



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
17106	OMNI Mixer Homogenizer 220V
17506	OMNI Macro Homogenizer 220V
MES-220	OMNI Macro ES Digital Programmable Homogenizer 220V
MESP-220	OMNI Macro ES Homogenizer Package 220V
17105	OMNI Mixer Homogenizer 115V
17505	OMNI Macro Homogenizer 115V
MES-115	OMNI Macro ES Digital Programmable Homogenizer 115V
MESP-115	OMNI Macro ES Homogenizer Package 115V

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

The OMNI Mixer/Macro/Macro ES 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
	CFR 47 FCC Part 15 Subpart B. Class A
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 61010-2-051:2015
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
For Canada and USA	CAN/CSA-C22.2 No. 61010-1:2012-05
	CAN/CSA-C22.2 No. 61010-2-010:2015-01
	UL 61010-1:2012-05
	UL 61010-2-010:2015-01

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

OMNI International, 935 Cobb Place Blvd., Kennesaw, GA 30144
An ISO 9001 Company

PN: DOC-17106