



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
06-021	OMNI Prep Homogenizer 115/220 V

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

The OMNI Prep Homogenizer 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
	CFR 47 FCC Part 15 Subpart B. Class A
<b>Directive 2014/35/EU</b> Standards used:	<b>Low Voltage Directive (LVD)</b>
	EN/IEC 61010-1:2010+A1:2016
	EN/IEC 61010-2-051:2015
<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