

SARTORIUS

Speak-Up Policy

Simplifying Progress

1. Purpose of the Speak-Up Policy	3
2. Scope of the Speak-Up Policy.....	3
3. Main Principles	4
3.1. Protection from Retaliation and Prohibition of Misuse.....	4
3.2. Confidentiality and Anonymity.....	5
3.3. Data Protection.....	6
(a) Informing Individuals Subject to Speak-Up Reports.....	6
(b) Retention of Records.....	6
4. Procedure for Submitting Reports and Follow-Up	7
4.1. Internal Reporting Channels and Submitting Reports.....	7
(a) Whistleblower Portal.....	7
(b) Telephone Hotline.....	7
(c) Email Inbox.....	7
(d) Surface Post.....	7
(e) In Person	8
(f) Country-specific Channels.....	8
4.2. External Reporting Channels	8
5. Handling of Complaints.....	8
6. Questions and concerns	9
7. Revision History.....	9
Annex 1: Internal Local Reporting Channels (country specific).....	10
Annex 2: External Reporting Channels.....	12

1. Purpose of the Speak-Up Policy

Sartorius AG and each company in which Sartorius AG directly or indirectly has effective control, i.e., holds a majority of shares or has management control ("**Sartorius**"), are subject to various laws and regulations as well as internal compliance policies worldwide. In addition, Sartorius has an ethical responsibility towards its customers, employees and business partners.

If applicable laws and regulations are violated, this can lead to considerable risks for Sartorius and its employees. It is therefore not only Sartorius's culture to ensure that all employees behave in accordance with applicable laws and regulations, but also to ensure that compliance violations are identified, investigated and prevented in the future.

Anyone inside or outside Sartorius (including, but not limited to, employees of direct or indirect suppliers) may file a complaint. Sartorius' employees as well as third parties such as suppliers, customers and other partners have access to various complaint channels.

Any indications of possible violations of laws and/or applicable regulations (e.g. the Sartorius Code of Conduct for Employees or the Code of Conduct for Business Partners) or of human rights or environmental risks or violations can be reported through the whistleblowing and complaints procedure.

To this end, Sartorius makes an effort to establish an open and reliable working environment where questions, doubts and concerns can be raised without fear of retaliation. As part of this effort, Sartorius issues this Speak-Up Policy ("**Policy**").

Any mandatory local laws prevail over this Policy.

2. Scope of the Speak-Up Policy

This Policy applies directly without further transposition to all employees of Sartorius, i.e. each and every one working at and for the company, including trainees, volunteers, officers, as well as members of the executive bodies, including any management board member or supervisory board member, each on group and local level. Also, this Policy applies to persons having self-employed status working for and behalf of Sartorius, as well as applicants and persons whose contracts of engagement have already ended or have not commenced yet. All persons referred to in this section are defined as "Employees" in the sense of this Policy.

This Policy is intended to encourage and enable all Employees to report established or soundly suspected breaches of applicable laws, standards and regulations and internal policies and guidelines that they become aware of in the

course of their employment. The reports can relate to various areas, such as environmental and human rights, bribery, corruption, data protection, ethical misconduct, antitrust law.

Sartorius expects Employees to report in particular the following in line with this Policy:

- Breaches of applicable national and international laws and regulations, in particular with regard to corruption, bribery, fraud, danger to health and safety (including excessive working hours), conflicts of interests, human rights violations (including labour rights, discrimination, harassment etc.), antitrust and unfair competition, tax and books and records requirements, money laundering and pollution of the environment (water, air, soil, etc.);
- In connection with human rights or the environment: (Alleged) Incidents in Sartorius' own business area as well as in the supply chain, e.g., health and safety in the workplace / depletion of natural resources / prohibition of child labor, forced labor and all forms of slavery / disregard of freedom of association, freedom of organization and the right to collective bargaining;
- Breaches of internal policies and regulations, in particular our Code of Conduct and Anti-Corruption Code; and
- Concerns regarding safety and quality of the products and services Sartorius is providing.

This Policy is applicable in addition to local speak-up policies to the extent in place. Employees can decide which option they want to use to raise their concerns.

