



Certificate of Compliance

Certificate: 1432240 (LR 34862)

Master Contract: 169502

Project: 70015996

Date Issued: November 20, 2014

Issued to: Beckman Coulter Inc.
250 S. Kraemer Blvd
Brea, CA 92822
USA
Attention: Jeffrey Gonzalez

The products listed below are eligible to bear the CSA Mark shown with adjacent indicators 'C' and 'US' for Canada and US or with adjacent indicator 'US' for US only or without either indicator for Canada only.



Ruo Xuan Lim

Issued by: Ruo Xuan Lim

PRODUCTS

CLASS 8721 85 - ELECTRICAL EQUIPMENT FOR LABORATORY USE - Certified to US Standards

CLASS 8721 05 - LABORATORY EQUIPMENT - Electrical

Avanti Centrifuge, J Series, Models and Ratings as follows:

Model	BCI P/N	Electrical Ratings
J-30I	363118	200/208/240V, 50/60 Hz, 24A
J-30I	363120	230V, 50Hz, 24A
J-30I with elutriation	363122	200/208/240V, 50/60 Hz, 24A
J-30I with elutriation	363124	230V, 50Hz, 24A
HP-30I	366418	200/208/240V, 50/60 Hz, 24A
HP-30I with elutriation	366422	200/208/240V, 50/60 Hz, 24A
J-HC with elutriation	367501	200/208/240V, 50/60 Hz, 24A
J-HC with elutriation	367502	230V, 50Hz, 24A
J26XP	393124	200/208/240V, 50/60 Hz, 24A
J26XP	393125	230V, 50Hz, 24A
J26XPI	393127	200/208/240V, 50/60 Hz, 24A



Certificate: 1432240 (LR 34862)

Master Contract: 169502

Project: 70015996

Date Issued: November 20, 2014

J26XPI	393128	230V, 50Hz, 24A
J26XP with elutriation	393130	200/208/240V, 50/60 Hz, 24A
J26XP with elutriation	393131	230V, 50Hz, 24A
J26XPI with elutriation	393133	200/208/240V, 50/60 Hz, 24A
J26XPI with elutriation	393134	230V, 50Hz, 24A
HP-J26XP	393136	200/208/240V, 50/60 Hz, 24A
HP-J26XPI	393137	200/208/240V, 50/60 Hz, 24A
HP-J26XP with elutriation	393138	200/208/240V, 50/60 Hz, 24A
HP-J26XPI with elutriation	393139	200/208/240V, 50/60 Hz, 24A
J-26S XP IVD	B06102	200/208/240V, 50/60 Hz, 24A
J-26S XP IVD	B06103	230V, 50Hz, 24A
J-26S XP IVD with elutriation	B06108	200/208/240V, 50/60Hz, 24A
J-26S XP IVD with elutriation	B06109	230V, 50Hz, 24A
JXN-26	B34183	200/208/240V, 50/60Hz, 24A
JXN-26	B37912	230V, 50Hz, 24A
JXN-26	B34182	220V, 50Hz, 12A (3 Phase)
JXN-26 IVD	B38619	200/208/240V, 50/60Hz, 24A
JXN-26 IVD	B38623	230V, 50Hz, 24A
JXN-26 IVD	B38618	220V, 50Hz, 12A (3 Phase)
JXN-30	B37913	230V, 50Hz, 24A
JXN-30	B34193	200/208/240V, 50/60Hz, 24A
JXN-30 (3 Phase)	B34192	220V, 50Hz, 12A
JXN-30 IVD	B38624	230V, 50Hz, 24A
JXN-30 IVD	B38621	200/208/240V, 50/60Hz, 24A
JXN-30 IVD (3 Phase)	B38620	220V, 50Hz, 12A

Notes:

The above models are Equipment Class I, Pollution Degree 2, Installation Category II.

APPLICABLE REQUIREMENTS

CAN/CSA-C22.2 No. 61010-1-04 - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements

UL Std. No. 61010-1 (2nd Edition) - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use - Part 1: General Requirements

CAN/CSA-C22.2 No. 61010-2-020:2009 - Safety requirements for electrical equipment for measurement, control and laboratory use - Part 2-020, Particular requirements for laboratory centrifuges

CAN/CSA-C22.2 No. 61010-2-101-04 - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use – Part 2-101: Particular Requirements for In Vitro Diagnostic (IVD) Medical Equipment – First Edition