

# LI-COR CUSTOMER PRIVACY POLICY

## CONTENTS

SCOPE OF THIS POLICY .....	1
A NOTE ON PRIVACY SHIELD (INTERNATIONAL DATA TRANSFERS FROM THE EEA OR SWITZERLAND) .....	1
CATEGORIES OF INFORMATION COLLECTED .....	2
PURPOSES FOR PROCESSING INFORMATION .....	4
DATA TRANSFERS TO THIRD PARTIES .....	4
SECURITY OF INFORMATION .....	6
ACCESS .....	6
ADDITIONAL NOTICES FOR EU PERSONAL DATA ONLY .....	6
DATA CONTROLLER .....	6
LEGAL BASIS FOR PROCESSING .....	6
YOUR RIGHTS .....	6
HOW LONG WE RETAIN YOUR DATA .....	7
USE OF THIRD-PARTY WEB SERVICES .....	7
NOTICE TO END USERS .....	8
CHILDREN'S INFORMATION .....	8
CHANGES TO THIS PRIVACY POLICY .....	8
QUESTIONS OR COMPLAINTS .....	9
E-MAIL OPT-OUT .....	9

## SCOPE OF THIS POLICY

This policy describes how LI-COR, Inc., d/b/a LI-COR Biosciences, and its subsidiaries, LI-COR Ltd and LI-COR GmbH, (collectively, “LI-COR”, “we”, “us”, or “our”) collect, use, and disclose information from you, or that you may provide, when you visit a LI-COR Website (your “Website Information”), your personal data that we may receive from the European Union (your “EU Personal Data”), and your personal data that we may receive in the United States (“US”) from the European Economic Area (your “EEA Personal Data”) or Switzerland (your “Swiss Personal Data”). Your Website Information, EU Personal Data, EEA Personal Data, and Swiss Personal Data are collectively referred to in this policy as your “Information.” Website Information may include your EU, EEA, or Swiss Personal Data. The term “personal data” as used in this policy refers to information relating to an identified or identifiable natural person that is protected as personal data under applicable data protection law, including without limitation the General Data Protection Regulation 2016/679 (“GDPR”). The term “LI-COR Website” includes all

websites operated by us, including without limitation trainingcenter.licor.com, training.licor.com, www.fluxsuite.com, www.fluxsuite.net, www.fluxsuite.org, www.eddypro.com, www.eddypro.net, www.eddypro.org, www.eco2s.com, www.eco2s.net, learn.lambdau.net, www.lambdau.net, tovi.io, and licor.com.

This policy does not apply to any third-party websites, applications, software, or products that integrate with any of our services or products or that are accessible on any LI-COR Website or any third-party products, services, or businesses. Additionally, a separate services agreement that you enter into with us may govern the delivery, access, and use of certain services we offer and provide additional or different terms with respect to those certain services.

If you do not agree with the terms of this policy, do not access or use the LI-COR Websites or other services or products we offer. This policy went into effect on the date set forth below. We may update this policy from time to time as described in the section of this policy called Changes to this Privacy Policy.

## A NOTE ON PRIVACY SHIELD (INTERNATIONAL DATA TRANSFERS FROM THE EEA OR SWITZERLAND)

LI-COR recognizes that the European Economic Area (the “EEA”) and Switzerland have established strict protections regarding the handling of EEA and Swiss Personal Data, including requirements to provide adequate protection for EEA and Swiss Personal Data transferred outside of the EEA or Switzerland. To provide adequate protection for certain EEA and Swiss Personal Data received by LI-COR, Inc. in the US, LI-COR, Inc. has elected to self-certify to the EU-US Privacy Shield Framework and Swiss-US Privacy Shield Framework administered by the US Department of Commerce (“Privacy Shield”). LI-COR, Inc. adheres to the Privacy Shield Principles of Notice, Choice, Accountability for Onward Transfer, Security, Data Integrity and Purpose Limitation, Access, and Recourse, Enforcement, and Liability. If there is any conflict between the terms of this policy and the Privacy Shield Principles, the Privacy Shield Principles shall govern.

