



**Salesforce's Processor Binding Corporate Rules  
for the  
Processing of Personal Data**

# Salesforce Processor BCR

## Table of Contents

<b><i>1. Introduction</i></b>	<b><i>4</i></b>
<b><i>2. Definitions</i></b>	<b><i>4</i></b>
<b><i>3. Scope and Application</i></b>	<b><i>5</i></b>
<b><i>4. Responsibilities Towards Customers</i></b>	<b><i>6</i></b>
<b>A. General Obligations</b>	<b>6</b>
<b>B. Transparency, Fairness, Lawfulness and Cooperation with Customers</b>	<b>6</b>
<b>C. Data Subject Rights</b>	<b>6</b>
i. Data Subject Requests	6
ii. Handling of Complaints	7
<b>D. Regulatory Inquiries and Complaints</b>	<b>7</b>
<b>E. Data Protection Impact Assessments</b>	<b>7</b>
<b>F. Records of Processing Activities</b>	<b>7</b>
<b><i>5. Description of Processing Operations and Transfers</i></b>	<b><i>7</i></b>
<b>A. Purpose Limitation</b>	<b>7</b>
<b>B. Nature of Personal Data Processed</b>	<b>8</b>
<b>C. Affected Data Subjects</b>	<b>8</b>
<b>D. Countries of location of the Salesforce Group Affiliate Sub-processors</b>	<b>9</b>
<b>E. Data Quality</b>	<b>9</b>
<b>F. Sub-processing</b>	<b>9</b>
i. Sub-processing Within the Salesforce Group	9
ii. Sub-processing by Third Parties	9
iii. Notification of New Sub-processors and Objection Rights	10
<b><i>6. Confidentiality and Security Measures</i></b>	<b><i>10</i></b>
<b>A. Confidentiality and Training</b>	<b>10</b>
<b>B. Data Security</b>	<b>10</b>
<b>C. Personal Data Incident Management and Notification</b>	<b>10</b>
<b>D. Audits</b>	<b>11</b>
i. Third-Party Audits and Certifications	11

## Salesforce Processor BCR

ii. Network of Privacy Personnel and Internal Verification	11
iii. Customer Audits	12
<b>7. <i>Third-Party Beneficiary Rights</i></b>	<b>12</b>
<b>A. Rights directly enforceable against the Salesforce Group</b>	<b>13</b>
<b>B. Rights enforceable against the Salesforce Group where the Data Subject is not able to bring a claim against the Customer acting as Controller</b>	<b>13</b>
<b>C. Modalities</b>	<b>14</b>
<b>8. <i>Liability and Enforcement</i></b>	<b>14</b>
<b>9. <i>Cooperation with Supervisory Authorities</i></b>	<b>15</b>
<b>10. <i>Local Law Requirements</i></b>	<b>15</b>
<b>11. <i>Data Protection Officer</i></b>	<b>16</b>
<b>12. <i>Salesforce Processor BCR and Applicable Law</i></b>	<b>16</b>
<b><i>Appendix A – Services to which the Salesforce Processor BCR applies</i></b>	<b>17</b>

# Salesforce Processor BCR

## 1. Introduction

Salesforce, Inc. and its affiliates are committed to achieving and maintaining customer trust. Integral to this mission is providing a robust security and privacy program that carefully considers data protection matters.

In accordance with the General Data Protection Regulation and, as applicable, the Swiss Data Protection Laws and Regulations, the Salesforce Processor BCR (as defined below) is intended to provide an adequate level of protection for Personal Data during international transfers within the Salesforce Group made on behalf of Customers and under their instructions.<sup>1</sup>

## 2. Definitions

- **Controller** means the entity which determines the purposes and the means of the processing of Personal Data.
- **Customer(s)** means (i) a legal entity with whom a member of the Salesforce Group has executed a contract to provide the Services (or a legal entity placing an order under such contract) and such contract incorporates by reference the Salesforce Processor BCR or (ii) a legal entity with whom a member of the Salesforce Group has executed a contract under which the legal entity is entitled to resell the Services to its end customers and such contract incorporates by reference the Salesforce Processor BCR.
- **Data Subject** means the identified or identifiable person to whom Personal Data relates.
- **General Data Protection Regulation** or **GDPR** means European Regulation 2016/679 of the European Parliament and of the Council of 27 April 2016 on the protection of natural persons with regard to the processing of personal data and on the free movement of such data and repealing European Directive 95/46/EC.
- **Personal Data** means any information relating to (i) an identified or identifiable natural person; and (ii) an identified or identifiable legal entity (where such information is protected similarly as personal data or personally identifiable information under Swiss Data Protection Laws and Regulations).
- **Processor** means the entity which processes Personal Data on behalf of the Controller.
- **Salesforce Group** means Salesforce, Inc. and its affiliate Sub-processors of Personal Data, listed as a 'Salesforce Affiliate' in the Infrastructure and Sub-processor documentation for each Service covered by the Salesforce Processor BCR, available [here](#).

