

Quanterix Corporation Data Privacy Framework Statement

This Data Privacy Framework Statement ("Privacy Framework Statement") describes how Quanterix Corporation (the "Company," "we," or "us") collect, use, and disclose certain personally identifiable information that we receive in the US from the European Union ("EU"), the United Kingdom ("UK") and Switzerland ("Personal Data").

This Privacy Framework Statement supplements our [Website Privacy Policy](#) and our Employee/Consultant Privacy Policy (together, the "Company Policies"), and unless specifically defined in this Privacy Framework Statement, the terms in this Privacy Framework Statement have the same meaning as the Company Policies.

We comply with the EU-U.S. Data Privacy Framework (EU-U.S. DPF) and the UK Extension to the EU-U.S. DPF, and the Swiss-U.S. Data Privacy Framework (Swiss-U.S. DPF) as set forth by the U.S. Department of Commerce. We have certified to the U.S. Department of Commerce that we adhere to the EU-U.S. Data Privacy Framework Principles (EU-U.S. DPF Principles) with regard to the processing of personal data received from the EU and the UK in reliance on the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF. We have certified to the U.S. Department of Commerce that we adhere to the Swiss-U.S. Data Privacy Framework Principles (Swiss-U.S. DPF Principles) with regard to the processing of personal data received from Switzerland in reliance on the Swiss-U.S. DPF. If there is any conflict between the terms in this Privacy Framework Statement and the EU-U.S. DPF Principles and/or the Swiss-U.S. DPF Principles, the Principles shall govern. To learn more about the Data Privacy Framework (DPF) Program, and to view our certification, please visit <https://www.dataprivacyframework.gov/>.

For purposes of enforcing compliance with the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF and the Swiss-U.S. DPF we are subject to the investigatory and enforcement authority of the US Federal Trade Commission (FTC).

Personal Data Collection and Use

Our Company Policies describe the categories of Personal Data that we may receive in the US as well as the purposes for which we use that Personal Data. The Company will only process Personal Data in ways that are compatible with the original purpose for which it was collected, or for purposes the individual later authorizes. Before we use your 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. The Company maintains reasonable procedures to help ensure that Personal Data is reliable for its intended use, accurate, complete, and current.

We may collect the following categories of sensitive Personal Data: certain demographic information that may include information regarding gender, nationality, and/or racial origin. When we collect sensitive Personal Data, we will obtain your express opt-in consent where the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF or the Swiss-U.S. DPF requires, including if we disclose your sensitive Personal Data to third parties, or before we use your sensitive Personal Data for a different purpose than we collected it for or than you later authorized.

Data Transfers to Third Parties

Third-Party Agents or Service Providers. We may transfer Personal Data to our third-party agents or service providers who perform functions on our behalf or on behalf of our customers as described in our Company Policies. Where required by the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF or the Swiss-U.S. DPF, we enter into written agreements with those third-party agents, resellers, licensees, and service providers requiring them to provide the same level of protection the DPF Program requires and limiting their use of the data to the specified services provided on our behalf or on behalf of our customers. We take reasonable and appropriate steps to ensure that third-party agents and service providers process Personal Data in accordance with our DPF Program obligations and to stop and remediate any unauthorized processing. Under certain

circumstances, we may remain liable for the acts of our third-party agents, resellers, licensees, or service providers who perform services on our behalf or on behalf of our customers for their handling of Personal Data that we transfer to them.

Disclosures for National Security or Law Enforcement

Under certain circumstances, we may be required to disclose your Personal Data in response to valid requests by public authorities, including to meet national security or law enforcement requirements.

Security

We maintain reasonable and appropriate security measures to protect Personal Data from loss, misuse, unauthorized access, disclosure, alteration, or destruction in accordance with the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF or the Swiss-U.S. DPF.

Access Rights

You may have the right to access the 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 EU-U.S. DPF and the UK Extension to the EU-U.S. DPF or the Swiss-U.S. DPF. These access rights may not apply in some cases, including where providing access is unreasonably burdensome or expensive under the circumstances, or where it would violate the rights of someone other than the individual requesting access. If you would like to request access to, correction, amendment, or deletion of your Personal Data, you can submit a written request to the contact information provided below. We may request specific information from you to confirm your identity. In some circumstances we may charge a reasonable fee for access to your information. If we have entered into a contractual agreement with a customer to administer assessments, we will refer your access request to our customer for response.

Questions or Complaints

In compliance with the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF and the Swiss-U.S. DPF, we commit to resolve DPF Principles-related complaints about our collection and use of your Personal Data. EU and UK individuals and Swiss individuals with inquiries or complaints regarding our handling of personal data received in reliance on the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF, and the Swiss-U.S. DPF should first contact us at:

Quanterix Corporation
900 Middlesex Turnpike, Building 1
Billerica, MA 01821
ATTN: Privacy
Phone: 617-301-9400
Email: privacy@quanterix.com

We will investigate and attempt to resolve any complaints or disputes regarding the use or disclosure of your Personal Data within 45 days of receiving your complaint.

In compliance with the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF and the Swiss-U.S. DPF, we commit to refer unresolved complaints concerning our handling of personal data received in reliance on the EU-U.S. DPF and the UK Extension to the EU-U.S. DPF and the Swiss-U.S. DPF to JAMS, an alternative dispute resolution provider based in the United States. If you do not receive timely acknowledgment of your DPF Principles-related complaint from us, or if we have not addressed your DPF Principles-related complaint to your satisfaction, please visit the [JAMS website](#) for more information or to file a complaint. The services of JAMS are provided at no cost to you.

Binding Arbitration

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 the Company and provided us 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, see [Annex I to the Data Privacy Framework](#).

Contact Us

If you have any questions about this Privacy Framework Statement or would like to request access to your Personal Data, please contact us as follows:

Quanterix Corporation
900 Middlesex Turnpike, Building 1
Billerica, MA 01821
ATTN: Privacy
Phone: 617-301-9400
Email: privacy@quanterix.com

Changes to This Policy

We reserve the right to amend this Privacy Framework Statement from time to time consistent with the DPF Program's requirements.

Effective Date: October 11, 2023