For purposes of enforcing compliance with the Privacy Shield, LI-COR, Inc. is subject to the investigatory and enforcement authority of the US Federal Trade Commission. For more information about the Privacy Shield, see the US Department of Commerce's Privacy Shield website located at: <https://www.privacyshield.gov>. To review LI-COR, Inc.'s representation on the Privacy Shield list, see the US Department of Commerce's Privacy Shield self-certification list located at: <https://www.privacyshield.gov/list>.

If you have a question or complaint related to LI-COR, Inc.'s participation in the EU-US or Swiss-US Privacy Shield, please contact us as provided in the Questions or Complaints section below. LI-COR, Inc. will investigate and attempt to resolve any complaints or disputes regarding the use or disclosure of your EEA or Swiss Personal Data within 45 days of receiving your complaint. For any unresolved complaints regarding EEA or Swiss Personal Data, LI-COR, Inc. has agreed to cooperate with ICDR/AAA EU-US Privacy Shield Program and ICDR/AAA Swiss-US Privacy Shield Program located in the United States. If you are unsatisfied with the resolution of your complaint regarding your EEA or Swiss Personal Data, you may contact ICDR/AAA Privacy Shield Program at <http://go.adr.org/privacyshield.html> for further information and assistance or to file a complaint. The services of ICDR/AAA are provided at no cost to you.

Further, with regard to EEA Personal Data, you may have the option to select binding arbitration for the resolution of your complaint under certain circumstances, provided you have taken the following steps: (1) raised your complaint directly with LI-COR, Inc. and provided LI-COR, Inc. the opportunity to resolve the issue; (2) made use of the independent dispute resolution mechanism identified above; and (3) raised the issue through the relevant data protection authority and allowed the US Department of Commerce an opportunity to resolve the complaint at no cost to you. For more information on binding arbitration for the EU-US Privacy Shield, please see the US Department of Commerce's Privacy Shield Framework: Annex I (Binding Arbitration), available at: <https://www.privacyshield.gov/article?id=ANNEX-I-introduction>. Regarding Swiss Personal Data, the binding arbitration option is expected to be implemented during the first annual review of the Swiss-US Privacy Shield by the Department of Commerce and Swiss government. Please visit <https://www.privacyshield.gov> for more information on the availability of the binding arbitration option for your Swiss Personal Data.

## CATEGORIES OF INFORMATION COLLECTED

LI-COR is a global leader in the design, manufacture, and marketing of high quality, innovative instruments, software, reagents, and integrated systems for plant biology, biotechnology, drug discovery, and environmental research. LI-COR'S global headquarters (LI-COR, Inc.) is located in the US, with subsidiaries in the United Kingdom (LI-COR Ltd) and Germany (LI-COR GmbH). Business operations are supported worldwide through a network of distributors and sales agents, although LI-COR provides services to individuals directly in many countries. Privacy and protection of your Information is of the highest importance to LI-COR. For the purposes described below under the section titled Purposes for Processing Information, LI-COR collects the following Information:

1. Information voluntarily provided either electronically, verbally, or in writing to us from individuals including without limitation first and last name, position or title, company/institution, lab information, publication information, addresses, phone numbers, and email addresses. Examples of this information include information you provide by filling in forms on a LI-COR Website, such as information provided at the time of registering to use a LI-COR Website, subscribing to a service, ordering a product, and requesting information, support, or further services from us. We may also collect financial information from you, which may include for example debit or credit card information, banking information, or other billing information, for billing purposes.
2. Order transaction and fulfillment details, which may include for example orders you place through a LI-COR Website or with one of our service representatives. You may be required to provide financial information before placing an order through a LI-COR Website or otherwise. Through e-commerce features on a LI-COR Website, we collect a session ID for the customer that is related to your customer account identity that allows us to associate your order with your account.
3. Contact information for sales enablement and marketing available from publicly available sources.
4. Instrument, software, and product information, which may include for example serial numbers of instruments sold and in a customer's possession, instrument configuration, software download and instrument update history, factory service history, and support history and requests, including communications with customers.
5. Research focus and potential application interest gathered from publicly available sources, provided

voluntarily by individuals, or gathered through automated systems including automated decision making systems which monitor individuals using our web services and LI-COR Websites as described in this policy.