---

<sup>1</sup> For clarity, a Customer (as defined in Section 2) may be a Controller or a Processor of Personal Data. Where a Customer is a Processor of Personal Data, the Salesforce Group shall process Personal Data as Sub-processors on behalf of the Controller. Instructions from the Controller regarding the processing Personal Data shall be given through the Processor.

## Salesforce Processor BCR

- **Salesforce Processor BCR** means Salesforce's Processor Binding Corporate Rules for the Processing of Personal Data, the most current version of which is available on Salesforce Group's website, currently located [here](#).
- **Services** means the online services provided to Customer by the Salesforce Group, as listed in **Appendix A**.
- **Sub-processor** means any Processor engaged by a member of the Salesforce Group.
- **Supervisory Authority** means an independent public authority which is established by an EU member state pursuant to Article 51 of the GDPR, and/or, as applicable, the Swiss data protection authority established under the Swiss Data Protection Laws and Regulations.
- **Swiss Data Protection Laws and Regulations** means the Swiss Federal Data Protection Act 1992 and its successor laws.

### 3. Scope and Application

The purpose of the Salesforce Processor BCR is to govern international transfers of Personal Data to and between members of the Salesforce Group, and to third-party Sub-processors (in accordance with written agreements with any such third-party Sub-processors) when acting as Processors and/or Sub-processors on behalf and under the documented instructions of Customers.

The Salesforce Processor BCR applies to Personal Data submitted to the Services by:

- Customers established in an EEA member state or Switzerland whose processing activities for the relevant data are governed by the GDPR or, as applicable, by the Swiss Data Protection Laws and Regulations; or
- Customers established in non-EEA member states for which the customer has contractually specified that the GDPR and implementing national legislation shall apply.

The Salesforce Group may update the Salesforce Processor BCR with approval from the Salesforce Group's appointed privacy leader, general counsel and compliance officer. All changes to the Salesforce Processor BCR shall be communicated to members of the Salesforce Group.

The Salesforce Group's appointed privacy leader shall be responsible for keeping a fully updated list of the members of the Salesforce Group and third-party Sub-processors and making appropriate notifications to Customers and the French data protection authority ("CNIL") in its capacity as competent Supervisory Authority for the Salesforce Processor BCR. The Salesforce Group shall not transfer Personal Data to a new member of the Salesforce Group until such member is appropriately bound by and complies with the Salesforce Processor BCR.

The Salesforce Group shall make the most current version of the Salesforce Processor BCR, including the members of the Salesforce Group, available [here](#). Significant changes to the Salesforce Processor BCR and/or the list of members of the Salesforce Group will be reported (a) in a timely fashion to Customers and (b) once per year to the relevant Supervisory Authorities via the CNIL in its capacity as competent Supervisory Authority for the Salesforce Processor BCR accompanied by a brief explanation of the changes.

## **Salesforce Processor BCR**

When the changes to the Salesforce Processor BCR affect the processing conditions, the Salesforce Group shall inform the Customer in such a timely fashion that Customer has the possibility to object to the change or to terminate the contract before the modification is made.

The categories of Personal Data, the types of processing and its purposes, the types of Data Subjects affected and the identification of the recipients in the third countries are set out in Section 5 below.

It shall be the responsibility of a Customer to apply the Salesforce Processor BCR to:

- All Personal Data processed for processor activities and that are submitted to EU and, as applicable, Swiss law; or
- All processing of Personal Data for processor activities within the Salesforce Group whatever the origin of the data.

### **4. Responsibilities Towards Customers**

#### **A. General Obligations**

The Salesforce Group and its employees shall comply with the Salesforce Processor BCR, process Personal Data only upon a Customer's documented instruction and shall have a duty to respect Customer's instructions regarding the data processing and the security and confidentiality of Personal Data, pursuant to the measures provided in the contracts executed with Customers.

The Salesforce Group shall immediately inform the Customer if in its opinion an instruction infringes the GDPR or other EU or EU member state law or, as applicable, Swiss data protection provisions.

#### **B. Transparency, Fairness, Lawfulness and Cooperation with Customers**

The Salesforce Group undertakes to be transparent regarding its Personal Data processing activities and to provide Customers with reasonable cooperation and assistance within a reasonable period of time to help facilitate their respective data protection obligations regarding Personal Data, to the extent Customer, in its use of the Services, does not have the reasonable ability to address such obligations.

#### **C. Data Subject Rights**

Members of the Salesforce Group act as Processors on behalf of Customers. As between the Salesforce Group and Customers, Customers have the primary responsibility for interacting with Data Subjects, and the role of the Salesforce Group is generally limited to assisting Customers as needed.

