

Decontamination Form

Any laboratory instrument used for clinical or research analysis maybe considered a biohazard and requires decontamination prior to handling. Universal precautions are suggested wherever applicable.

Before shipping any instrument back to Berthold Detection Systems the luminometers have to be cleaned in the following way:

1. Wash the housing with a moist cloth, if necessary use a mild detergent or 70% Isopropanol.
2. Use 70% Isopropanol for sterilizing the injection lines after 10 minutes incubation time. Rinse with distilled water thoroughly.
3. Empty the injection lines.
4. Open the access door, remove and clean the reflector, then clean the measurement chamber with the measurement window carefully. Decontaminate the instruments surfaces, measurement chamber and sample holder with 70% Isopropanol according to the description in the chapter "Maintenance". Install the reflector again according to the description.
5. Use cotton tipped applicators for hard to reach areas, if necessary.

Berthold Detection Systems only accepts instruments for repair together with a filled out decontamination form. Thank you.

Instrument Type: _____

Serial Number: _____

- ☐ I confirm that the specified instrument was decontaminated according to the above described decontamination procedure
- ☐ I confirm that the above specified instrument had no contact to any hazardous material.

Company	
Contact Person	
Position	
Telephone	

Date: _____

Signature: _____

Please copy and fax to:

Berthold Detection Systems, Bleichstr. 56-68, 75173 Pforzheim,
Tel.: +49-(0)7231-9206-0, Fax -50