6. Information provided to us from our affiliates and third-party distributors or sales agents for the purpose of business operations, order fulfillment, service, support, marketing, and to otherwise support our business. We will obtain assurances from our affiliates, including our subsidiaries or parent entity, and our third-party distributors or sales agents that permission has been granted from the individual for the exchange of personal information to LI-COR where such permission is required by applicable law. If LI-COR has knowledge that personal information was disclosed to LI-COR without any required permission, LI-COR will take reasonable steps to prevent or stop the use or disclosure of this information immediately.
7. Automatically collected information from individuals using our web services and LI-COR Websites, including browser type and version, browser language, operating system, device manufacturer and model, device settings, unique device identifiers, crash data, network information, statistics on page views, traffic to and from our user interfaces, browsing activity including address of website visited prior to using the LI-COR Website or services, date and time the services were used, ad data, internet protocol (IP) address, language preferences, and cookie data, and inferred information about a customer including country, company, city, state, postal code, metropolitan area, and phone area code. This may include information about your online activities over time (behavioral tracking). This information may include personal information or we may maintain or associate it with personal information we collect in other ways or receive from third parties. Like many websites, LI-COR Websites and our web services may use a “cookie” identifier to obtain this automatic information. The cookies placed by the LI-COR server are readable only by LI-COR, and cookies cannot access, read, or modify any other data on a computer. This information is used to provide a personalized and relevant experience for our website users and to track general statistics on traffic. The “help” portion of the toolbar on most browsers will direct users on how to prevent the browser from accepting new cookies, how to configure the browser to tell the user when they receive a new cookie, or how to fully disable cookies. If you disable cookies,

some features of the LI-COR Website or our web services may not work.

8. Response to email communication including when an individual “opens” or “clicks”, if supported by the individual’s program.
9. LI-COR provides data management and processing software solutions through LI-COR Websites, web services, instruments, and downloaded software that allow users to enter or collect certain data to be processed by the solution at the direction of the user (“Customer Data”). Those who voluntarily use these solutions agree to the processing, transfer, and storage of such Customer Data. Any Customer Data that is EEA or Swiss Personal Data may be transferred to a server in the US according to this policy. LI-COR has taken steps to secure Customer Data. Customer Data is encrypted in transit, and passwords are encrypted at rest for the protection of our users. LI-COR only uses, accesses, processes, or discloses Customer Data in accordance with instructions provided by the user, the user’s consent, the applicable services agreement, and this policy or as required by applicable law, legal process, or regulation.
10. Information provided by customers, our subsidiaries or parent entity, distributors, or sales agents in order to provide technical support or service for instrumentation or other products or services or guidance on the use of the instrumentation or other products or services. This may include data collected or processed by the instrument user or information automatically collected by instruments, software, or other products and services for support and product and service improvement purposes.

Generally, you are not under a statutory or contractual requirement to provide any EU Personal Data to us. However, it may be necessary for you to provide certain EU Personal Data to us in order to purchase products or services from us or for us to provide certain services, such as warranty or repair services or web services, and without such EU Personal Data we may not be able to sell or provide such products or services to you. For example, we may need your address in order to return a repaired instrument to you or you may need to provide certain EU Personal Data to us when subscribing to our web service that allows us to validate that your license remains active in order for you to continue to use such service.

As noted above, we may receive some Information from third parties, including our subsidiaries, parent entity, distributors, and sales agents, as well as publicly available sources. We protect Information obtained from third parties according to the practices described in this policy, plus any

additional restrictions imposed by the source of the Information. We may combine this data with Information we collect.

