



## PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

| Part Number   | Description  |
|---------------|--|
| 12-500        | OMNI THQ - Digital Tissue Homogenizer  |
| 12-200-220    | OMNI THQ - Digital Tissue Homogenizer  |
| 12-200-220-05 | OMNI THQ - Digital Tissue Homogenizer with 5 x 75 mm Stainless Steel Generator Probe   |
| 12-200-220-07 | OMNI THQ - Digital Tissue Homogenizer with 7 x 110 mm Stainless Steel Generator Probe  |
| 12-200-220-10 | OMNI THQ - Digital Tissue Homogenizer with 10 x 110 mm Stainless Steel Generator Probe |

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

The OMNI THQ - Digital Tissue 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><br>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><br>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><br>Standards used: | <b>The Restriction of Hazardous Substances Directive (RoHS)</b>   |
|   | EN 63000:2018   |
| <b>Directive 2012/19/EU</b><br>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