##### **i. Data Subject Requests**

The Salesforce Group shall promptly notify Customer if the Salesforce Group receives a request from a Data Subject to exercise the Data Subject's right of access, right to rectification, restriction of processing, erasure ("right to be forgotten"), data portability, object to the processing, or its right not to be subject to an automated individual decision making ("Data Subject Request"). Taking into account the nature of the processing, the Salesforce Group shall assist Customer by appropriate technical and organizational measures, insofar as this is possible, for the fulfilment of Customer's obligation to respond to a Data Subject Request under the GDPR or, as applicable, an equivalent obligation under Swiss Data Protection Laws and Regulations. In addition, to the extent Customer, in its use of the Services, does not have the ability to address a Data Subject Request, the Salesforce Group shall upon Customer's request provide commercially reasonable efforts to assist Customer in responding to such Data Subject Request, to the

## **Salesforce Processor BCR**

extent the response to such Data Subject Request is required under the GDPR or, as applicable, an equivalent obligation under Swiss Data Protection Laws and Regulations. To the extent legally permitted, Customer shall be responsible for any costs arising from the Salesforce Group's provision of such assistance.

### **ii. Handling of Complaints**

The Salesforce Group's Privacy department shall be responsible for handling complaints related to compliance with the Salesforce Processor BCR.

Data Subjects may lodge a complaint about processing of their respective Personal Data that is incompatible with the Salesforce Processor BCR by contacting the relevant Customer or the Salesforce Group's Privacy department at the email address [privacy@salesforce.com](mailto:privacy@salesforce.com). The Salesforce Group shall without undue delay communicate the complaint to the Customer to whom the Personal Data relates without obligation to handle it (except if it has been agreed otherwise with Customer).

Customers shall be responsible for responding to all Data Subject complaints forwarded by the Salesforce Group except in cases where a Customer has disappeared factually or has ceased to exist in law or become insolvent. Where the Salesforce Group is aware of such a case, it undertakes to respond directly to Data Subjects' complaints within one (1) month, including the consequences of the complaint and further actions Data Subjects may take if they are unsatisfied by the reply (such as lodging a complaint before the relevant Supervisory Authority). Taking into account the complexity and number of requests, this period of one (1) month can be extended by two (2) further months in which case the Salesforce Group will inform the Data Subjects accordingly.

### **D. Regulatory Inquiries and Complaints**

The Salesforce Group shall, to the extent legally permitted, promptly notify a Customer if the Salesforce Group receives an inquiry or complaint from a Supervisory Authority in which that Customer is specifically named. Upon a Customer's request, the Salesforce Group shall provide the Customer with cooperation and assistance in a reasonable period of time and to the extent reasonably possible in relation to any regulatory inquiry or complaint involving the Salesforce Group's processing of Personal Data.

### **E. Data Protection Impact Assessments**

Upon Customer's request, the Salesforce Group shall provide Customer with reasonable cooperation and assistance needed to fulfil Customer's obligation under the GDPR (or, as applicable, under the Swiss Data Protection Laws and Regulations) to carry out a data protection impact assessment related to Customer's use of the Services, to the extent Customer does not otherwise have access to the relevant information, and to the extent such information is available to the Salesforce Group. The Salesforce Group shall provide reasonable assistance to Customer in the cooperation or prior consultation with the Supervisory Authority in the performance of its tasks relating to this Section 4 E. of the Salesforce Processor BCR to the extent required under the GDPR (or, as applicable, under the Swiss Data Protection Laws and Regulations).

### **F. Records of Processing Activities**

As required by data protection laws and regulations, the Salesforce Group shall maintain a record of all categories of processing activities carried out on behalf of each Customer.

### 5. Description of Processing Operations and Transfers

#### A. Purpose Limitation

The Salesforce Group shall only process Personal Data on behalf of and in accordance with Customer's documented instructions for the following purposes: (i) processing in accordance with a Customer's instructions set forth in the Customer's contract with a member of the Salesforce Group including with regard to transfers of personal data to a third country (unless the Salesforce Group is legally required to do so by EU or EU member state law, or, as applicable, Swiss law, in which case prior information will be provided by the Salesforce Group to Customer unless such information is legally prohibited); and (ii) processing initiated by the Customer in its use of the Services. If the Salesforce Group cannot comply with such purpose limitation, a member of the Salesforce Group shall promptly notify the relevant Customer, and such Customer shall be entitled to suspend the transfer of Personal Data and/or terminate the applicable order form(s) in respect to only those Services which cannot be provided by the Salesforce Group in accordance with such Customer's instructions. On the termination of the provision of such Services, the Salesforce Group and third-party Sub-processors shall, at the choice of the Customer, return the Personal Data to the Customer and/or delete the Personal Data as set forth in the applicable customer contract and upon request from Customer, the Salesforce Group shall certify that it has done so. The only exception to this is if the law applicable to the Salesforce Group and its third-party Sub-processors requires the Salesforce Group and its third-party Sub-processors to retain the data that has been transferred in which case the Salesforce Group will inform the Customer and warrant that it will guarantee the confidentiality of the Personal Data transferred and will not actively process the Personal Data transferred anymore.

#### B. Nature of Personal Data Processed

The Salesforce Processor BCR will apply to Personal Data submitted by Customers to the Services. The Salesforce Group's Customers determine what Personal Data, if any, is submitted to the Services under the conditions set out in the contract.

The following types of Personal Data are oftentimes submitted to the Services.

