



CERTIFICATE OF ACCREDITATION

ANSI National Accreditation Board
11617 Coldwater Road, Fort Wayne, IN 46845 USA

This is to certify that

Sartorius Corporation
545-1 Johnson Avenue
Bohemia, NY 11716

has been assessed by ANAB and meets the requirements of international standard

ISO/IEC 17025:2017

while demonstrating technical competence in the field of

CALIBRATION

Refer to the accompanying Scope of Accreditation for information regarding the types of activities to which this accreditation applies

ACT-1252

Certificate Number



ANAB Approval

Certificate Valid Through: 12/29/2020
Version No. 015 Issued: 01/31/2019



This laboratory is accredited in accordance with the recognized International Standard ISO/IEC 17025:2017. This accreditation demonstrates technical competence for a defined scope and the operation of a laboratory quality management system (refer to joint ISO-ILAC-IAF Communiqué dated April 2017).



SCOPE OF ACCREDITATION TO ISO/IEC 17025:2017

Sartorius Corporation

545-1 Johnson Avenue
Bohemia, New York 11716
Ilhyun Joung
631-254-4249 ext 8469

CALIBRATION

Valid to: **December 29, 2020**

Certificate Number: **ACT-1252**

Mass and Mass Related

Parameter/Equipment	Range	Expanded Uncertainty of Measurement (+/-)	Reference Standard, Method, and/or Equipment
Ultra-Micro Balances ^{1,2}	Up to 2.1 g (0.000 1 mg)	2.7 µg + 2.5 µg/g	Euramet cg-18 and ASTM Class1 Weights
Micro Balances ^{1,2}	Up to 31 g (0.001 mg)	5.7 µg + 0.42 µg/g	
Semi-Micro Balances ^{1,2} / Speedcal	Up to 220 g (0.01 mg)	33 µg + 0.59 µg/g	
Analytical Balances ^{1,2}	Up to 520 g (0.000 1 g)	200 µg + 0.44 µg/g	
Precision Balances ^{1,2}	Up to 5 200 g (0.001 g)	1.8 mg + 0.000 42 mg/g	
Top Loading Balances ^{1,2}	Up to 14.2 kg (0.01 g)	18 mg + 0.001 4 mg/kg	
Top Loading Balances ^{1,2}	Up to 70.2 kg (0.1 g)	1.7 g + 0.002 5 mg/kg	
Top Loading Balances ^{1,2}	Up to 150 kg (0.1 g)	1.7 g + 1.2 g /kg	
Top Loading Balances ^{1,2}	Up to 70.2 kg (1 g)	2.4 g + 0.024 g/kg	



Mass and Mass Related

Parameter/Equipment	Range	Expanded Uncertainty of Measurement (+/-)	Reference Standard, Method, and/or Equipment
Moisture Analyzers ^{1,2} Balance	Up to 200 g (0.001 g) Up to 100 g (0.000 1 g)	1.8 mg 200 µg	Partial calibration using Euramet cg-18 and ASTM Class1 Weights
Volume ³ Pipettes, Liquid Handling Devices	Up to 10 µl (10 to 20) µl (20 to 50) µl (50 to 100) µl (100 to 200) µl (200 to 300) µl (300 to 500) µl (500 µl to 1) ml (1 to 2) ml (2 to 5) ml (5 to 10) ml	0.096 µl 0.29 % of reading 0.12 % of reading 0.06 % of reading 0.036 % of reading 0.028 % of reading 0.024 % of reading 0.021 % of reading 0.021 % of reading 0.020 % of reading 0.020 % of reading	ISO 8655-6 : Gravimetric method using Micro Balance, Semi-Micro Balance, Analytical Balances

Calibration and Measurement Capability (CMC) is expressed in terms of the measurement parameter, measurement range, expanded uncertainty of measurement and reference standard, method, and/or equipment. The expanded uncertainty of measurement is expressed as the standard uncertainty of the measurement multiplied by a coverage factor of 2 ($k=2$), corresponding to a confidence level of approximately 95%.

Notes:

1. On-site calibration service is available for this parameter, since on-site conditions are typically more variable than those in the laboratory, larger measurement uncertainties are expected on-site than what is reported on the accredited scope.
2. Range Number in parentheses is the instrument resolution.
3. Expanded uncertainty for volume in this Scope do not include contribution due to process repeatability, which will be included in the expanded uncertainty in the calibration certificates.
4. This scope is formatted as part of a single document including Certificate of Accreditation No. ACT-1252.



Vice President