



Certificate of Compliance

Certificate: 80103918

Master Contract: 169502

Project: 80103918

Date Issued: 2022-01-27

Issued To: Beckman Coulter Inc. - Brea
250 S. Kraemer Blvd
M/S C1.NW.02
Brea, California, 92822
United States

Attention: Bret Widdifield

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.

Issued by: Tony Nguyen
Tony Nguyen



PRODUCTS

CLASS - C872106 - ELECTRICAL LABORATORY EQUIPMENT Electrical Laboratory Equipment

CLASS - C872186 - ELECTRICAL LABORATORY EQUIPMENT Certified to US Standards

Next Generation Library Prep System, Rated 100-240V, 50/60 Hz, 10 A.

Models:

Biomek Ngenius Dx Next Generation Library Prep System (C62704, IVD)

Biomek Ngenius Next Generation Library Prep System (C62703, Non-IVD)

Notes:

1. The above models are detachable cord connected, Equipment Class I, Pollution Degree 2, and Overvoltage Category II.



Certificate: 80103918
Project: 80103918

Master Contract: 169502
Date Issued: 2022-01-27

2. Mode of operation: Continuous
3. Environmental Conditions: Normal: 19°C to 25°C, 2000 m max, 20-60% R.H. (non-condensing)

APPLICABLE REQUIREMENTS

CAN/CSA C22.2 No. 61010-1-12, UPD1: 2015, UPD2: 2016, AMD1: 2018

Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements

CSA C22.2 No. 61010-2-081:19 - Safety requirements for electrical equipment for measurement, control and laboratory use — Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes

CSA-C22.2 No. 61010-2-101:19 - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use - Part 2-101: Part 2-101: Particular requirements for in vitro diagnostic (IVD) medical equipment

UL 61010-1, 3rd Edition (2012), AMD1: 2018

Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements

UL 61010-2-081, 3rd Edition - Safety requirements for electrical equipment for measurement, control and laboratory use — Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes

UL 61010-2-101, 3rd edition (2019) - Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use - Part 2-101: Part 2-101: Particular requirements for in vitro diagnostic (IVD) medical equipment

MARKINGS

The manufacturer is required to apply the following markings:

- Products shall be marked with the markings specified by the particular product standard.
- Products certified for Canada shall have all Caution and Warning markings in both English and French.

Additional bilingual markings not covered by the product standard(s) may be required by the Authorities Having Jurisdiction. It is the responsibility of the manufacturer to provide and apply these additional markings, where applicable, in accordance with the requirements of those authorities.



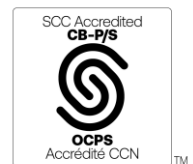
Certificate: 80103918
Project: 80103918

Master Contract: 169502
Date Issued: 2022-01-27

The products listed are eligible to bear the CSA Mark shown with adjacent indicators 'C' and 'US' for Canada and US (indicating that products have been manufactured to the requirements of both Canadian and U.S. Standards) or with adjacent indicator 'US' for US only or without either indicator for Canada only.

Notes:

Products certified under Class C872106, C872186 have been certified under CSA's ISO/IEC 17065 accreditation with the Standards Council of Canada (SCC). www.scc.ca





Supplement to Certificate of Compliance

Certificate: 80103918

Master Contract: 169502

*The products listed, including the latest revision described below,
are eligible to be marked in accordance with the referenced Certificate.*

Product Certification History

Project	Date	Description
80103918	2022-01-27	C/US certification of Next Generation Library Prep System, model Biomek Ngenius Dx (C62704, IVD) and Biomek Ngenius (C62703, Non-IVD).