- First and last name
- Title
- Position
- Employer
- Contact information (company, email, phone, physical business address)
- ID data
- Professional life data
- Personal life data
- Location data

These types of Personal Data oftentimes relate to the following categories of data subjects:

- Prospects, customers, business partners and vendors of Customer (who are natural persons)
- Employees or contact persons of Customer's prospects, customers, business partners and vendors
- Employees, agents, advisors, freelancers of Customer (who are natural persons)
- Customer's users authorized by Customer to use the Services



## **Salesforce Processor BCR**

Customers are allowed to submit sensitive personal data or special categories of Personal Data to some Services under the conditions set out in the contract.

### **C. Affected Data Subjects**

The Salesforce Group does not choose or determine the categories of Data Subjects that relate to the Personal Data submitted to the Services. The Salesforce Group's Customers solely determine the Data Subjects whose Personal Data is submitted to the Services.

### **D. Countries of location of the Salesforce Group Affiliate Sub-processors**

The countries where the Salesforce affiliate Sub-processors of Personal Data are located are listed in the Infrastructure and Sub-processor documentation for each Service covered by the Salesforce Processor BCR, available [here](#).

### **E. Data Quality**

Customers have access to, and control of, Personal Data in their use of the Services. To the extent a Customer, in its use of the Services, does not have the ability to anonymize, correct, amend, update or delete Personal Data, as required by applicable laws, the Salesforce Group shall comply with any request by a Customer in a reasonable period of time and to the extent reasonably possible to facilitate such actions by executing any measures necessary to comply with the law, in a reasonable period of time and to the extent reasonably possible to the extent the Salesforce Group is legally permitted to do so. The Salesforce Group will, to the extent reasonably required for this purpose, inform each member of the Salesforce Group to whom the Personal Data may be stored of any anonymization, rectification, amendment, update or deletion of such data. If any such anonymization, correction, amendment, update or deletion request is applicable to a third-party Sub-processor's processing of Personal Data, the Salesforce Group shall communicate such request to the applicable third-party Sub-processor(s).

### **F. Sub-processing**

#### **i. Sub-processing Within the Salesforce Group**

As set forth in applicable contracts with Customers, members of the Salesforce Group may be retained as Sub-processors of Personal Data, and depending on the location of the Salesforce Group member, processing of Personal Data by such Sub-processors may involve transfers of Personal Data. The Salesforce Processor BCR extends to all members of the Salesforce Group. The Salesforce Processor BCR is incorporated by reference into the Salesforce Group's code of conduct which is available [here](#).

#### **ii. Sub-processing by Third Parties**

As set forth in applicable contracts with Customers, members of the Salesforce Group may retain third-party Sub-processors, and depending on the location of the third-party Sub-processor, processing of Personal Data by such Sub-processors may involve transfers of Personal Data. Such third-party Sub-processors shall process Personal Data only: (i) in accordance with the Customer's instructions set forth in the Customer's contract with a member of the Salesforce Group; or (ii) if processing is initiated by the Customer in its use of the Services. The current list of third-party Sub-processors engaged in processing Personal Data, including a description of their processing activities, is available in the Infrastructure and Sub-processor documentation for each Service covered by the Salesforce Processor

## **Salesforce Processor BCR**

BCR, available [here](#). Such third-party Sub-processors have entered into written agreements with a member of the Salesforce Group in accordance with the applicable requirements of Articles 28, 29, 32, 45, 46 and 47 of the GDPR, or, as applicable, corresponding provisions of the Swiss Data Protection Laws and Regulations, as well as the relevant sections of the Salesforce Processor BCR as applicable to the third-party Sub-processor's processing activities.

### **iii. Notification of New Sub-processors and Objection Rights**

As set forth in applicable contracts with Customers, the Salesforce Group shall provide Customers with prior notification before a new Sub-processor begins processing Personal Data. Within thirty (30) days of receiving such notice, a Customer may object to Salesforce Group's use of a new Sub-processor by notifying the Salesforce Group in accordance with the provisions set forth in the Customer's contract. In the event Customer objects to a new Sub-processor, as permitted in the preceding sentence, the Salesforce Group will use reasonable efforts to make available to Customer a change in the Services or recommend a commercially reasonable change to Customer's configuration or use of the Services to avoid processing of Personal Data by the objected-to new Sub-processor without unreasonably burdening the Customer. If the Salesforce Group is unable to make available such change within a reasonable period of time, which shall not exceed sixty (60) days, Customer may terminate the applicable order form(s) with respect only to those Services which cannot be provided by the Salesforce Group without the use of the objected-to new Sub-processor by providing written notice to the Salesforce Group. The Salesforce Group will refund Customer any prepaid fees covering the remainder of the term of such order form(s) following the effective date of termination with respect to such terminated Services, without imposing a penalty for such termination on Customer.

## **6. Confidentiality and Security Measures**

### **A. Confidentiality and Training**

The Salesforce Group shall ensure that its personnel engaged in the processing of Personal Data are informed of the confidential nature of the Personal Data, have executed written confidentiality agreements and have received appropriate training on their responsibilities. Additionally, the Salesforce Group shall ensure that its personnel responsible for the development of the tools used to process Personal Data have received appropriate training on their responsibilities. The Salesforce Group shall also ensure that its personnel engaged in the processing of Personal Data are limited to those personnel who require such access to perform the Salesforce Group's obligations under applicable contracts with Customers.