## **PURPOSES FOR PROCESSING INFORMATION**

LI-COR collects, uses, and processes your Information for purposes of:

1. Supporting marketing, sales, ordering, and shipping of products and services by us or our third-party distributors or sales agents.
2. Supporting the service, support, provision, updating, improvement, and maintenance of our products and services and development of current and future products and services.
3. Providing you with the support and services you request, including for example providing customer support and processing Customer Data according to your instructions.
4. Communication with our subsidiaries or parent entity, third-party distributors, or sales agents, or with you, including without limitation for promotional communications. If you contact us, we may use your Information to respond.
5. Communication and advertising of LI-COR products and services.
6. For the operation of our business, including without limitation for billing, account management, and other administrative and operational purposes.
7. With regard to Website Information specifically:
  - a. Presenting a LI-COR Website and contents to you.
  - b. Allowing you to participate in interactive features on a LI-COR Website.
  - c. Notifying you about changes to a LI-COR Website or any products or services we offer or provide through it.
8. Carrying out our obligations and enforcing our rights arising from any contracts entered into between you and us, including without limitation for billing and collection.
9. Fulfilling any other purpose for which you provide it or for which you consent.
10. Investigating and helping prevent fraud and security issues and abuses.
11. As required by applicable law, legal process, or regulation.

If Information is aggregated or de-identified so it is no longer reasonably associated with an identified or

identifiable natural person, LI-COR may disclose or use it for any purpose as allowed by applicable law.

LI-COR will only process EU, EEA, and Swiss Personal Data in ways that are compatible with the purpose that LI-COR collected it for, or for purposes the individual later authorizes. Before we use your EU, EEA, or Swiss Personal Data for a purpose that is materially different than the purpose we collected it for or that you later authorized, we will provide you with the opportunity to opt out. You may also opt out from disclosures of your EU, EEA, or Swiss Personal Data to unaffiliated third-party data controllers. Opt-out requests can be sent to the contact listed under the Questions or Complaints section below. LI-COR maintains reasonable procedures to help ensure that EU, EEA, and Swiss Personal Data is reliable for its intended use, accurate, complete, and current.

Individuals may choose to opt out of correspondence or communication from LI-COR. Those who opt out will only receive communication necessary for safe and accurate operation of instruments in their possession or as a direct result of a product purchase or warranty registration. If we have sent you a promotional email, you may send us a return email asking to be omitted from future email distributions. Individuals may select their email opt-out preferences at [www.licor.com/email-preference](http://www.licor.com/email-preference) or by contacting LI-COR as set forth under the E-mail Opt-Out section below. Further, individuals may contact LI-COR as set forth below under Questions or Complaints and request to opt out of all correspondence and communication from LI-COR, subject to the limitations set forth above.

## **DATA TRANSFERS TO THIRD PARTIES**

To the extent allowed by applicable law, we may disclose Information to a third party: (i) as described in this section, (ii) with your consent, or (iii) where such third party is a buyer or other successor in the event of a merger, acquisition, divestiture, restructuring, reorganization, dissolution, or other sale or transfer of some or all of our assets or equity, whether as a going concern or as part of bankruptcy, liquidation, or similar proceeding, in which Information held by us is among the assets transferred.

You have the right to opt out of having Information disclosed to an unaffiliated third-party data controller or used for some reason other than that for which the data was originally collected. In order to opt out of such disclosures, please email [privacy@licor.com](mailto:privacy@licor.com) or contact LI-COR as provided in the Questions or Complaints section below.

## Subsidiaries / Parent

We may disclose your Information to our subsidiaries and parent entity. If LI-COR, Inc. transfers your EEA or Swiss Personal Data to one of its subsidiaries, LI-COR, Inc. will take steps to ensure that your EEA or Swiss Personal Data is protected with the same level of protection that the Privacy Shield requires.

