

PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

FT-IR Spectrometer, Model Frontier

This PerkinElmer product conforms to the regulations stipulated in the CE Mark requirements for the EMC Directive (2014/30/EU), the Low Voltage Directive (2014/35/EU), and the RoHS 2 Directive (2011/65/EU as amended by (EU) 2015/863), and the corresponding UK Statutory Instruments (SI) for the UKCA Mark, the Electromagnetic Compatibility Regulations 2016 (SI 2016 No. 1091), the Electrical Equipment (Safety) Regulations 2016 (SI 2106 No. 1101), and the Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012 (SI 2012 No. 3032):

EN 55011:2009 + A1:2010 Group 1, Class A, EMC -- RF Characteristics of ISM Equipment
EN 61326-1:2013, EMC -- Requirements for Electrical Equipment for Laboratory Use
EN 61000-4-2:2009, EMC -- Electrostatic Discharge Requirements
EN 61000-4-3: 2006 + A1:2008 + A2:2010, EMC -- Radiated Electromagnetic Field Requirements
EN 61000-4-4: 2012, EMC -- Electrical Fast Transient/Burst Requirements
EN 61000-4-5:2006, EMC -- Surge Immunity Requirements
EN 61000-4-6:2009, EMC -- Conducted Disturbances (induced by RF fields) Requirements
EN 61000-4-11:2004, EMC -- Voltage Dips, Short Interruptions, Voltage Variations Requirements
EN 61000-3-2:2006 + A2:2009, EMC -- Harmonic Current Emissions
EN 61000-3-3:2013, EMC -- Voltage Fluctuations and Flicker
EN 61010-1:2010, Safety Requirements for Electrical Equipment for Laboratory Use
EN 60825-1:2014, Safety of laser products - Part 1: Equipment classification and requirements
EN 63000:2018, Technical Documentation for the assessment of electrical and electronic products with respect to the RoHS



CAN/CSA C22.2 No. 61010-1-12 + GII + GI2 (R2017), Safety Requirements for Electrical Equipment for Laboratory Use
UL 61010-1:2012 R4.16, Safety Requirements for Electrical Equipment for Laboratory Use
ICES-003, Issue 6, Class A, Radiated and Conducted Emissions
FCC Part 15, Class A, For Radiated and Conducted Emissions
FDA CDRH 21 CFR 1000-1050 Laser Products

AS/NZS CISPR 11:2011
Korean Radio Waves Act, Article 58-2, Clause 3

This product also conforms to the Waste Electrical and Electronic Equipment Directive (WEEE, 2012/19/EU), and the Regulation on Registration, Evaluation, Authorisation and Restriction of Chemicals (REACH, EC 1907/2006), and the corresponding UK Statutory Instruments (SI), the WEEE Regulations 2013 (SI 2013 No. 3113), and the REACH Enforcement Regulations 2008 (SI 2008 No. 2852).

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

Signed for and on behalf of:



Richard Covell
Compliance Engineering Manager

July 22, 2021

PerkinElmer, 710 Bridgeport Avenue, Shelton, CT 06484 USA
[in the EU: DANI ANALITICA S.R.L., Part of PerkinElmer Group, Viale Brianza, 87 20093 Cologno Monzese MI - ITALY]
[in the UK: PerkinElmer, Chalfont Road, Buckinghamshire, Seer Green, HP9 2FX, United Kingdom]
An ISO 9001 Company