### **B. Data Security**

The Salesforce Group shall maintain appropriate administrative, technical and physical measures for protection of the security (including protection against unauthorized or unlawful processing and against accidental or unlawful destruction, loss or alteration or damage, unauthorized disclosure of, or access to, Personal Data), confidentiality and integrity of Personal Data, as set forth in applicable contracts with Customers. The Salesforce Group shall implement technical and organizational measures which at least meet the requirements of the GDPR or, as applicable, Swiss Data Protection Laws and Regulations and any existing particular measure specified in the contract with the Customer. The Salesforce Group regularly monitors compliance with these measures. The Salesforce Group will not materially decrease the overall security of the Services during a Customer's applicable subscription term.

## Salesforce Processor BCR

### C. Personal Data Incident Management and Notification

In the event a member of the Salesforce Group becomes aware of the accidental or unlawful destruction, loss, alteration, unauthorized disclosure of, or access to Personal Data, transmitted, stored or otherwise processed by the Salesforce Group or its Sub-processors (a “Personal Data Incident”) the Salesforce Group will without undue delay after becoming aware notify affected Customers. The Salesforce Group shall make reasonable efforts to identify the cause of such Personal Data Incident and take those steps as the Salesforce Group deems necessary and reasonable in order to remediate the cause of such a Personal Data Incident to the extent the remediation is within the Salesforce Group’s reasonable control. The obligations herein shall not apply to incidents that are caused by Customer or Customer’s users.

### D. Audits

The Salesforce Group shall maintain an audit program to help ensure compliance with the Salesforce Processor BCR, including the following third-party audits and certifications, internal verification and audits by Customers. The audit program covers all aspects of the Salesforce Processor BCR, including methods for ensuring non-compliance is addressed.

#### i. Third-Party Audits and Certifications

The Security, Privacy and Architecture Documentation, available [here](#) describes the third-party audits and certifications applicable to each Service. The scope of these third-party audits and certifications is set forth in the corresponding audit reports and certificates which the Salesforce Group shall make available to its Customers upon request. Where the Salesforce Group has obtained the following third-party audits and certifications, the Salesforce Group agrees to maintain these, or their successors.

- **ISO 27001 certification:** The Salesforce Group is subject to an information security management system (ISMS) in accordance with the ISO 27001 international standard. Members of the Salesforce Group have achieved ISO 27001 certification for their ISMS from an independent third party.
- **SSAE 18 Service Organization Control (SOC) reports:** The Salesforce Group’s information security control environment applicable to the Services undergoes an independent evaluation in the form of SSAE 18 Service Organization Control (SOC) reports, which are available to Customers upon request.

#### ii. Network of Privacy Personnel and Internal Verification

The Salesforce Group has appointed a network of privacy personnel responsible for overseeing and ensuring compliance with the Salesforce Group’s data protection responsibilities at a local and global level, including compliance with this Salesforce Processor BCR, advising management on data protection matters, liaising with data protection authorities, and handling data protection-related complaints. Each member of the Salesforce Group shall be assigned such a member of network of privacy personnel. Such privacy personnel are primarily responsible for privacy-related matters and report to the Salesforce Group’s appointed privacy leader (who reports to the Salesforce Group’s general counsel) and benefit from the support of the Salesforce Group’s senior management. The Salesforce Group’s appointed privacy leader is responsible for the Salesforce Group’s compliance with applicable privacy and data protection laws and leads the Salesforce Group’s network of privacy personnel. The Salesforce Group’s network of privacy personnel have regional responsibility for the Salesforce Group’s compliance with applicable privacy and data protection laws.

## **Salesforce Processor BCR**

The Salesforce Group's compliance department shall conduct an annual assessment of the Salesforce Group's compliance with the Salesforce Processor BCR, which is provided to the Salesforce Group's appointed privacy leader, compliance officer and Salesforce, Inc.'s board of directors. Such an assessment shall include any necessary corrective actions, timeframes for completing such corrective actions, and follow up by Salesforce's compliance department to ensure such corrective actions have been completed.

The Supervisory Authority competent for the Customer may upon request have access to the results of such annual assessment and may carry out a data protection audit of any member of the Salesforce Group, if required.

### **iii. Customer Audits**

Upon a Customer's written request, and subject to appropriate confidentiality obligations, the Salesforce Group shall make available to the Customer (or such Customer's independent, third-party auditor that is not a competitor of the Salesforce Group) information regarding the Salesforce Group's and third-party Sub-processors' compliance with the data protection controls set forth in this Salesforce Processor BCR.

With respect to the Salesforce Group's compliance with the data protection controls set forth in the Salesforce Processor BCR, the Salesforce Group shall make available third-party certifications and audits set forth in the contract to the extent Salesforce makes them generally available to its customers.