3. Main Principles

3.1. Protection from Retaliation and Prohibition of Misuse

Sartorius firmly believes that effective reporting structures and an open and transparent culture help protect the Company, its employees, customers and business partners from damage and reputational harm.

Hence, Sartorius is committed to a policy of non-retaliation. Therefore, Employees and any third parties who are awarded protection by applicable laws do not have to fear any negative consequences of discriminatory or disciplinary nature for making a report or raising questions, doubts and concerns. Sartorius does not tolerate any kind of threats, detrimental behaviour or acts of retaliation. Such conduct is itself a breach of our values and can result in disciplinary actions.

This policy of non-retaliation also applies if reporters are not absolutely certain about the facts, provided that they act in good faith. Acting in good faith means that one has reasonable grounds to suspect that, in light of the circumstances available at the time of reporting, the information reported is true.

However, intentionally misusing the reporting system and this Policy for abusive practice (i.e., filing reports in bad faith) is strictly prohibited. An Employee who knowingly reports incorrect information is not protected by this Policy and may face disciplinary, criminal and civil sanctions.

Further, this Policy does not guarantee protection from disciplinary action if the reporter was involved in the reported breach. Nevertheless, the fact of having reported said breach might be considered as a mitigating factor in any disciplinary proceedings.

3.2. Confidentiality and Anonymity

Confidentiality of the reporting as well as the accused persons and any third parties mentioned in their reports will be preserved as much as possible.

Sartorius' reporting system allows anonymous reports to extent this is permitted by applicable laws and regulations, in particular by data protection and labor laws and regulations. Despite this, it may be possible that the role as a reporting person or the accused person or any third party might become apparent to others during later investigations. In any case, Sartorius will make every effort to keep the identity of the reporter and the reports confidential unless:

- The reporter has given consent for disclosure;
- Disclosure is necessary for compliance with a legal obligation to which Sartorius is subject;
- Disclosure is required for Sartorius to effectively investigate the report and take related measures;
- Disclosure to law enforcement authorities upon their request.

Reporters should keep in mind that investigating a report and therefore addressing the raised concerns is usually facilitated if one does not submit the report anonymously and Sartorius can contact the reporter in case of any questions.

3.3. Data Protection

Sartorius attaches great importance to data protection and privacy as well as data security requirements. Sartorius therefore adheres to applicable data protection laws and regulations as a minimum standard also in relation to this Policy.

(a) Informing Individuals Subject to Speak-Up Reports

Sartorius will inform the accused persons whose behavior has been reported after receiving the respective complaint in compliance with applicable data protection laws and regulations.

Accused persons will generally be informed of the alleged facts and of the person or unit in charge of handling the report within one (1) month.

However, when such information may jeopardize the effectiveness of the investigation, the protection of evidence or the reporting process, it must only be provided if and when those risks no longer exist. This requires a case-by-case analysis.

In any case, the identity of the reporting person will not be disclosed unless this is strictly necessary for compliance with a legal obligation to which Sartorius is subject or for other legitimate purposes and furthermore permitted by applicable data protection and labor laws and regulations.

(b) Retention of Records

Records of reports and the related investigations will be stored in accordance with applicable laws as long as this is necessary for the achievement of Sartorius' legitimate interests. Legitimate interests include the establishment, exercise or defense of legal claims as well as the operation of internal reporting channels and an effective compliance management in line with legal obligations to which Sartorius is subject.

Records of your report(s) will, as a rule, be kept for three (3) years after completion of the reporting procedure under this Policy. Records may be kept longer to meet legal obligations (including the German Act on the Protection of Persons reporting Misconduct – Hinweisgeberschutzgesetz ("HinSchG")) as long as it is necessary and proportionate and then be deleted in accordance with applicable data protection laws and regulations.

In line with this, records that relate to pending disputes or investigations are not deleted as long as data storage is necessary for achieving the purposes of such proceedings. In such cases, the information will be stored for the duration of the respective dispute or investigation or even longer, if required and permitted under applicable data protection laws and regulations.

