

IH-Centrifuge L and IH-Incubator L

L-Range Instruments













IH-Centrifuge L

Innovation

- Only one instrument for both IH-Cards and tubes
- Head for post-transportation IH-Card centrifugation with adapted program (5 minutes at 85 g)
- Panel PC touch screen
- Automatic head type detection and selection of the appropriate settings
- Multi-language software

Main Features

- Only one centrifuge for tubes and IH-Cards
- Automatic detection of the head type
- Automatic selection of the appropriate program
- Head for post-transportation IH-Card centrifugation with adapted program (5 minutes at 85 g)
- 28-tube capacity
- Panel PC touch screen with easy-to-use interface (icon based)
- 4 compatible centrifuge heads
- Multi-language software
- FDA 510(k)-cleared device

Technical Data

Performance	
RCF maximum	1,000 g (for tubes)
RCF	85 g (for IH-Cards)
Dimensions	
Width	380 mm (15.0 in)
Height	230 mm (9.1 in)
Depth	500 mm (19.7 in)
Weight, not including rotors	13 kg (28.7 lb)

Ordering Information

Catalog #	Description
009301	IH-Centrifuge L , 1 unit
009410	Head 24 IH-Cards , 1 unit
009415	Head 28 tubes , 1 unit
009413	Head 12 tubes , 1 unit
009414	Head 2 racks , 1 unit











IH-Incubator L

Innovation

- Only one instrument for both IH-Cards and tubes
- Incubation zones are independent and can be deactivated
- Panel PC touch screen
- Multi-language software

Main Features

- Incubator for both tubes and IH-Cards
- Independent incubation areas (IH-Cards and/or tubes)
- Up to 36 IH-Card and 24 tube capacity
- Panel PC touch screen with easy-to-use interface (icon based)
- Multi-language software
- Batch-specific timers (3 for IH-Cards, 2 for tubes)
- Standard program: 15 minutes at 37°C
- Programmable incubation time
- FDA 510(k)-cleared device

Technical Data

Dimensions

Width	380 mm (15.0 in)
Height	230 mm (9.1 in)
Depth	500 mm (19.7 in)
Weight, not including rotors	13 kg (28.7 lb)

Ordering Information

Catalog #	Description
009303	IH-Incubator L , 1 unit

Visit bio-rad.com/ih-manual-workplace for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.
© 2026 Bio-Rad Laboratories, Inc.



Bio-Rad
Laboratories, Inc.

Clinical Diagnostics Group	Website www.bio-rad.com/diagnostics	U.S. 1 800 224 6723	Australia 61 2 9914 2800	Austria 00 800 00 24 67 23	Belgium 00 800 00 24 67 23
	Brazil 55 11 3065 7550	Canada 1 800 361 1808	China +86 21 6169 8500	Czech Republic 00 800 00 24 67 23	Denmark 00 800 00 24 67 23
	Finland 00 800 00 24 67 23	France 00 800 00 24 67 23	Germany 00 800 00 24 67 23	Greece 30 210 7774396	Hong Kong 852 2789 3300
	Hungary 00 800 00 24 67 23	India 91 124 4029300	Israel 00 800 00 24 67 23	Italy 00 800 00 24 67 23	Japan 81 3 6361 7070
	Luxembourg 00 800 00 24 67 23	Mexico 52 55 5488 7670	The Netherlands 00 800 00 24 67 23	New Zealand 64 9415 2280	Norway 00 800 00 24 67 23
	Poland 00 800 00 24 67 23	Portugal 00 800 00 24 67 23	Republic of Korea 82 080 007 7373	Russia 7 495 721 14 04	Singapore 65 6415 3170
	South Africa 27 11 442 8508	Spain 00 800 00 24 67 23	Sweden 00 800 00 24 67 23	Switzerland 00 800 00 24 67 23	Taiwan 886 2 2578-7189
	Thailand 662 651 8311	United Arab Emirates 971 4 818 7300	United Kingdom 00 800 00 24 67 23		