



Biosciences

LI-COR Biosciences GmbH

Siemensstraße 25 A
61352 Bad Homburg

Fax: +49 6172 17177-99
Phone: +49 6172 17177-71

LI-COR Decontamination-Certificate

Please fill in this form and return with the instrument!

Name:

Company/University:

Department/Institute:

Street:

City:

Country:

Phone:

Fax:

Type of instrument:

Serial-no:

Possible Contamination (please specify):

Radioactive:

Biological

Chemical:

Other substances:

Amount of contamination:

Equipment / instrument decontaminated?

Yes:

No:

Method of decontamination:

Please contact LI-COR by phone:

+49 6172 171 77 22
or **+49 6172 171 77 71**

if you can't decontaminate the equipment / instrument, or if the activity is more than 5 cps.

Signature

Place, Date

Position