4. Procedure for Submitting Reports and Follow-Up

4.1. Internal Reporting Channels and Submitting Reports

Employees are encouraged to submit reports of suspected irregularities, concerns, doubts and suspicions, at any time.

When submitting a report, Sartorius asks reporters to submit a description of their report. Reporters should note that the more information they provide, the better Sartorius can understand and investigate the information reported. Therefore, reporters should submit all relevant information on the facts in question, including on the people involved and relevant background information as well as supporting documents of other evidence that could be helpful in investigating the report.

When submitting information, reporters are free to choose whether they like to disclosure their identity or remain anonymous, provided that this is permitted by applicable laws.

Sartorius has implemented various complaint channels that are available at all times and in many languages.

(a) Whistleblower Portal

Sartorius has established the following online whistleblower portal ([Compliance | Sartorius](#)). The portal allows for written reports in various different languages. The portal is available at all times and allows to complain anonymously.

(b) Telephone Hotline

Reports can also be made by telephone. There is a toll-free number (00800 22 44 22 11) and a chargeable number (0049 551 50 41 99 41) where reporters can make a complaint orally.

(c) Email Inbox

Reporters can also contact the Compliance Team in Legal Affairs & Compliance directly by sending an e-mail to compliance@sartorius.com, speaking with them in person or by telephone describing your concern.

(d) Surface Post

Reports can also be sent by conventional surface post to the following address:

Sartorius AG
Attention: Chief Compliance Officer
Otto-Brenner-Str. 20
37079 Göttingen

(e) In Person

Reports can also be in person to the Chief Compliance Officer.

Of course, reporters are free to make a report also to their direct supervisor or a member of the line management, who will then pass the report on to the channels listed above.

(f) Country-specific Channels

In certain countries additional local reporting channels for receipt and handling of reports are available. The reporting channels and further information can be found on the Sartorius website and below in **Annex 1**. Sartorius' central reporting channels are only considered as supplementary channels for reports on the respective local entities.

4.2. External Reporting Channels

Sartorius encourages everyone to submit reports directly and internally at Sartorius as such compliance violations can jeopardize our core business values and high standards of ethics and integrity.

However, reporting persons are permitted to report certain violations to the competent national authorities. Reporting persons may refer to the competent national authority's website for information on when and how to report according to applicable laws and regulations. Employees can find further information on the respective external channels for the respective country under **Annex 2**.

5. Handling of Complaints

Complaints are handled by Corporate Compliance. Corporate Compliance is

- Impartial,
- Independent, i.e. not bound by instructions,
- Committed to confidentiality,
- Specially trained, and
- Equipped with the necessary resources to handle the complaint appropriately.

Corporate Compliance will handle the complaint independently, objectively and impartially.

After receipt of a report, Corporate Compliance will follow-up with the reporter and keep the reporter informed about the status of the review of the report in accordance with applicable laws.

6. Questions and concerns

If you have any questions about this Policy, please contact your Regional Compliance Officer or Corporate Compliance. They will provide you with all the support and information needed. The same applies if you identify possible violations of this Policy.

7. Revision History

Version	Effective Date	Description of Revision
1.0	13.12.2023	New Policy
[...]	[...]	[...]

Annex 1: Internal Local Reporting Channels (country specific)

Country	Internal Local Reporting Channels
<p>Belgium:</p> <p>Xpress Biologics SA</p>	<p>Reports can be made via the following channel(s)</p> <p>Whistleblower Portal Compliance Sartorius</p> <p>Telephone Hotline 0032 487 24 70 89</p> <p>Email Inbox compliance@xpress-biologics.com</p>
<p>Slovenia:</p> <p>Sartorius BIA Separations, separacijske tehnologije, d.o.o.</p>	<p>Reports can be made via the following channel(s)</p> <p>Person of Trust Kim Matavž Stefanovski</p> <p>Whistleblower Portal Compliance Sartorius</p> <p>Telephone Hotline 00386 59 699 506</p> <p>Email Inbox zaupnik@biaseparations.com</p>

Sweden:

Sartorius Stedim Data Analytics AB

Reports can be made via the following channel(s)

Whistleblower Portal

[Compliance | Sartorius](#)

Telephone Hotline

0046 90 184850

0046 73 682 4850

0046 76 145 7378

Email Inbox

WhistleblowSSDA@Sartorius.com

Annex 2: External Reporting Channels

As laid out above in clause 4.2 Sartorius encourages everyone to submit reports directly and internally at Sartorius as such compliance violations can jeopardize our core business values and high standards of ethics and integrity.

