		160	500	1,600	5,000	10,000	1,6000	30,000	50,000
200 kg		300	1,000	3,000	10,000	20,000	30,000	60,000	100,000
500 kg		800	2,500	8,000	25,000	50,000	80,000	160,000	250,000
1,000 kg		1,600	5,000	16,000	50,000	100,000	160,000	300,000	500,000
2,000 kg			10,000	30,000	100,000	200,000	300,000	600,000	1,000,000
5,000 kg			25,000	80,000	250,000	500,000	800,000	1,600,000	2,500,000



# Weight Set Combinations

Range	Contents	mg	g	kg	Range	Contents	mg	g	kg
<b>1 mg – 5 g</b>	1	●	●		<b>1 mg – 5 kg</b>	1	●	●	●
Total	2	●○	●○		Total	2	●○	●○	●○
contents:	5	●	●		contents:	5	●	●	●
11.11 g	10	●			11,111.11 g	10	●	●	
16 units	20	●○			28 units	20	●○	●○	
	50	●				50	●	●	
	100	●				100	●	●	
	200	●○				200	●○	●○	
	500	●				500	●	●	
<b>1 mg – 100 g</b>	1	●	●		<b>1 g – 1 kg</b>	1		●	●
Total	2	●○	●○		Total	2		●○	
contents:	5	●	●		contents:	5		●	
211.11 g	10	●	●		2,110 g	10		●	
21 units	20	●○	●○		13 units	20		●○	
	50	●	●			50		●	
	100	●	●			100		●	
	200	●○				200		●○	
	500	●				500		●	
<b>1 mg – 200 g</b>	1	●	●		<b>1 g – 5 kg</b>	1		●	●
Total	2	●○	●○		Total	2		●○	●○
contents:	5	●	●		contents:	5		●	●
611.11 g	10	●	●		11,110 g	10		●	
23 units	20	●○	●○		16 units	20		●○	
	50	●	●			50		●	
	100	●	●			100		●	
	200	●○	●○			200		●○	
	500	●				500		●	
<b>1 mg – 1 kg</b>	1	●	●	●	<b>1 g – 10 kg</b>	1		●	●
Total	2	●○	●○		Total	2		●○	●○
contents:	5	●	●		contents:	5		●	●
2,111.11 g	10	●	●		21,110 g	10		●	●
25 units	20	●○	●○		17 units	20		●○	
	50	●	●			50		●	
	100	●	●			100		●	
	200	●○	●○			200		●○	
	500	●	●			500		●	

● Weight without marking  
 ○ Weight with marking

\* 1 kg and up: gloves included







# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Weender Landstrasse 94–108  
37075 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan

Phone +33.1.70.62.50.00  
Fax +33.1.64.59.76.39

### Austria

Sartorius Austria GmbH  
Franzosengraben 12  
1030 Vienna

Phone +43.1.7965760.0  
Fax +43.1.7965760.24

### Belgium

Sartorius Belgium N.V.  
Leuvensesteenweg, 248/B  
1800 Vilvoorde

Phone +32.2.756.06.71  
Fax +32.2.253.45.95

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki

Phone +358.9.755.951  
Fax +358.9.755.95.200

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi

Phone +3623.457.227  
Fax +3623.457.147

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11

Phone +353.1.8089050  
Fax +353.1.8089388

### Italy

Sartorius Italy S.r.l.  
Viale A. Casati, 4  
20835 Muggiò (MB)

Phone +39.039.4659.1  
Fax +39.039.4659.88

### Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001  
Fax +31.30.60.52.917  
[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn

Phone +48.61.6473830  
Fax +48.61.6473839

### Russian Federation & Ukraine

LLC "Sartorius ICR"  
and LLC "Biohit"  
Uralskaya str. 4, Lit. B  
199155, Saint-Petersburg

Phone +7.812.327.5.327  
Fax +7.812.327.5.323

LLC "Biohit"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918  
Fax +380.50.623.3162

### Spain & Portugal

Sartorius Spain S.A.  
Offices in Madrid:  
C/ Isabel Colbrand, 10–12, of. 70  
28050 Madrid

Phone Spain +34.902.123.367  
Phone Portugal +351.800.855.800  
Fax Spain +34.91.358.84.85  
Fax Portugal +351.800.855.799

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00  
Fax +41.44.746.50.50

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ

Phone +44.1372.737159  
Fax +44.1372.726171

## America

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
Fax +1.631.254.4253

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 – SP- Brasil

Phone +55.11.4362.8900  
Fax + 55.11.4362.8901

### Canada

Sartorius Canada Inc.  
2179 Dunwin Drive #4  
Mississauga, ON L5L 1X2

Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

### Mexico

Sartorius de México S.A. de C.V.  
Circuito Circunvalación Poniente  
No. 149  
Ciudad Satélite  
53100, Estado de México  
México

Phone +52.5555.62.1102  
Fax +52.5555.62.2942

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7–11 Rodeo Drive  
Dandenong South Vic 3175

Phone +61.3.8762.1800  
Fax +61.3.8762.1828

### China

Sartorius Scientific  
Instruments (Beijing) Co., Ltd.  
33 Yu An Road, Airport Industrial Park  
Zone B,  
Shunyi District, Beijing 101300,  
P.R.China

Phone +86.10.8042.6300  
Fax +86.10.8042.6486

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwung Tong  
Kowloon, Hong Kong

Phone +852.2774.2678  
Fax +852.2766.3526

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India

Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### Japan

Sartorius Stedim Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300  
Fax +81.3.4331.4301

### Malaysia

Sartorius Malaysia Sdn. Bhd  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Singapore Pte. Ltd  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528

Phone +65.6872.3966  
Fax +65.6778.2494

### South Korea

Sartorius Korea Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, BunDang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700  
Fax +82.31.622.5799

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310

Phone +66.2643.8361-6  
Fax +66.2643.8367



◀ [www.sartorius.com](http://www.sartorius.com)