



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
29-060-450-120	OM500 Ultrasonic Homogenizer 110 VAC, 50/60 Hz
29-060-450-220	OM500 Ultrasonic Homogenizer 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 OM500 Sonicator 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
	EN 55011:2016, CFR 47 FCC Part 18, ICES-003 Issue 5
	EN 61000-3-2:2018, EN 61000-3-3:2013 3rd Edition
<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 61000-4-2:2009, EN 61000-4-3:2006 + A2:2010
	EN 61000-4-4:2012, EN 61000-4-5:2014, EN 61000-4-6:2014
	EN 61000-4-11:2004 + A1:2017
<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
<b>For Canada and USA</b>	CAN/CSA-C22.2 No. 61010-1:2012-05
	CAN/CSA-C22.2 No. 61010-2-010:2015-01
	CAN/CSA-C22.2 No. 61010-2-081:2015-11
	UL 61010-1:2012-05
	UL 61010-2-010:2015-01
	UL 61010-2-081:2015-11

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