With respect to third-party Sub-processors' compliance with the data protection controls set forth in the Salesforce Processor BCR, the Salesforce Group shall provide the requesting Customer a report of the Salesforce Group's audits of third-party Sub-processors and/or a report of third party auditors' audits of third-party Sub-processors that will have been provided by those third-party Sub-processors to the Salesforce Group.

Customer acknowledges and agrees to exercise its audit right by hereby instructing the Salesforce Group and the Salesforce Group's third party Sub-processors to carry out the audit as described in this Section 6.D (iii).

Customer has the right to change at any moment its instruction regarding the exercise of its audit right by sending the relevant member of the Salesforce Group a notice in writing.

If Customer changes its instruction and thereby requests to exercise its audit right directly, Customer shall reimburse the Salesforce Group for any time expended by the Salesforce Group or its third-party Sub-processors for any on-site audit carried out by the Customer. Before any such on-site audit commences, the requesting Customer and the Salesforce Group or its third party Sub-processors as appropriate shall mutually agree upon the scope, timing, and duration of the audit in addition to the reimbursement rate for which the Customer shall be responsible. All reimbursement rates shall be reasonable, taking into account the resources expended by the Salesforce Group or its third-party Sub-processors.

As set forth in applicable contracts with Customers, a Customer who performs an audit in accordance with this Section must promptly provide the Salesforce Group with information regarding any non-compliance discovered during the course of an audit.

Nothing in this Section affects any Supervisory Authority's or Data Subject's rights under the Salesforce Processor BCR.

### 7. Third-Party Beneficiary Rights

#### A. Rights directly enforceable against the Salesforce Group

Data Subjects may directly enforce the following elements of the Salesforce Processor BCR against the Salesforce Group as third party beneficiaries:

- a. Duty to respect the instructions from the Customer acting as Controller regarding the Data Processing including for data transfers to third countries located outside the EEA;
- b. Duty to implement appropriate technical and organizational security measures and duty to notify any security breach to the Customer acting as Controller;
- c. Duty to respect the conditions when engaging a Sub-processor either within or outside the Salesforce Group;
- d. Duty to cooperate with and assist the Customer acting as Controller in complying and demonstrating compliance with the law such as for answering requests from Data Subjects in relation to their rights;
- e. Provide an easy access to the Salesforce Processor BCR;
- f. Right to complain through internal complaint mechanisms;
- g. Duty to cooperate with the Supervisory Authority;
- h. Liability, compensation and jurisdiction provisions; and
- i. National legislation preventing respect of the Salesforce Processor BCR.

#### B. Rights enforceable against the Salesforce Group where the Data Subject is not able to bring a claim against the Customer acting as Controller

Data Subjects may directly enforce against the Salesforce Group the following elements of the Salesforce Processor BCR as third-party beneficiaries in those limited situations where a Data Subject is unable to bring a claim against the relevant Customer because such Customer has factually disappeared or ceased to exist in law or become insolvent unless a successor entity has been appointed to assume the legal obligations of the Customer by contract or by operation of law:

- Duty to respect the Salesforce Processor BCR;
- Creation of third party beneficiary rights for Data Subjects;
- Liability of salesforce.com France S.A.S for paying compensation and to remedy breaches to the Salesforce Processor BCR;
- Burden of proof on salesforce.com France S.A.S to demonstrate that the member of the Salesforce Group outside of the EU or the external Sub-processor is not liable for any violation of the rules which has resulted in the Data Subject claiming damages;
- Easy access for the Data Subjects to access the Salesforce Processor BCR and in particular information about their third party beneficiary rights and on the means to exercise those rights;
- Existence of a complaint handling process for the Salesforce Processor BCR;
- Duty for the Salesforce Group to cooperate with the Supervisory Authorities;
- Duty for the Salesforce Group to cooperate with the Controller;
- Description of the privacy principles;
- List of entities bound by the Salesforce Processor BCR; and
- Transparency requirement where national legislation prevents the Salesforce Group from complying with the Salesforce Processor BCR.

## Salesforce Processor BCR

### C. Modalities

The Data Subjects' rights as mentioned under sections A and B above shall cover the judicial remedies for any breach of the third party beneficiary rights guaranteed and the right to obtain redress and where appropriate, receive compensation for any damage.

In particular, Data Subjects shall be entitled to lodge a complaint before:

- the competent Supervisory Authority (with a choice for the Data Subject to choose between the Supervisory Authority of the EU Member State of his/her habitual residence, place of work or place of alleged infringement, or, if the Data Subject is based in Switzerland, the Swiss Federal Data Protection and Information Commissioner; and
- the competent court of the EU Member State (with a choice for the Data Subject to act before the courts where the Customer or the Salesforce Group has an establishment or where the Data Subject has his or her habitual residence, or, if the Data Subject is based in Switzerland, the competent Swiss court.

Where the Salesforce Group and the Customer involved in the same processing are found responsible for any damage caused by such processing, the Data Subject shall be entitled to receive compensation for the entire damage directly from the Salesforce Group.

The Salesforce Processor BCR are made available to Data Subjects [here](#).

