



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

| Part Number | Description |
|-------------|----------------|
| 19-050A | Bead Ruptor 12 |

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

The Bead Ruptor 12 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 61326-1:2021 / IEC 61326-1:2020 |
| | 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
May 2024