

March 2010

## **SAFETY REGULATIONS FOR SERVICING LABORATORY INSTRUMENTS**

Dear Customer

When working with biological and/or radioactive materials and the instruments used thereby, it cannot be ruled out that service personnel could be exposed to health risks by coming into contact with residues of hazardous substances, especially when the instrument has been used for making measurements using radioactive, infectious or toxic substances.

The current regulations and laws, as well as the extended guidelines and norms, stipulate that we, as a manufacturer of measuring systems, observe more stringent safety regulations in order to ensure the safety of our employees.

Moreover, the environmental regulations issued by the environmental protection and industrial inspection board as well as the quality assurance system DIN/ISO 9001, which was awarded to BERTHOLD TECHNOLOGIES GmbH & Co. KG in 1990, also have to be observed.

Therefore, prior to allowing any repair work in your laboratory or before returning the instrument to us we would ask you, either to carefully clean, disinfect or decontaminate the instrument or components to be serviced, or confirm that the instrument or components have not come into contact with any hazardous substances.

On the reverse side you will find the „Confirmation on Decontamination/Declaration of Safety“ which should be filled out and attached to the shipping papers together with your repair order, or handed out directly to our service technicians in your laboratory.

**We are unable to commence repair work without a declaration that the instrument has been decontaminated. Should the declaration not be received within three weeks, for safety reasons we have to return the instruments unrepaired freight forward.**

For radiation protection instruments and detectors returned without a decontamination declaration we offer to test for radioactive contamination at the price of 100,00 Euro.

For further questions, please do not hesitate to contact directly our Service Department.

Yours sincerely

**BERTHOLD TECHNOLOGIES GmbH & Co. KG**

p.t.o.

**Important: please affix this confirmation outside the shipping box**

## Confirmation on Decontamination / Declaration of Safety

(PLEASE USE CAPITAL LETTERS !!!)

instrument / component: _____	serial no.: _____
<input type="checkbox"/> I hereby confirm that the instrument or component specified above has not been used with any hazardous or contagious samples or reagents	
instrument or component has come into contact with:	
<input type="checkbox"/> radioactive substances Isotope(s): _____ _____	means of decontamination applied: _____ _____
<input type="checkbox"/> chemical reagents specify: _____ R and S rules: _____	means of decontamination applied: _____ _____
<input type="checkbox"/> contagious biological material specify: _____ _____	means of decontamination applied: _____ _____
indicate security level of the laboratory the instrument has been used in <input type="checkbox"/> S1 <input type="checkbox"/> S2 <input type="checkbox"/> S3 <input type="checkbox"/> S4	
<input type="checkbox"/> I hereby confirm that the instrument or component specified above was not contaminated with any of the above mentioned substances / reagents / agents	
<input type="checkbox"/> I hereby confirm that the instrument or component specified above was decontaminated / cleansed using the appropriate method	
date: _____	signature: _____
name: _____	address: _____
title: _____	_____
e-mail: _____	_____
phone: _____	fax: _____

DC0000DV2 Rev.05