

Environmental Policy of Bio-Rad Laboratories GmbH, Munich, Germany

As a manufacturer of life science and diagnostic products, Bio-Rad Laboratories wants to look after its corporate responsibility for the protection of people and the environment. The protection of environmental and human health are part of ensuring corporate efficiency. Thus, we want to integrate them into our corporate policy and direct the actions of our employees towards these essential goals.

It is therefore our duty to reduce any effects on the environment to a minimum. This concept shall enter the company's philosophy in addition to the pursuit for safe and profitable production and continuous improvement of our products.

1. Advancement of environmental consciousness in our employees

All managers at Bio-Rad Laboratories support the awareness of environmental protection of their employees through model behavior, goal-oriented instructions and information. Effective environmental protection is based on the commitment and motivation of our coworkers.

2. Environmental protection provisions for new procedures and products

Environmental and occupational protection at Bio-Rad Laboratories already starts at the development and planning of new products and procedures.

3. Safe facilities and procedures

Bio-Rad Laboratories continuously reviews and evaluates the effects of existing operations, procedures and facilities. All environmental impact is reduced to an extent which can be reached with economically justifiable applications of the best technology available. Provisions are made and developed continuously, so that in the case of a production failure, damages caused by accident on people and the environment are kept to a minimum through optimal reaction.

4. Fending off environmental and human health hazards

The people in charge at Bio-Rad Laboratories take on the necessary precautions to avoid environmental hazards due to products, waste and processes. In the case of deviations from planned procedures, possible damaging effects on the protection and health of employees and residents are excluded through effective corrective measures.

5. Evaluation of environmental protection activities

Control procedures at Bio-Rad Laboratories ensure that the environmental policy is carried out on location and the environmental goals are attained. Systematic environmental audits ensure a continuous improvement of environmental protection. The management continuously informs itself of the status of the environmental protection management system.

6. Dialog with the public

Bio-Rad Laboratories wants to gain additional acceptance for products and the company by active information of the public about environmental and environment-based activities. Suggestions from the public will be accepted and reviewed with regard to their practicability.

7. Customer advice about environmental aspects

Customers of Bio-Rad Laboratories will be informed about the safe and environmentally conserving handling and environmentally compatible disposal of products and their packaging.

8. Cooperation with government agencies

Bio-Rad Laboratories conducts a cooperative interaction with governmental agencies for the coordination and improvement of environmental protection activities. This ensures that processes are maintained up to the latest standard and the compliance with applicable regulations and governmental requirements as minimum standard.

9. Evaluation of our contractors

We are committed that our vendors and service contractors apply the same standards regarding environmental protection and protection like our company itself. We want to persuade our contractual partners to comply with environmentally compatible processes by using a special evaluation procedure.

10. Efficient environmental protection organization

The management of Bio-Rad Laboratories has determined the responsibilities for the creation of an environmental protection organization, to ensure and constantly improve

environmental protection at our location. All responsibilities and processes are clearly defined in the framework of this organization.

11. Continuous improvement of environmental protection

Bio-Rad Laboratories wants to constantly advance the achieved status of operational environmental protection. In the foreground is the constant improvement of all environmental protection activities, including all technical and organizational procedures and processes, provided they seem to be economically meaningful.



**Bio-Rad
Laboratories GmbH**

*D-80901 München · Postfach 45 01 33
D-80939 München · Heidemannstraße 164
Telefon (089) 318 84-0
Telefax (089) 318 84-100
Internet: www.bio-rad.com*