### 8. Liability and Enforcement

Salesforce's contracts with Customers shall include a reference to the Salesforce Processor BCR and the Salesforce Processor BCR shall form part of those contracts. These contracts shall comply with Article 28 of the GDPR.

In accordance with such contracts, Customers shall have the right to enforce the Salesforce Processor BCR against any member of the Salesforce Group, for breaches they caused including judicial remedies and the right to receive compensation. Moreover, Customers shall have the right to enforce the Salesforce Processor BCR against salesforce.com France S.A.S. in case of: (i) a breach of the Salesforce Processor BCR or of the contract by members of the Salesforce Group established outside of the EU; or (ii) a breach by external Sub-processors established outside the EU of their sub-processing agreement with the Salesforce Group.

salesforce.com France S.A.S accepts responsibility for and agrees to take the necessary actions to remedy the acts of other members of the Salesforce Group established outside of the EU and third-party Sub-processors for breaches of the Salesforce Processor BCR or breaches caused by third-party Sub-processors established outside the EU and to pay compensation for any damages resulting from a violation of the Salesforce Processor BCR.

salesforce.com France S.A.S accepts liability as if the violation had taken place by salesforce France S.A.S in France instead of the member of the Salesforce Group outside of the EU or the third party Sub-processor established outside the EU. Salesforce France S.A.S may not rely on a breach by a Sub-processor (internal or external to the Salesforce Group) of its obligations in order to avoid its own liabilities.

With regard to Data Subjects, salesforce.com France S.A.S has the burden of proof to demonstrate that the member of the Salesforce Group outside of the EU or the third party Sub-processor is not liable for any violation of the rules which has resulted in the Data Subject claiming damages.

## **Salesforce Processor BCR**

With respect to Customer, to the extent a Customer can demonstrate that Customer has suffered damages and establishes facts showing that it is likely that such damages have occurred because of the Salesforce Group's breach of the Salesforce Processor BCR salesforce.com France S.A.S. shall be responsible for providing that the Salesforce member outside of the EU – or the third-party Sub-processor – was not responsible for the breach of the Salesforce Processor BCR giving rise to the damages or that no such breach took place.

If salesforce.com France S.A.S. can prove that the Salesforce member outside of the EU is not responsible for the act leading to the damages suffered by Customer or the Data Subject, salesforce.com France S.A.S. may discharge itself from any responsibility.

### **9. Cooperation with Supervisory Authorities**

The Salesforce Group shall cooperate with Supervisory Authorities with jurisdiction over the Salesforce Group or competent for Customers, reply to any requests they make within a reasonable time frame and abide by the advice and recommendations of the relevant EU member state regarding the interpretation and application of the Salesforce Processor BCR.

Upon request and subject to duties of confidentiality, the Salesforce Group shall provide relevant EU member state Supervisory Authorities with jurisdiction over the Salesforce Group or competent for Customers: (i) a copy of the Salesforce Group's annual assessment of compliance with the Salesforce Processor BCR and/or other documentation reasonably requested; and (ii) the ability to conduct an onsite audit of the Salesforce Group's architecture, systems and procedures relevant to the protection of Personal Data.

### **10. Local Law Requirements**

As set forth in applicable contracts with Customers, the Salesforce Group shall comply with applicable law in its processing of Personal Data. Where applicable law requires a higher level of protection for Personal Data than provided for in the Salesforce Processor BCR, the local applicable law shall take precedence.

Where the Salesforce Group reasonably believes that applicable existing or future enacted or enforceable law prevents it from fulfilling its obligations under the Salesforce Processor BCR or the instructions of a Customer, it shall promptly notify the Salesforce Group's Privacy department in addition to affected Customers, the Supervisory Authority competent for the Customer and the Supervisory Authority competent for Salesforce. In such a case, the Salesforce Group shall use reasonable efforts to make available to the affected Customers a change in the Services or recommend a commercially reasonable change to the Customers' configuration or use of the Services to facilitate compliance with applicable law without unreasonably burdening Customers. If the Salesforce Group is unable to make available such change within a reasonable period of time, Customers may terminate the applicable order form(s) in respect to only those Services which cannot be provided by the Salesforce Group in accordance with applicable law by providing written notice to the member of the Salesforce Group with whom the Customer has contracted. Such Customer shall receive a refund of any prepaid fees for the period following the effective date of termination for such terminated Services.

In accordance with applicable contracts with Customers, the Salesforce Group shall communicate any legally binding request for disclosure of Personal Data by a law enforcement authority or state security

## **Salesforce Processor BCR**

body to the impacted Customer unless the Salesforce Group is prohibited by law from providing such notification.

To the extent the Salesforce Group is prohibited by law from providing such notification, the Salesforce Group shall: (i) review each request on a case-by-case basis; (ii) use best efforts to request that the confidentiality requirement be waived to enable the Salesforce Group to notify the appropriate Supervisory Authority competent for the Customer and the CNIL in its capacity as competent Supervisory Authority for the Salesforce Processor BCR; and (iii) maintain evidence of any such attempt to have a confidentiality requirement waived

On an annual basis, the Salesforce Group shall provide the appropriate Supervisory Authorities competent for impacted Customers and the CNIL with general information about the types of legally binding requests for disclosure of Personal Data the Salesforce Group receives by law enforcement authorities.

