

COPY

(English Translation)



Certificate of Accreditation

To Sartorius Japan K. K.

IAJapan hereby accredits the following laboratory as a calibration laboratory based on the Measurement Law as it meets the requirements of relevant international standards. This laboratory also meets the requirements for Mutual Recognition Arrangements (MRA) of ILAC and APLAC.

Accreditation No. JCSS0288

Name of Laboratory

Sartorius Japan K. K.

Service Department, Liquid Handling Service

Address of Laboratory

1-2-34 Ichigaya Sadohara-cho Shinjuku-ku Tokyo 162-0842,
Japan

Accreditation Scope

Volume (as attached)

Accreditation Criterion ISO/IEC 17025:2005

Date of Initial Accreditation: 2012-03-29

Latest Date of Issue: 2014-10-21

Ichiro Fujima

Chief Executive, IAJapan

National Institute of Technology and Evaluation

-
- International Accreditation Japan (IAJapan) is a laboratory accreditation body which has signed MRAs of ILAC(International Laboratory Accreditation Cooperation) and APLAC (Asia Pacific Laboratory Accreditation Cooperation).
 - MRA requirements are, in addition to relevant international standards and guides, requirements for participation in proficiency testing programmes, surveillance and reassessment, and the policy for the traceability of measurement for MRA purpose.
 - This accreditation demonstrates technical competence for a defined scope and the operation of a laboratory quality management system. The management system requirements in ISO/IEC 17025:2005 meet the principles of ISO 9001:2008 and are aligned with its pertinent requirements.

General Field of Calibration : Volume

Date of Initial Accreditation of the Field : 2012-03-29

Permanent Laboratory/On-site Calibration : Permanent Laboratory

Type of Service		Calibration Scope	CMC (Level of Confidence Approximately 95 %)
Volumetric Apparatus	Pipette	From 0.2 μ L up to 10 μ L	0.05 μ L
		More than 10 μ L up to 120 μ L	0.24 μ L
		More than 120 μ L up to 150 μ L	0.30 μ L
		More than 150 μ L up to 250 μ L	0.47 μ L
		More than 250 μ L up to 1000 μ L	1.9 μ L
		More than 1000 μ L up to 1200 μ L	2.4 μ L
		More than 1200 μ L up to 2500 μ L	5.6 μ L
		More than 2500 μ L up to 5 mL	13 μ L
		More than 5 mL up to 10 mL	34 μ L