## Third-Party Data Agents or Service Providers

LI-COR may disclose your Information to third-party service providers, distributors, and sales agents (“Third-Party Partners”) to support LI-COR’s business, including for example to support the sales and service of our products and services, the administration of our business, our global and domestic operations, logistics, and other functions and purposes related to LI-COR’s business, and to fulfill the purposes for which you provided the Information or any other purposes disclosed by us when you provided the Information. Functions performed on our behalf by Third-Party Partners include for example consulting, cloud storage services, customer information management software services, marketing software services, ecommerce software services, sales and support, and professional services. Third-Party Partners are not under the direct control of LI-COR, however LI-COR will take reasonable steps to ensure that these companies comply with Privacy Shield Principles (if applicable), applicable data protection law, and this policy. The privacy and control of Information is of the utmost concern to LI-COR. If LI-COR determines that a Third-Party Partner is using Information inconsistent with Privacy Shield Principles (if applicable), applicable data protection law, or this policy, LI-COR will immediately take steps to protect the Information, and prevent or stop any misuse or disclosure.

Where required by the Privacy Shield (if applicable) or other applicable data protection law, we enter into written agreements with Third-Party Partners requiring them to provide the same level of protection the Privacy Shield (if applicable) or other applicable data protection law requires and limiting their use of EU, EEA, and Swiss Personal Data to the specified services provided on our behalf. We take reasonable and appropriate steps to ensure that Third-Party Partners process EU, EEA, and Swiss Personal Data in accordance with our Privacy Shield obligations (if applicable) and other obligations under applicable data protection law and to stop and remediate any unauthorized processing. Under certain circumstances, we may remain liable for the acts of our Third-Party Partners who perform services on our behalf for their handling of EU, EEA, and Swiss Personal Data that we transfer to them.

## Third-Party Data Controllers

In some cases we may transfer your Information to unaffiliated third-party data controllers, which may include unaffiliated distributors and sales agents. These third parties do not act as our agents or service providers and are not performing functions on our behalf. We may transfer your Information to third-party data controllers for the purpose of processing the sale of our products and services and providing you with additional information about our products and any related services and support that are available from such third-party data controllers. We will only provide your EU, EEA, and Swiss Personal Data to unaffiliated third-party data controllers where you have not opted out of such disclosures, or in the case of sensitive EU, EEA, and Swiss Personal Data, where you have opted-in if the Privacy Shield or other applicable data protection law requires consent. We enter into written contracts with any unaffiliated third-party data controllers requiring them to provide the same level of protection for EU, EEA, and Swiss Personal Data the Privacy Shield or other applicable data protection law requires. We also limit their use of your EU, EEA, and Swiss Personal Data so that it is consistent with any consent you have provided and with the notices you have received.

## Disclosures for National Security or Law Enforcement

Under certain circumstances, we may be required to disclose your Information in response to valid requests by public authorities, including without limitation to meet national security or law enforcement requirements.

## Disclosures Required by Law

If we receive a request for your Information, we may disclose your Information if we reasonably believe disclosure is in accordance with or required by any applicable law, regulation, or legal process.

## Aggregated or De-identified Data

We may disclose aggregated or de-identified Information to any third party for any purpose as allowed by applicable law.

## Other Disclosures

As allowed by applicable law, we may share your Information with third parties to enforce our contracts or policies, investigate and prevent fraud or security issues, or otherwise protect and defend our rights, property, or safety or that of third parties.



## SECURITY OF INFORMATION

All Information collected and stored by LI-COR is maintained in a secure corporate data center and/or otherwise subject to reasonable and appropriate access protection. Further, LI-COR maintains reasonable and appropriate security measures to help protect your EU, EEA, and Swiss Personal Data from loss, misuse, unauthorized access, disclosure, alteration, or destruction in accordance with the Privacy Shield (if applicable) and other applicable data protection law. These measures include appropriate technical and organizational measures to ensure a level of security appropriate to the risk. While LI-COR strives to secure Information, even strict security measures do not guarantee the security of any information, and LI-COR cannot guarantee that the Information it seeks to protect will be protected in all circumstances, including those beyond its reasonable control.

