

Dojindo EU GmbH Privacy Policy

Effective from 01. 09. 2021

Introduction

Dojindo EU Privacy Policy describes how Dojindo EU GmbH treats your personal data provided to and collected by Dojindo EU GmbH, Leopoldstr. 254, 80807 Munich, Germany (“we”, “us”, “our”, “Dojindo”) 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 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.



What is covered by this privacy policy

This document covers our processing of your personal data, online via our website, as well as offline in case you contact us. The website are as follows:

<https://www.dojindo.eu.com/>

Legal grounds of our processing

We will only process your personal data if there is a legal basis to do so. Under GDPR, there are 4 obligations that we use.

1. Consent – If you agree to our processing. This will be clear when we ask for it. You are also able to withdraw your consent at any time.
2. Contractual obligation  – If we have a contract with you, we will need to store and use some information about you.
3. Legal obligation § – If there is a law forcing us to process your personal data, we have to comply.
4. Balance of interests (Legitimate interest)  – Legitimate interest allows us to process personal data, if it is necessary for our legitimate interests and your interests or fundamental rights and freedoms do not outweigh our interest. We will inform below if we do such processing and what interests we pursue. You can also find information about your rights

We have marked all processing activities described below with these symbols, so that you can easily see and understand what our legal grounds are.

Deletion of your personal data §

We will generally delete your personal information as soon as we no longer require it for the purposes for which we collected it. If any legal obligation stops us from deleting your information for a certain time, we will stop access to this information, store it securely and delete it after our legal obligation is over.

Steuernummer/Tax number:143/130/21379, VAT ID: DE 281981280

Amtsgericht/District court: München (HRB 197296) Geschäftsführer/Managing director: Mr. Mutsuo Yoshida

Bankverbindung/Bank information: MUFG Bank (Europe) N.V. Germany Branch

Kontonr/Account number: 222961 BLZ: 300 107 00 BIC:BOTKDE33 IBAN: DE15 3001 0700 0000 2229 61

In case of processing based on your consent, we will delete information as soon as you revoke your consent or when the agreed upon processing has finished.

This Website

When using our website, certain personal information will be transferred to our servers automatically. This is necessary to provide you with the website. For example, our servers need to know your IP-Address to know where to send information.

Informational use

If you simply use our website to get information, we will automatically receive and store the following information about you:

Content data

- Any text you enter or files you upload

Usage data

- Websites visited
- Date and time of your request

Metadata

- Your IP-Address
- Your computers time zone
- Your computers date and time
- The website you requested
- Which website you came from
- Information about your browser (version, language, fonts installed)
- Information about your operating system (version, language)
- Potentially other information your browser sends via the HTTP protocol.

This information is necessary for us to present you our website and show all our products and services. We have a legitimate interest to present our company with a modern looking website. As we only collect necessary information, there is no technical way for us to limit the kinds of information we receive.

We use an external hosting provider to serve the website. The provider receives and stores the content data, usage data and metadata on our behalf.

We use Cloudflare as our CDN-provider to load some pictures and elements of this website as well as protect our website from DDoS attacks and malicious use. The CDN-provider will receive usage data and metadata of your visit. By using a CDN, we can more efficiently provide you with the information you requested and keep our website secure. The legal basis for this use is our legitimate interest. Cloudflare is a service offered by Cloudflare Inc., 101 Townsend St, San Francisco, CA 94107, USA. You can review their privacy policy here: <https://www.cloudflare.com/privacypolicy>

Contact

If you contact us, we will receive and store the information you have entered additional to the information described above.

Steuernummer/Tax number:143/130/21379, VAT ID: DE 281981280

Amtsgericht/District court: München (HRB 197296) Geschäftsführer/Managing director: Mr. Mutsuo Yoshida

Bankverbindung/Bank information: MUFGBank (Europe) N.V. Germany Branch

Kontonr/Account number: 222961 BLZ: 300 107 00 BIC:BOTKDE33 IBAN: DE15 3001 0700 0000 2229 61

We will use the information you have entered to respond to your message and take any actions you may have requested.