Transfers of Personal Data by the Salesforce Group to any public authority cannot be massive, disproportionate and indiscriminate in a manner that would go beyond what is necessary in a democratic society.

### **11. Data Protection Officer**

The Salesforce Group has appointed a data protection officer in accordance with its legal obligation for the Salesforce Group who can be reached at [privacy@salesforce.com](mailto:privacy@salesforce.com)

### **12. Salesforce Processor BCR and Applicable Law**

Where national law applicable to the Salesforce Group requires a higher level of protection for Personal Data than what is set out in the Salesforce Processor BCR, then that national applicable law will take precedence over the Salesforce Processor BCR.

In any event the Salesforce Group shall process Personal Data in accordance with the national law applicable to the Salesforce Group.



## **Salesforce Processor BCR**

### **Appendix A – Services to which the Salesforce Processor BCR applies**

The Salesforce Processor BCR applies to the services branded as the following:

- Accounting Subledger
- Admissions Connect
- Advertising Studio (formerly branded as Social.com and Active Audiences)
- Audience Studio and Data Studio (formerly branded as Krux or Salesforce DMP)
- Automotive Cloud
- B2B Commerce and B2B Commerce on Lightning Experience (formerly branded as CloudCraze)
- B2C Commerce (formerly branded as Commerce Cloud or Demandware)
- Chatter
- Consumer Goods Cloud
- Customer Data Cloud (formerly branded as Salesforce Data Cloud)
- Customer Data Platform (formerly branded as CDP formerly branded as Customer 360 Audiences)
- Customer 360 Data Manager
- Datorama
- Datorama Reports for Marketing Cloud
- Digital Process Automation
- D2C Commerce (formerly branded as Salesforce Commerce for B2C and B2B2C Commerce)
- Education Cloud
- Einstein Bots
- Einstein Conversation Insights
- Einstein Copy Insights
- Einstein Discovery Classic (formerly branded as Einstein Discovery and BeyondCore)
- Einstein Engagement Scoring
- Einstein Prediction Builder
- Einstein Relationship Insights
- Einstein Vision and Language
- Einstein Vision for Social Studio
- Emergency Program Management
- Employee Productivity
- Enablement
- Enhanced Messaging
- Evergage (services branded or sold as Evergage, Data Science Workbench, and Data Warehouse)
- ExactTarget
- Experience Cloud (formerly branded as Community Cloud)
- Financial Services Cloud
- foundationConnect
- Government Cloud Plus
- Grants Management
- Headless Browser Service
- Health Cloud
- Heroku
- High Velocity Sales
- Intelligent Form Reader (IFR)
- Interaction Studio

## Salesforce Processor BCR

- IoT Cloud
- IoT Explorer
- Lightning Platform (including Force.com)
- LiveMessage
- Loyalty Management
- Manufacturing Cloud
- Marketing Cloud Einstein (formerly branded as Predictive Intelligence)
- Marketing Cloud for Nonprofits
- Messaging
- Messaging for In-App and Web
- Microsoft Teams Integration
- MuleSoft
- Net Zero Cloud (formerly branded as Sustainability Cloud)
- Non-Profit Cloud
- Nonprofit Cloud Case Management
- Non-Profit Cloud for Grantmaking
- Order Management
- Pardot and Pardot Einstein
- Privacy Center
- Public Sector Solutions
- Safety Cloud
- Sales Cloud and Sales Cloud Einstein
- Sales Enablement (formerly branded as myTrailhead)
- Salesforce Anywhere (formerly branded as Quip)
- Salesforce Connect
- Salesforce Contracts
- Salesforce CPQ and Salesforce Billing (together formerly branded as Salesforce Quote to Cash)
- Salesforce Inbox
- Salesforce Maps (Map Anything)
- Salesforce Order Management
- Salesforce Private Connect
- Salesforce Sales Planning
- Salesforce Slack Integration Proxy
- Salesforce Web3
- Salesforce.org Elevate
- Salesforce.org Insights Platform: Data Integrity
- Service Cloud and Service Cloud Einstein
- Service Cloud Voice
- Shift Management
- Site.com
- Slack
- Social Studio
- Student Success Hub (including the former Salesforce Advisor Link).
- Subscription Management
- Tableau CRM (formerly branded as Einstein Analytics, Analytics Cloud or Wave Analytics)
- Tableau Online
- Vlocity Services comprising Vlocity Communications, Vlocity Media and Entertainment, Vlocity Energy & Utilities, Vlocity Insurance, Vlocity Health, Vlocity Public Sector, Communications Cloud, Energy & Utility Cloud, and Media Cloud (all formerly branded as Vlocity Managed Packages)

## **Salesforce Processor BCR**

- WDC
- Workplace Command Center
- Workforce Engagement Management (WEM)