LI-COR limits access to Information to only those persons in LI-COR who have legitimate and specific business reasons. Only employees who need Information to perform a specific job are granted access to it. Any employee granted access to Information is trained and kept up to date on our privacy and security practices. During training they are made aware of their responsibilities to protect the security, confidentiality, and integrity of Information. LI-COR will conduct regular and random audits of its privacy practices to verify adherence to this policy. Any employee that LI-COR determines is in violation of this policy will be subject to disciplinary action up to and including termination of employment.

## ACCESS

Subject to applicable exemptions, you may have the right to access the EEA or Swiss Personal Data that LI-COR, Inc. holds about you and to request that LI-COR, Inc. correct, amend, or delete it if it is inaccurate or processed in violation of the Privacy Shield. If you would like to request access to, correction, amendment, or deletion of your EEA or Swiss Personal Data held by LI-COR, Inc., you can submit a written request to LI-COR to the contact set forth below in the Questions or Complaints section. We may request specific information from you to confirm your identity.

## ADDITIONAL NOTICES FOR EU PERSONAL DATA ONLY

This section provides additional notices to you with respect to your EU Personal Data.

## DATA CONTROLLER

A data controller decides how your EU Personal Data is processed and for what purposes.

LI-COR, Inc. is a data controller of your EU Personal Data and may be a data processor of your EU Personal Data in some situations on behalf of its subsidiaries and a data processor of your EU Personal Data that is Customer Data.

Further, LI-COR Ltd and LI-COR GmbH are data controllers of your EU Personal Data with respect to operations outside of the US and may be data processors of your EU Personal Data in some situations on behalf of their affiliates and data processors of your EU Personal Data that is Customer Data.

Contact information for each entity is provided below under the Questions or Complaints section.

## LEGAL BASIS FOR PROCESSING

In order to process your EU Personal Data, including any sensitive EU Personal Data, we may rely on one or more of the following legal bases:

- consent to the processing;
- processing is necessary for the performance of a contract to which you are a party;
- processing is necessary in order to take steps at your request prior to entering into a contract;
- compliance with a legal obligation of the European Union or a member state of the European Union to which we are subject;
- processing is necessary to comply with your vital interests; and
- for the purposes of our legitimate interests.

Where processing is based on our legitimate interests, those legitimate interests include operating our business, offering and supporting our products and services, and promoting our products and services.

## YOUR RIGHTS

You may exercise any of the rights described in this section by submitting a written request to LI-COR to the contact set forth below in the Questions or Complaints section. We may request specific information from you to confirm your identity. Your rights with respect to your EU Personal Data may include, subject to any exemptions provided by applicable law:

- **Right to Access and Correction:** You may have the right to access your EU Personal Data that we hold about you and to request that we correct, amend, or delete it if it is inaccurate or processed in violation of the Privacy Shield (if applicable) or other applicable data protection law.
- **Right to Withdraw Consent:** Where you have provided your consent to the processing of your EU Personal Data by LI-COR you may withdraw your consent at any time by sending a communication to the LI-COR contact listed below specifying which consent you are withdrawing. Please note that the withdrawal of your consent does not affect the lawfulness of any processing activities based on such consent before its withdrawal.
- **Right to Restrict Use:** You may have the right to restrict our use of your EU Personal Data, including for example when you object to the processing of your EU Personal Data on the basis of legitimate interests or you object to the accuracy of your EU Personal Data we hold.
- **Right to Complain about Processing:** You may have the right to lodge a complaint about our processing of your EU Personal Data with a supervisory authority in the EU.
- **Right to Object to Processing:** Applicable data protection law may entitle you to require us not to process your EU Personal Data for certain specific purposes (including profiling or automated decision making) where such processing is based on our legitimate interests. If you object to such processing, pursuant to applicable data protection law we will no longer process your EU Personal Data for these purposes unless we can demonstrate compelling legitimate grounds for such processing or such processing is required for the establishment, exercise, or defense of legal claims. You further have an absolute right to object to our processing your EU Personal Data for direct marketing purposes, and you can exercise this right by submitting a written request to LI-COR to the contact set forth below in the Questions or Complaints section.
- **Right to be Forgotten:** You may be entitled in certain situations to request that we erase your EU Personal Data and close your accounts. Please note that if you request the erasure of your EU Personal Data, we may retain and use your EU Personal Data to the extent necessary to comply with our legal obligations or if required for the establishment, exercise, or defense of legal claims. For example, for our tax, legal reporting, and auditing obligations.
- **Right to Data Portability:** In some cases, you may have the right to require that we transmit your EU Personal Data to another data controller.

