



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
CF-350-120	CF-350 Centrifuge 110 VAC, 50/60 Hz
CF-350-220	CF-350 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-350 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.

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