

EU declaration of conformity

Product model:

DropletQuant

Manufacturer and contact authorized to compile the relevant technical documents:

Trinean nv
 Dulle Grietlaan 17/3
 9050 Gentbrugge
 Belgium
www.trinean.com

This declaration of conformity is issued under the sole responsibility of the manufacturer.

The product described above is in conformity with the relevant Union harmonisation legislation:

2014/35/EU Low Voltage Directive
 2014/30/EU Electromagnetic Compatibility Directive
 2011/65/EU Restriction of Hazardous Substances (RoHS)

In order to achieve conformity, following harmonized standards were consulted:

IEC61010-1:2010 (Third Edition)
 Safety requirements for electrical equipment for measurement, control, and laboratory use.

IEC61010-2-081:2015 (Second Edition)
 Safety requirements for electrical equipment for measurement, control, and laboratory use.

EN61326-1 (2013) Electrical equipment for measurement, control, and laboratory use

EN55011 (2009) + am1	EN61000-4-5 (2014)
EN61000-3-2 (2014) + am1 + am2	EN61000-4-6 (2014)
EN61000-4-2 (2009)	EN61000-4-8 (2010)
EN61000-4-3 (2006) + am1 + am2	EN61000-4-11 (2004)
EN61000-4-4 (2012)	

Signed for and on behalf of:

Trinean nv
 At Gentbrugge 3rd of Februari, 2017



Philippe Stas
 CEO