



Biosciences

LI-COR Biosciences UK Ltd.

St. John's Innovation Centre
Cowley Road
Cambridge CB4 0WS

Fax: +44 1223 422 105
Phone: +44 1223 422 104

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:

+44 1223 422 104
or **+49 6172 17 177 71**

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

Signature

Place, Date

Position