However, reporting persons are permitted to report certain violations to the competent national authorities. Reporting persons may refer to the competent national authority's website for information on when and how to report according to applicable laws and regulations. In addition, external reporting can be made to the institutions, bodies, offices or agencies of the European Union outlined in the table below.

If you choose to report to an external reporting channel you can find further information for your respective country of residence and the European Union as follows:

Country/ EU	External Reporting Channels
European Union	<ul style="list-style-type: none"> ▪ European Commission: Whistleblower (europa.eu) ▪ European Anti-Fraud Office (OLAF): Fraud Reporting Home (europa.eu) ▪ European Maritime Safety Agency (EMSA): Home - EMSA - European Maritime Safety Agency (europa.eu) ▪ European Union Aviation Safety Agency (EASA): EASA European Union Aviation Safety Agency (europa.eu) ▪ European Securities and Markets Authority (ESMA): European Securities and Markets Authority (europa.eu) ▪ European Medicines Agency (EMA): European Medicines Agency (europa.eu)
Austria	<ul style="list-style-type: none"> ▪ Federal Office for the Prevention of and Fight against Corruption: https://www.bak.gv.at/601/ ▪ Auditor Oversight Authority: APAB : Whistleblower-Service ▪ Balance Sheet Accounting Authority: WhistleB, Whistleblowing Centre ▪ Federal Competition Authority: WhistleB, Whistleblowing Centre ▪ Financial Market Authority: Einführung (bkms-system.net) ▪ Money Laundering Reporting Office: Geldwäschemeldestelle (usp.gv.at) ▪ Austrian Chamber of Notaries: WhistleB, Whistleblowing Centre ▪ Austrian Bar Associations: Die Rechtsanwaltskammern - oerak.at ▪ Chamber of tax advisors and auditors: WhistleB, Whistleblowing Centre ▪ Head of the federal disciplinary authority: Einführung (bkms-system.net)
Belgium	<ul style="list-style-type: none"> ▪ Service Federal Public Economy, SMEs, Self-Employed and Energy: Homepage FOD Economie (fgov.be) ▪ Federal Public Service Finance: FOD Financiën (belgium.be)

	<ul style="list-style-type: none"> ▪ Federal Public Service Public Health, Food Chain Safety and Environment: Whistleblowing procedure FPS Public Health (belgium.be) ▪ Federal Public Service Mobility and Transport: mobiliteit.belgium.be/nl ▪ Federal Public Service Employment, Labor and Social Dialogue: Homepage Federale Overheidsdienst Werkgelegenheid, Arbeid en Sociaal Overleg (belgie.be) ▪ Public Programming Service Integration Social, Fight against Poverty, Social Economy and Political of Large Cities: POD Maatschappelijke Integratie POD Maatschappelijke Integratie, Armoedebestrijding, Sociale Economie en Grootstedenbeleid (mi-is.be) ▪ Agency Federal Nuclear Control: FANC - Federaal Agentschap voor Nucleaire Controle Samen beschermen (fgov.be) ▪ Federal Agency for Medicines and Health Products: FAGG Uw geneesmiddelen en gezondheidsproducten (fagg-afmps.be) ▪ Federal Agency for the Safety of the Food Chain: FAVV : Federaal Agentschap voor de Veiligheid van de Voedselketen (favv-afsa.be) ▪ Belgian Competition Authority: Report a competition issue? Belgian Competition Authority ▪ Data Protection Authority: Startpagina burger Gegevensbeschermingsautoriteit ▪ Financial Services and Markets Authority: News & Warnings FSMA ▪ National Bank of Belgium: Welkom bij de Nationale Bank van België nbb.be ▪ Belgian Audit Oversight Board: College van toezicht op de bedrijfsrevisoren (CTR) CTR (ctr-csr.be) ▪ authorities referred to in section 85 of the Act of 18 September 2017 on the prevention of money laundering and financing terrorism and limiting the use of cash: Online melding (ctif-cfi.be) ▪ National Committee for the supply and distribution of drinking water ▪ Belgian Institute for Postal Services and telecommunications: Consumenten BIPT ▪ National Institute for Sickness and Invalidity Insurance: Onthaal - RIZIV (fgov.be) ▪ National Social Insurance Institute for the Self-Employed: RSVZ Sociale Zekerheid Zelfstandige Ondernemers ▪ National Office Employment: Welcome The National Employment Office (rva.be) ▪ National Social Security Office: Home RSZ ▪ Information Service and Social Research: Homepage Sociale Inlichtingen-en Opsporingsdienst (belgie.be) ▪ Autonomous Anti-Fraud Coordination Service (CAF) ▪ Navigation Control: mobiliteit.belgium.be/nl/scheepvaart
Finland	<ul style="list-style-type: none"> ▪ The Office of the Chancellor of Justice: https://oikeuskansleri.fi/en/whistleblower-protection <p>Please note before reporting:</p>

