



DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

Privacy Policy

Effective Date:10/29/2018

Overview

Dojindo Molecular Technologies, Inc. Privacy Policy describes how Dojindo Molecular Technologies, Inc. treats your personal data provided to and collected by Dojindo Molecular Technologies, Inc., 30 West Gude Drive, Suite 260, Rockville, MD 20850 USA in accordance with this policy, as well as applicable laws, rules and, regulations. This Privacy Policy does not apply to third-party websites where the website of Dojindo Molecular Technologies may be linked. Personal data includes the information generally obtained such as your name, address, phone number, email, IP address, name of organization you belong to, history of purchased products and inquiry of our products and services.

1. The Purpose of Data Collection

The collected data is used to deliver our products and provide services you requested. We may also use collected data to analyze our marketing activities and improve our services.

2. The Way of Data Collection

Information given by you

We may collect the data when you decide to provide it to us, such as by communication through phone, fax, email, and meeting, or entering your information on a contact form or questionnaire.

Information automatically collected

In some cases, data (such as IP Address, browser information, operating system, web server logs) is collected automatically by IT systems (including tracking technologies such as Cookies and Google Analytics) when you visit the website. Cookies are necessary to communicate through the system to provide better services. For more information about the IT system, please review [Appendix](#) attached.



DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

3. Data Sharing and Transfer

We do not share, transfer or disclose your personal data to third parties without your express consent unless we are required by public authorities under the law.

4. Your Right for Your Personal Data

You have the right to your personal data in accordance with data protection regulations if the exercise of your request is legally accepted, such as;

- to request or access information about your stored data
- to correct or/and delete your data
- to transfer your data (when you request the direct transfer of data to another responsible party, "data portability")
- to object or/and restrict to the processing of data
- to withdraw your consent to the use of your data
- to lodge complaint with a supervisory authority

If you have any questions or comments, please contact us anytime at

Dojindo Molecular Technologies, Inc.
30 West Gude Drive, Suite 260
Rockville, MD 20850 USA

Telephone: 301-987-2667

Email: info@dojindo.com

5. The Period Your Data is Kept

We store data as long as it is legally necessary or necessary to fulfill the service requested by you, or as long as it has been agreed upon in a declaration of consent. We erase your personal data in compliance with data protection regulations according to your request anytime if not prohibited by law or contract.



DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

6. Data Security

The site uses SSL/TLS for security, in order to protect the transmission of confidential content including your inquiry sent to us.

7. Use of this Website by Children

We do not knowingly collect or use any personal data directly from children under the age of 16 unless the prior, verifiable consent of their legal representative is given. Upon request, their legal representatives have the right to view the information provided by the children and/or to request deletion of such information.

8. Our Privacy Policy Change

Dojindo Molecular Technologies may revise Privacy Policy from time to time, and any such changes will be immediately reflected on our website with a new effective date. Any personal data collected will be handled in accordance with the updated Privacy Policy.



DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

Appendix: General Information

What is Cookie?

Cookies are used to remember the information such as for logging in websites if the log-in process is required, recording which pages were accessed, or items in the shopping cart in an online store or to record the user's browsing activity. For example, when you visit a website's login page, the server of the website sends a cookie, small piece of data, and your computer stores it. When you log in, the server remembers the cookie and allow you to access the webpages. Cookies can be used to remember information about you in order to show you relevant content that you accessed several times. Many websites use cookies for providing information you may be interested in based on your preferences. You select your preferences by entering your information in a web form and submitting the form to the server. The server encodes the preferences in a cookie and sends the cookie back to the browser. This way, every time you access a page on the website, the server can personalize the page according to your preferences.

Type of Cookies

There are several cookies used for certain purposes. Session cookie is used for navigation of webpages and be deleted by closing the browser. Persistent cookies are used as ways to compile records of browsing history. The lifespan of persistent cookie/ tracking cookie is set arbitrarily. These cookies are reset if the expiration times is reached or you delete the cookie manually. Secure cookie can only be transmitted by an encrypted connection, HTTPS. Third-party cookies are used to build up a browsing history of the user across all the websites that have advertisements from the advertiser. Currently, we are using session cookies and persistent cookies such as to maintain the website ("Necessary Cookie") or to optimize the website according to your choice ("Functionality Cookie") or for the analytical purpose ("Analytical Cookie") * on our website. There are several websites are available to learn more about cookies. Followings are examples of the websites describe cookies.

*Necessary Cookie is needed for the operation of the website and to give you all the necessary functions (e.g., to identify trusted web traffic)

*Functionality Cookie is used to record information about choices you have made, in order to



DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

customize the website for you.

*Analytical Cookie is used for internal purposes such as the better understanding of how visitors use our website. This helps us to make an improvement on our website and offer better experience on our website.

For more information, you can visit websites below,

- <https://developer.mozilla.org/en-US/docs/Web/HTTP/Cookies>
- <https://computer.howstuffworks.com/cookie.htm>
- <https://www.epic.org/privacy/internet/cookies/>
- <https://support.mozilla.org/en-US/kb/cookies-information-websites-store-on-your-computer>

Disclaimer: We are not responsible for the information provided from the websites indicated above or you may need to seek for additional information.

Use of Google Analytics

This website uses Google Analytics, a web analytics service provided by Google Inc. ("Google"). Google Analytics uses "cookies" which allow an analysis of users' browsing habits. The information generated by the cookie about your use of the website is usually transmitted to a Google server in the USA and stored there.

IP anonymization

We have also extended the code "anonymizeIP" on this website to Google Analytics. This guarantees the masking of your IP address so that all data is collected anonymously.

How to Prevent/Disable Cookies or Opt-out from Automatic Data Collection

Prevent or Disable Cookies

You can prevent the storage of cookies and collecting data by a corresponding setting of your browser software. However, some pages may not be accessible or do not work properly without necessary cookies.

DOJINDO
DOJINDO MOLECULAR TECHNOLOGIES, INC.

*30 West Gude Drive, Suite 260, Rockville, MD 20850
Phone: 301-987-2667 Fax: 301-987-2687*

The opt-out cookies (disable cookies) are only valid in this browser you are currently using and only for our website and are stored on your device.

For suggestions on how to disable cookies, please visit following websites.

<https://www.wikihow.com/Disable-Cookies>

<https://support.google.com/accounts/answer/61416?co=GENIE.Platform%3DDesktop&hl=en>

Google Analytics Opt-out

If you do not want your data used by Google Analytics, you can opt-out (prevent it) by downloading and installing the Google Analytics browser plugin available at the following link:

<https://tools.google.com/dlpage/gaoptout?hl=en>.

Disclaimer: We are not responsible for the information provided from the websites indicated above or you may need to seek for additional information.