



PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

Part Number	Description
25-010	Bead Ruptor 4

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

The Bead Ruptor 4 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 EN/IEC 61010-2-081:2015
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
	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