



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
CF-260-120	CF-260 Centrifuge 110 VAC, 50/60 Hz
CF-260-220	CF-260 Centrifuge 220 VAC, 50/60 Hz

OMNI International declares under our sole responsibility that the products listed above conform with the following standards or other normative documents.

The CF-260 Centrifuge will meet the CE and UKCA specifications when applied in accordance with the manufacturer's instructions. This product also conforms to the corresponding UK Statutory Instruments (SI) requirements.

Directive 2014/30/EU Standards used:	Electromagnetic Compatibility Directive (EMC) EN 61326-1:2013 / IEC 61326-1:2012
Directive 2014/35/EU Standards used:	Low Voltage Directive (LVD) EN/IEC 61010-1:2010+A1:2016, EN 61010-1:2010/A1:2019 EN/IEC 61010-2-020:2017
Directive 2015/863/EU Standards used:	The Restriction of Hazardous Substances Directive (RoHS) EN 63000:2018
Directive 2012/19/EU Standards used:	Waste Electrical and Electronic Equipment Directive (WEEE) EN 50419:2006

This declaration of conformity is issued under the sole responsibility of OMNI International.

Signed for and on behalf of:

Paul H. Roberts
Senior Product Manager
November 2023