

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. JCSS0089

Name of Laboratory

Lab Products & Services, JCSS Calibration Service
Sartorius Japan K.K.

Address of Laboratory

1-2-34 Ichigaya Sadohara-cho Shinjuku-ku Tokyo, JAPAN

Accreditation Scope Mass (as attached)

Accreditation Criterion ISO/IEC 17025:2005

Date of Initial Accreditation : 2000-12-21

Latest Date of Issue : 2013-10-17

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.

COPY

General Field of Calibration : Mass

Date of Initial Accreditation of the Field : 2000-12-21

Permanent Laboratory/On-site Calibration : Permanent Laboratory

Type of Service		Calibration Scope	CMC(Level of Confidence Approximately 95 %) (Conventional Mass)	
Weight	Weight	20 kg	12 mg	
		10 kg	5.3 mg	
		5 kg	1.5 mg	
		2 kg	0.7 mg	
		1 kg	0.30 mg	
		500 g	0.15 mg	
		200 g	0.06 mg	
		100 g	0.030 mg	
		50 g	0.019 mg	
		20 g	0.015 mg	
		10 g	0.012 mg	
		5 g	0.010 mg	
		2 g	0.008 mg	
		1 g	0.006 mg	
		500 mg	0.0048 mg	
		200 mg	0.0037 mg	
		100 mg	0.0030 mg	
		50 mg	0.0024 mg	
		20 mg	0.0019 mg	
		10 mg	0.0017 mg	
		5 mg	0.0017 mg	
		2 mg	0.0017 mg	
		1 mg	0.0017 mg	
		Deadweight	More than 10 kg up to 20 kg	100 mg
			More than 5 kg up to 10 kg	50 mg
			More than 2 kg up to 5 kg	25 mg
			More than 1 kg up to 2 kg	10 mg
			More than 500 g up to 1 kg	5 mg
	More than 200 g up to 500 g		2.5 mg	
	More than 100 g up to 200 g	1.0 mg		
	More than 50 g up to 100 g	0.5 mg		
	More than 20 g up to 50 g	0.30 mg		
	More than 10 g up to 20 g	0.25 mg		
More than 5 g up to 10 g	0.20 mg			
More than 2 mg up to 5 g	0.15 mg			

Permanent Laboratory/On-site Calibration : On-site Calibration

Type of Service		Calibration Scope	CMC(Level of Confidence Approximately 95 %) (Conventional Mass)
Weight	Weight	20 kg	300 mg
		10 kg	150 mg

COPY

General Field of Calibration : Mass

Date of Initial Accreditation of the Field : 2003-12-10

Permanent Laboratory/On-site Calibration : Permanent Laboratory and On-site Calibration

Type of Service		Calibration Scope	CMC (Level of Confidence Approximately 95 %)	
			Permanent	Laboratory On-Site Calibration
Scale	Non-Automatic Electronic Weighing Instruments	1 mg	2.6 µg	2.6 µg
		2 mg	2.6 µg	2.6 µg
		3 mg	5.2 µg	5.2 µg
		4 mg	5.2 µg	5.2 µg
		5 mg	2.6 µg	2.6 µg
		6 mg	5.2 µg	5.2 µg
		7 mg	5.2 µg	5.2 µg
		8 mg	7.8 µg	7.8 µg
		9 mg	7.8 µg	7.8 µg
		10 mg	2.6 µg	2.6 µg
		15 mg	5.2 µg	5.2 µg
		20 mg	2.9 µg	3.0 µg
		30 mg	5.5 µg	5.5 µg
		40 mg	5.9 µg	5.9 µg
		50 mg	3.7 µg	3.7 µg
		60 mg	6.3 µg	6.3 µg
		70 mg	6.6 µg	6.6 µg
		80 mg	9.2 µg	9.2 µg
		90 mg	9.5 µg	9.5 µg
		100 mg	4.6 µg	4.6 µg
		120 mg	7.5 µg	7.5 µg
		150 mg	8.3 µg	8.3 µg
		200 mg	5.7 µg	5.7 µg
		250 mg	9.4 µg	9.4 µg
		300 mg	11 µg	11 µg
		350 mg	14 µg	14 µg
		400 mg	12 µg	12 µg
		450 mg	15 µg	15 µg
		500 mg	7.4 µg	7.4 µg
		550 mg	11 µg	11 µg
600 mg	12 µg	12 µg		
650 mg	16 µg	16 µg		
700 mg	13 µg	13 µg		
750 mg	17 µg	17 µg		
800 mg	18 µg	18 µg		
850 mg	22 µg	22 µg		
900 mg	19 µg	19 µg		
1 g	9.2 µg	9.3 µg		
More than 1 g up to 1.5 g	17 µg	17 µg		
More than 1.5 g less than 2 g	23 µg	24 µg		

COPY

	2 g	13 µg	13 µg
	More than 2 g less than 3 g	26 µg	26 µg
	3 g	22 µg	22 µg
	4 g	25 µg	25 µg
	5 g	16 µg	17µg
	6 g	25 µg	26 µg
	7 g	28 µg	29 µg
	8 g	37 µg	39 µg
	9 g	40 µg	42 µg
	10 g	19 µg	23 µg
	15 g	34 µg	39 µg
	20 g	24 µg	35 µg
	25 g	39 µg	51 µg
	30 g	43 µg	57 µg
	40 g	0.049 mg	0.071 mg
	50 g	0.035 mg	0.073 mg
	More than 50 g less than 100 g	0.081 mg	0.14 mg
	100 g	0.056 mg	0.14 mg
	More than 100 g less than 150 g	0.096 mg	0.19 mg
	150 g	0.088 mg	0.21 mg
	More than 150 g less than 200 g	0.15 mg	0.28 mg
	200 g	0.11 mg	0.28 mg
	More than 200 g up to 300 g	0.21 mg	0.45 mg
	More than 300 g up to 400 g	0.26 mg	0.59 mg
	More than 400 g up to 500 g	0.31 mg	0.73 mg
	More than 500 g up to 600 g	0.37 mg	0.87 mg
	More than 600 g up to 2000 g	1.8 mg	3.3 mg
	More than 2000 g up to 5000 g	3.2 mg	7.4 mg
	More than 5000 g up to 8000 g	15 mg	24 mg
	More than 8000 g up to 14 kg	31 mg	44 mg
	More than 14 kg up to 35 kg	0.16 g	0.22 g
	More than 35 kg up to 64 kg	0.50 g	0.55 g
	More than 64 kg up to 150 kg	2.0 g	2.9 g
	More than 150 kg up to 300 kg	4.0 g	5.9 g
	More than 300 kg up to 600 kg	33 g	54 g