We do not use the information for other purposes, unless you agree with it, and delete it as soon as we have completed your request, unless a legal obligation forces us to keep the information for longer.

If you contact us, we usually receive the following information from you:

- E-Mail: We need this to respond to you
- Name: We use this to personalize our message and to see if you had contacted us before
- Company: We use this to personalize our message and to see if your company had contact with us before
- Your message

In some cases, we will not be able to respond to your message if information is missing.

Our basis for this processing is our contractual obligation if your message concerns purchasing or services offered by Dojindo.

If you have a general request, our legal basis is legitimate interest, as we want to give you the best service we can.

Cookies on this website / 

Cookies are small text files that any website can write into your browser. They only contain text and cannot contain malware. Cookies may be set by ourselves (first party cookies), or by third parties. If we set a cookie in your browser, other websites cannot read it. If other third parties set a cookie in your browser, any other website which uses the same third party can read the information stored in these cookies.

This allows for tracking functionality across websites. If we use third party cookies, we will always ask for your consent before placing these cookies.

We use Cookiebot as our consent management platform (CMP) to manage your preferences on cookie selection. Cookiebot is a service offered by Cybot, Havnegade 39, 1058 Copenhagen, Denmark. You can review their privacy policy here: <https://www.cybot.com/privacy-policy/>

Below you can see a list of all first party cookies which we set:

Name	Content	Purpose	Duration
_ga	Unique User ID	Analysing your use of the website; Targeting advertisements to you	2 years
_gid	Unique User ID	Analysing your use of the website; Targeting advertisements to you	1 day
_gat	Unique User ID	Analysing your use of the website; Targeting advertisements to you	1 minute
SecureSessionKey	Unique Code	Verifying and storing your preferences and session	3 months

SessionKey	Unique Code	Verifying and storing your preferences and session	3 months
SessionID	Unique Session ID & Date and Time of Access	Verifying and storing your preferences and session	3 months
__cfduid	Trust token	Allowing your access to the site after an anti-bot check has been successfully run	1 month
ASP.NET_SessionId	Unique Session ID	Unused	Session only
cookieconsent_status	Cookie choices	Remembering your cookie choices	1 year

Third Parties

Tools /

On our websites, we may embed tools and contents loaded from the servers of third party service providers. Examples of this may be graphics, videos, social-media buttons, or other functionalities. Whenever you access a website which embeds such contents, your IP-address will be transmitted to these third parties.

Occasionally these tools might include so called web beacons, invisible graphics that are used to gather further statistics. If you notice any such web beacon, please notify us and we will strive to remove it from our websites.

Some third-party contents will also set cookies in your browser to store certain information on your usage of the tools and contents.

Occasionally, the third party might connect your information gathered on our websites to other information they have collected in order to build a profile on you.

Google Analytics

This website uses Google Analytics, a service for web analysis provided by Google Ireland Ltd., Gordon House, Barrow Street, Dublin 4, Ireland ("Google"). Google Analytics uses cookies to enable analysis of your usage of our site.

Google uses the information to analyse your usage of our site, to compile reports on website-activities and to provide further services in relation to internet usage and website usage to us. It is also used to determine your interests and serve relevant advertisements to you.

You can prevent the collection of cookie-generated data in relation to your usage of this site (including your IP-address) by Google, as well as the processing of such data by downloading and installing the opt-out plugin from the following address: <https://tools.google.com/dlpage/gaoptout>

We use Google Analytics with the extension “_anonymizeIp()”. IP-addresses are only processed after being shortened on servers within the EU/EEA, making it impossible to relate the collected data to any individual. Only in rare exemptions the full IP-address will be transmitted to the US and shortened there.

