





1.12537.0001

UV lamp 254 nm

Safety instructions

	<ul style="list-style-type: none">• Please read these instructions carefully before using the instrument.• The protection provided by this equipment may be compromised if it is used incorrectly.• Obey the instructions that are given to you. Do not perform any work on the inside of the system.• Keep these instructions for future reference.
	<ul style="list-style-type: none">• The instrument is emitting UV light, which is a potential hazard.• Do not expose eyes to rays from this lamp.• Protective glasses must be worn.
	<ul style="list-style-type: none">• Certificate of conformation No. NSE-E07101536: the device has been tested for electromagnetic compatibility and found in compliance with the council directive EMC 2004/108/EC.
	<ul style="list-style-type: none">• Do not dispose the instrument or the batteries in normal garbage.• Dispose the instrument and the batteries in a recycling bin suitable for the type of product.

UV lamp

Ultraviolet (UV) light is an electromagnetic radiation, with wavelenths below 400 nm. The UV lamp emits visible light, but also ultraviolet (UVC and UVB) wavelengths around 254 nm, which are used to detect fluorescence in laboratory applications such as chromatography, microbiology or other areas. In contrary to lamps with wavelength filters and very narrow spectrum, this lamp has a broader spectrum for quick analyses or screening purposes.

Additionally a small flash lamp is built in, allowing orientation in dark rooms.

Instructions for use

1. Open the lid of the instrument and insert 4 AA batteries (switch in middle position)
2. Before switching on, ensure that your eyes are not exposed to the UV light.
3. The switch allows the following selections: Left = UV lamp, Middle = Off, Right = Flash lamp

Specifications

Item number	1.12537.0001
Wavelength	A) Ultraviolet spectrum around 254 nm for laboratory applications, B) Visible light / Flash lamp
Output	4 Watt
Dimensions	16.1 x 2.4 x 5.5 cm
Weight	99.8 grams
Power supply	AA 1.5V Battery x 4 pcs (6V)

Handling and storage

The instrument is made for indoor use in clean and dry laboratory environment. Do not use any liquid or aerosol cleaning agent or solvent on or near the instrument.

The lamp is battery powered, 4 AA/Mignon/R6 batteries (1.5V) are included.

Do not charge non-rechargeable batteries.

Guarantee

The UV lamp is guaranteed for six months from the date of delivery.

Faults or functional disturbances brought about by incorrect handling are excluded.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

© 2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck, Supelco, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly available resources.

Merck Life Science KGaA,
64271 Darmstadt, Germany,
Tel. +49(0)6151 72-2440
www.sigmaaldrich.com

MERCK