## HOW LONG WE RETAIN YOUR DATA

We keep your EU Personal Data for no longer than reasonably necessary for the purposes described herein, including without limitation providing our products and services and fulfilling the transactions you have requested, or for other essential purposes such as complying with our legal obligations, resolving disputes, and enforcing our agreements.

We use the following criteria to determine how long to retain your EU Personal Data:

- How long such data is reasonably needed to offer our products and services and operate our business, including without limitation improving and maintaining our products and services and maintaining our business and financial records. This is a key consideration underlying our data retention policies for your EU Personal Data.
- Whether our customer reasonably expects we will maintain that data. This will lead us to maintaining the data until the customer requests deletion.
- If the data is of a sensitive type, a shortened retention time is generally more appropriate.
- Whether we have announced a specific data retention period that would apply to such data.
- Whether we are subject to a legal, contractual, or other similar obligation to retain the data.

Because these needs can vary for different data types in the context of different products or services, actual retention periods can vary significantly.

## USE OF THIRD-PARTY WEB SERVICES

LI-COR may also allow certain third parties, such as Google, Inc. (“Google”) or Microsoft, Inc. (“Microsoft”) to place tracking technologies on LI-COR Websites using their web services. We use such web services as Google Analytics, Google Adwords, and Google Remarketing, all provided by Google, and Bing Ads Connect, provided by Microsoft, among others. LI-COR and these third-party vendors use cookies together to inform, optimize, and serve advertisements based on a user’s previous visit to a LI-COR Website. This information may be used to report how the use of these advertisements and the interactions with them are related to visits to LI-COR Websites. Data collected from this advertising may be used to provide a more personalized experience for users. The “help” portion of the toolbar on most browsers will direct users on how to prevent the browser from accepting new cookies, how to configure the browser to tell the user when they receive a new cookie, or how to fully disable cookies. If you disable cookies, some

features of the LI-COR Website or our web services may not work. For more information about targeting and advertising cookies and how you can opt out, you can visit the Network Advertising Initiative's opt-out page at <http://optout.networkadvertising.org/?c=1#!/>, the Digital Advertising Alliance's opt-out page at <http://optout.aboutads.info/?c=2#!/>, or <http://youonlinechoices.eu>. Please note we do not control any of these opt out links and are not responsible for their availability or accuracy.

Google Analytics uses cookies to help analyze how users use LI-COR Websites. Through Google Analytics, we collect information through your browser or device such as browser type and version, operating system, device manufacturer and model, and network. In most cases, information generated by the cookie is transferred to a server of Google in the US and stored there. LI-COR uses IP anonymization, where the IP address of individuals located within the EEA or Switzerland is truncated before being sent to a US-based server. Google Analytics ensures that the IP address of the user is not associated with any other data held by Google. Users may opt-out of Google Analytics by following the instructions and downloading the tool available at: <https://tools.google.com/dlpage/gaoptout> (you may have to have cookies enabled), or customize their Google advertisement settings at: <http://www.google.com/settings/ads>.