Steuernummer/Tax number:143/130/21379, VAT ID: DE 281981280

Amtsgericht/District court: München (HRB 197296) Geschäftsführer/Managing director: Mr. Mutsuo Yoshida

Bankverbindung/Bank information: MUFG Bank (Europe) N.V. Germany Branch

Kontonr./Account number: 222961 BLZ: 300 107 00 BIC:BOTKDE33 IBAN: DE15 3001 0700 0000 2229 61

We use Google Analytics to learn more about the usage of our website in order to continuously improve it. By referencing the statistical reports generated, we can make our offerings more interesting to our users.

The legal grounds for this use of Google Analytics is your active consent which you have given when visiting our website.

You can read more about Google Analytics and the Google data protection below:

Terms of use: <https://www.google.com/analytics/terms/en.html>

Overview of privacy: <https://support.google.com/analytics/answer/6004245>

Privacy Policy: <https://policies.google.com/privacy>

Fonts

We use external fonts on our website to create a unified experience and deliver a modern looking website. These fonts are loaded from third parties to speed up loading processes in your browser. Our legal ground is our legitimate interest to provide our website with fonts securely, efficiently and maintenance free.

We use the Raleway font from Google.

When you load our website, Google will learn which website you accessed and what your IP-Address is. The information collected will only be used to show the embedded fonts as well as keep statistics on the popularity of fonts.

For more information please reference the following links:

FAQ:

https://developers.google.com/fonts/faq#what_does_using_the_google_fonts_api_mean_for_the_privacy_of_my_users

We also use Font Awesome, a service provided by Fonticons, Inc., 307 S. Main St., Suite 202, Bentonville, AR 72712, USA (<https://fontawesome.com>) to consistently present fonts and symbols, including emoticons. You can access their privacy policy here: <https://fontawesome.com/privacy>. Font Awesome will collect information on which icon is downloaded and when, in order to optimize the content delivery networks, find technical errors, defend against attacks, compile statistics on icon popularity and compile statistics on the kinds of software and computers used.

YouTube

On some of our webpages, we embed videos from the platform YouTube, offered by Google Ireland Ltd., Gordon House, Barrow Street, Dublin 4, Ireland ("Google"). When you access a webpage where a video is embedded, Google will become aware of your visit to the website, including your IP-Address, date and time of your visit and video that is embedded. To store additional information about your interaction with the video (for example, when you paused), Google will store a cookie in your browser.

Our legal grounds for embedding videos is our legitimate interest in the modern communication and promotion of our products.

To change your setting for ads and turning of personalized ads, please follow the link below:
<https://adssettings.google.com/authenticated>

Your Rights

You have the right of access regarding your personal data. You always have the right to ask us for copies of your personal information, as well as additional information about our processing. In some cases, we may ask for additional verification of your identity to prevent fraud.

You also have the right to rectification or erasure or to restriction of processing of your personal data, insofar no legal limits apply.

You have the right to object our data processing based on our legitimate interest on grounds arising from your special circumstances. We ask you to detail these grounds when you object so that we may evaluate if we still have a compelling legitimate interest that overrides your interests, rights, and freedoms.

You also have may have the right to data portability. This right applies to information you have provided to us. You have the right to ask that we transfer the information you gave us to another organization or give it to you. This right only applies if we process your information on grounds of your consent, or when the processing is based on your contract with us.

You also have the right to lodge a complaint against our processing of your data with a supervisory authority. A list of all German supervisory authorities can be accessed here:

https://www.bfdi.bund.de/DE/Infothek/Anschriften_Links/anschriften_links-node.html

We are based in Bavaria, so the “Bayerisches Landesamt für Datenschutzaufsicht”

(<https://www.lida.bayern.de/de/kontakt.html>) is competent for handling complaints against us.

Contact us

You can contact us for any questions regarding the privacy policy or our use of your personal data via the contact information below.

Dojindo EU GmbH
Leopoldstr.254
80807, Munich, Germany
Tel +49 89 3540 4805
Fax +49 89 3540 4806
Email: info@dojindo.eu.com