	<p>In Finland, reports to the competent national authorities can be made if the reporting person cannot use the internal reporting channel, the reporting person has a legitimate reason to believe that their internal report has not resulted in measures within the time prescribed or that it is ineffective, or the reporting person has a legitimate reason to believe that they may face retaliation due to their report. Reports to the competent national authority can be made in writing through an electronic form (only in Finnish and Swedish), post or email, or verbally.</p>
France	<ul style="list-style-type: none"> ▪ Criminal authorities ▪ Defender of rights: The Defender of Rights Défenseur des Droits (defenseurdesdroits.fr) ▪ Specific administrative authorities depending on the area concerned (see complete list of the administrative authorities is annexed to the French Decree No. 2022-1284 of October 3, 2022).
Germany	<ul style="list-style-type: none"> ▪ External Federal Reporting Office as part of the Federal Ministry of Justice: Startseite - Externe Meldestelle (Bundesamt für Justiz Schlichtungsverfahren) ▪ German Federal Cartel Office in the event of violations of the law against restraints of competition: Bundeskartellamt - Hinweise auf Kartellverstöße - Hinweise auf Kartelle, Marktmissbrauch und sonstige Verstöße ▪ Federal Financial Supervisory Authority in the event of violations of the financial services supervision act: https://www.bafin.de/DE/DieBaFin/Hinweisgeberstelle/hinweisgeberstelle_node.html
Hungary	<ul style="list-style-type: none"> ▪ Directorate-General for Audit of European Funds: https://eutaf.hu/ ▪ Integrity Authority: https://integritashatosag.hu/ ▪ Hungarian Competition Authority: https://www.gvh.hu/ ▪ Public Procurement Authority: https://www.kozbeszerzes.hu/ ▪ Hungarian Energy and Public Utility Regulatory Authority: https://www.mekh.hu/ ▪ Hungarian National Bank: https://www.mnb.hu/ ▪ Hungarian National Authority for Data Protection and Freedom of Information: https://www.naih.hu/ ▪ National Media and Infocommunications Authority: https://nmhh.hu/ ▪ Hungarian Atomic Energy Authority: https://www.haea.gov.hu/ ▪ Supervisory Authority for Regulated Services: https://sztfh.hu/
Ireland	<ul style="list-style-type: none"> ▪ The Office of the Protected Disclosures Commissioner: OPDC Home Protected Disclosures Commissioner OPDC.ie
Italy	<ul style="list-style-type: none"> ▪ National Anti-Corruption Authority: Autorità Nazionale Anticorruzione <p>Please note before reporting:</p>