## **NOTICE TO END USERS**

Many of our products and services are intended for use by organizations and are administered to you by your organization. We are not responsible for the privacy or security practices of your organization, which may differ from those set forth in this policy. If your organization is administering your use of our products and services, please direct your privacy inquiries to your administrator. Your use of our products and services may be subject to your organization's policies, if any.

## **CHILDREN'S INFORMATION**

No LI-COR Website, instrument, software, or other service or product that we offer is intended for children under 16 years of age. No one under age 16 may provide any personal data to us without parental consent. We do not knowingly collect personal data from children under 16. If we learn we have collected or received personal data from a child under 16 without verification of parental consent, we will delete that information. If you believe we might have any information from or about a child under 16, please send an email to [privacy@licor.com](mailto:privacy@licor.com) or mail your request to the address set forth below under Questions or Complaints.

## **CHANGES TO THIS PRIVACY POLICY**

We reserve the right to amend this policy from time to time. Any such amendment that affects EU, EEA, and Swiss Personal Data will be consistent with the Privacy Shield's requirements and other applicable data protection law. Unless otherwise provided by applicable law, use of Information we collect is subject to the policy in effect at the time such Information is collected.

If we make changes to this policy, the revised policy will be posted on <http://www.licor.com>. If we make material changes to this policy regarding your personal data, we will provide additional notice, for example via e-mail, prominently posting a notice of such changes on LI-COR Websites, or through the affected services. If you disagree with the changes to this policy, you should cease using our products and services. The date this policy was last revised is identified below. Please periodically check this date to determine when the policy was last updated.

You are responsible for ensuring we have an up-to-date active and deliverable e-mail address for you and for periodically visiting <http://www.licor.com> and this policy to check for any changes.



## QUESTIONS OR COMPLAINTS

You can direct any questions or complaints about this policy or the use or disclosure of your Information to the following contact:

*Daniel E. Hile*  
*Chief Operating Officer*  
*LI-COR, Inc.*  
*4647 Superior Street*  
*Lincoln, NE 68504 USA*  
*privacy@licor.com*  
*+1-402-467-3576*

Individuals located outside of the US may also contact our EU affiliate as follows with any questions or complaints:

*Glen McMurdo*  
*European Managing Director*  
*LI-COR GmbH*  
*Siemensstrasse 25 A*  
*D-61352 Bad Homburg*  
*Germany*  
*+49 (0) 6172 17 17 771*

Individuals who wish to contact LI-COR Ltd should contact our European Managing Director at LI-COR GmbH at the above address.

Our data protection representative is the following:

*LI-COR GmbH*  
*Siemensstrasse 25 A*  
*D-61352 Bad Homburg*  
*Germany*  
*+49 (0) 6172 17 17 771*

## E-MAIL OPT-OUT

If you wish to change or remove your contact information from our e-mail database, you may select your e-mail opt-out preferences at [www.licor.com/email-preference](http://www.licor.com/email-preference), send an email to [privacy@licor.com](mailto:privacy@licor.com) or mail your request to:

*Director of Marketing Communications*  
*LI-COR Biosciences*  
*4647 Superior Street*  
*Lincoln, NE 68504*  
*USA*

Effective Date: May 25, 2018

Last Revised: Thursday, May 24, 2018



**LI-COR Biosciences**  
**Global Headquarters**

4647 Superior Street  
Lincoln, Nebraska 68504  
USA

© LI-COR, Inc. 2018  
977-14752 v7 • 05/18

LI-COR® is a registered trademark of LI-COR, Inc. All brand and product names are trademarks or registered trademarks of their respective owners.

The LI-COR board of directors would like to take this opportunity to return thanks to God for His merciful providence in allowing LI-COR to develop and commercialize products, through the collective effort of dedicated employees, that enable the examination of the wonders of His works.

“Trust in the LORD with all your heart and do not lean on your own understanding. In all your ways acknowledge Him, and He will make your paths straight.”