

General Information

To ensure the safety of personnel and the environment, it is extremely important that both the device and its accessories be thoroughly disinfected before any maintenance or servicing, as well as before leaving the laboratory. Please ensure that the procedure described below complies with your national, regional, and local safety regulations. After completing the disinfection and decontamination procedure, fill out the form on the next page and sign it. If the device has been contaminated with a hazardous substance, it must also be countersigned by a responsible person, a supervisor, a safety officer, or a bio-safety officer. Subsequently, submit the certificate to your contact person at BMG LABTECH GmbH. If the equipment is being shipped, also include a printed copy with the shipment.

Preparation steps

Before decontaminating your item:

- It is advisable to consult the Operating Manual (Chapter "Servicing"), if applicable, as reference for which decontamination product(s) may be suitable when cleaning your BMG LABTECH product.
- The decontamination should be done by authorized personnel only in a well ventilated location.
- It is recommended to wear Personal Protective Equipment (PPE) such as a laboratory coat, gloves, safety glasses and any other PPE as recommended by your local safety protocols and the type of the hazards handled.
- Consult the disinfectant's Safety Data Sheet (SDS) before using it.
- Follow strictly the safety instructions of your laboratory for decontamination.

Use of disinfectant

Only propriety disinfectant or alcohol should be used for the purpose of disinfection. Use suitable disinfectants for the (bio-)hazard you handled. Please consult your lab host for any specific cleaning materials for decontamination. Here are some examples of disinfectants (BMG LABTECH assumes no responsibility for their effectiveness):

- Decon laboratories products
- Microcide SQ
- Steris Spor Klenz
- 1% Virkon solution or equivalent spray solution
- 70% Ethanol
- 70% Isopropanol

Procedure using liquid disinfectants

All parts of the instrument, which could potentially get into contact with patient sera or positive samples, must be handled as if they were hazardous. Please follow all the instructions carefully for a successful disinfection:

- 1) Press the in/out button briefly to move the plate carrier out.
- 2) Disconnect the instrument from the main power supply.
- 3) Remove the USB cable from the connector.
- 4) Clean all outside surfaces of the instrument and the plate carrier carefully with cotton wool, which has been soaked in disinfectant.
- 5) Open all lids, doors and slide doors and move out the plate carrier. Place the instrument in a large plastic bag along with the cotton wool that has been soaked in disinfectant. Ensure that the wool does not touch the instrument.
- 6) Close and seal the bag.
- 7) Keep the instrument in the plastic bag for at least 24 hours.
- 8) After exposure time has lapsed, remove the instrument from the plastic bag and clean all outside surfaces of the instrument with cotton wool that has been soaked in alcohol solution.
- 9) Repeat the procedure for disinfection on any accessories, which will be returned with the instrument.
- 10) Fill the form on the next page and signed it.

