



# CERTIFICATE OF ACCREDITATION

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

This is to certify that

**Sartorius Canada Inc.**  
**1173 North Service Road West, Unit D-4**  
**Oakville, ON L6M 2V9**

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-1252.01  
Certificate Number

  
ANAB Approval

Certificate Valid Through: 12/29/2020  
Version No. 005 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 Canada Inc.**

1173 North Service Road West, Unit D-4  
Oakville, ON L6M 2V9  
Ilhyun Joung  
631-254-4249 ext 8469

**CALIBRATION**

Valid to: **December 29, 2020**

Certificate Number: **ACT-1252.01**

**Mass and Mass Related**

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



Mass and Mass Related

Parameter/Equipment	Range <sup>2</sup>	Expanded Uncertainty of Measurement (+/-)	Reference Standard, Method, and/or Equipment
Moisture Analyzers <sup>1</sup>  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

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. This scope is formatted as part of a single document including Certificate of Accreditation No. ACT-1252.01.



Vice President