	<p>In Italy, reporting via external reporting channels is only permitted if one of the following conditions applies:</p> <ul style="list-style-type: none"> ▪ The reporting channel has not been activated by the company or does not meet the requirements set by the Italian Transposing Law; ▪ The reporting person has already submitted a report which the Company has failed to follow up on; ▪ The reporting person has a legitimate reason to believe that the Company would not follow up effectively on his/her internal report, or that the company would apply retaliatory measure against him/her as a result of his/her internal report or ▪ The reporting person has a legitimate reason to believe that the relevant violation may cause imminent or obvious harm to the public interest.
Netherlands	<ul style="list-style-type: none"> ▪ Authority for Consumers & Markets (<i>Autoriteit Consument en Markt</i> (ACM)): https://www.acm.nl/nl/bescherming-klokkenluiders ▪ Dutch Data Protection Authority (<i>Autoriteit persoonsgegevens</i> (AP)): https://autoriteitpersoonsgegevens.nl/nl/zelf-doen/gebruik-uw-privacyrechten/klacht-melden-bij-de-ap. ▪ The Dutch Whistleblowers Authority (<i>Huis voor klokkenluiders</i> (HvK)): https://www.huisvoorklokkenluiders.nl/onderzoek-door-het-huis ▪ Health and Youth Care Inspectorate (<i>Inspectie gezondheidszorg en jeugd</i> (IGJ)): https://www.igi.nl/onderwerpen/melden-over-uw-werkgever-klokkenluidersregeling ▪ Dutch Healthcare Authority (Nederlandse Zorgautoriteit (NZa)): https://www.nza.nl/contact/ons-meldpunt/meldpunt-misstanden-bij-zorgverzekeraars-of-zorgaanbieders
Slovenia	<ul style="list-style-type: none"> ▪ The Slovenian Commission for the Prevention of Corruption (CPC): Commission for the Prevention of Corruption GOV.SI ▪ The Slovenian Securities Market Agency (ATVP): https://www.a-tvp.si/eng. ▪ The Slovenian Competition Protection Agency (AVK): Slovenian Competition Protection Agency: Home (varstvo-konkurence.si) ▪ The Slovenian Insurance Supervision Agency (AZN): Home (a-zn.si) ▪ Slovenian Traffic Safety Agency: STSA - Slovenian Traffic Safety Agency - Agencija za varnost prometa - AVP (avp-rs.si) ▪ Agency for Public Oversight of Auditing: Zunanja prijava kršitve predpisa v delovnem okolju po ZZPri - ANR ▪ Bank of Slovenia: Reporting (bsi.si) ▪ National Commission of review for the Review of Public Procurement Procedures: National Review Commission for Reviewing Public Procurement Procedures GOV.SI ▪ Financial Administration of the Republic of Slovenia: FINANCIAL ADMINISTRATION OF THE REPUBLIC OF SLOVENIA (gov.si)

	<ul style="list-style-type: none"> ▪ Market Inspectorate of the Republic of Slovenia: Market Inspectorate GOV.SI ▪ Office for the Prevention of Money Laundering: Office for Money Laundering Prevention GOV.SI ▪ Information Commissioner: Information Commissioner GOV.SI ▪ Information Security Inspection: Government Information Security Office GOV.SI ▪ Radiation and Nuclear Safety Inspection: Radiation and Nuclear Safety Inspection GOV.SI ▪ Radiation Protection Inspection: Slovenian Radiation Protection Administration ▪ Inspection for Food Safety, Veterinary and Plant Protection: Inspection for Food Safety, Veterinary Sector and Plant Protection GOV.SI ▪ Labour Inspectorate of the Republic of Slovenia: Labour Inspectorate GOV.SI ▪ Public Sector Inspectorate: About the Public Sector Inspectorate GOV.SI ▪ Inspectorate of the Republic of Slovenia for the Environment and Spatial Planning: Natural Resources and Spatial Planning Inspectorate GOV.SI ▪ Slovenian Agency for Medicinal Products and Medical Devices: JAZMP - Agency for Medicinal Products and Medical Devices of the Republic of Slovenia ▪ Health Inspectorate of the Republic of Slovenia: Health Inspectorate GOV.SI ▪ Slovenian Sovereign Holding: SDH - Home <p>Please note before reporting:</p> <p>In Slovenia, reporting via external reporting channels is only permitted if one of the following conditions applies:</p> <ul style="list-style-type: none"> ▪ An internal reporting channel has not been established; ▪ A report could not be handled internally in an effective way; ▪ The reporting person believes there is the risk of retaliation in case they report via the internal channel.
Sweden	<ul style="list-style-type: none"> ▪ Swedish Work Environment Authority: https://www.av.se/om-oss/visselblasarlagen/extern-rapporteringskanal/# ▪ National Board of Housing, Building and Planning: https://www.boverket.se/sv/om-boverket/visselblasning/ ▪ National Electrical Safety Board: https://www.elsakerhetsverket.se/yrkespersoner/tillverka-och-salja-elprodukter/sla-larm-om-missforhallanden/ ▪ Swedish Economic Crime Authority: https://www.ekobrottsmyndigheten.se/visselblsarfunktion-eu-medel/ ▪ Swedish Estate Agents Inspectorate: https://fmi.se/det-har-ar-fmi/kontakta-oss/visselblasning-om-penningtvatt-eller-finansiering-av-terrorism/ ▪ Swedish Financial Supervisory Authority:

<https://www.finansinspektionen.se/sv/om-fi/kontakta-oss/visselblasare/>

- Public Health Agency of Sweden:
<https://www.folkhalsomyndigheten.se/livsvillkor-levnadsvanor/andts/regler-for-tillverkning-handel-och-hantering/visselblasning-tobaksområdet/>
- Swedish Agency for Marine and Water Management:
<https://www.havochvatten.se/om-oss-kontakt-och-karriar/om-oss/visselblasarfunktion.html>
- Swedish Authority for Privacy Protection:
<https://www.imy.se/privatperson/utfora-arenden/visselblasning/>
- Inspectorate of Strategic Products:
<https://isp.se/om-isp/visselblasning-till-isp/hur-rapportering-ska-ske/>
- Health and Social Care Inspectorate:
<https://www.ivo.se/om-ivo/kontakta-oss/visselblasning/extern-rapporteringskanal/>
- Swedish Chemicals Agency: <https://www.kemi.se/om-kemikalieinspektionen/kontakta-oss/extern-kanal-for-visselblasning>
- Swedish Consumer Agency:
<https://www.konsumentverket.se/om-konsumentverket/var-verksamhet/visselblasning/extern-kanal-for-visselblasning/>
- Swedish Competition Authority:
<https://www.konkurrensverket.se/tipsa-oss/visselblasarfunktion/>
- Swedish Food Agency:
<https://www.livsmedelsverket.se/om-oss/kontakt/visselblasning--rapportera-om-missforhallanden>
- Swedish Medical Products Agency:
<https://www.lakemedelsverket.se/sv/om-lakemedelsverket/kontakta-oss/visselblasning#hmainbody4>
- County Administrative Boards:
www.lansstyrelsen.se
- Swedish Civil Contingencies Agencies:
<https://www.msb.se/sv/om-msb/kontakta-oss/visselblasning--rapportera--om-missforhallanden/visselblasning--rapportera--om-missforhallanden-inom-området-produktsakerhet-och-produktoverensstammelse/>
- Swedish Environmental Protection Agency:
<https://www.naturvardsverket.se/om-oss/kontakt/visselblasning/rapportera-missforhallanden-inom-vissa-tillsynsomraden/>
- Swedish Post and Telecom Authority:
<https://www.pts.se/sv/bransch/radio/Radioutrustning/visselblas-spegel-3/>
- The Government Offices:
<https://www.regeringen.se/om-webbplatsen/rapportera-missforhallanden-om-statsstod/>

- Swedish Inspectorate of Auditors:
<https://www.revisorsinspektionen.se/tillsyn/rapportering-om-missforhallanden/>
- Swedish Tax Agency:
<https://www.skatteverket.se/omoss/varverksamhet/styrningochuppfoljning/skattekontroller/rapporteraommissforhallandeninomskatteomradet.4.1df9c71e181083ce6f636e5.html?q=visselbl%C3%A5sning>
- Swedish Forest Agency:
<https://www.skogsstyrelsen.se/kontakt/visselblasarfunktion/>
- Swedish Gambling Authority:
<https://www.spelinspektionen.se/lagar--forordningar/penningtvatt/visselblasarfunktion/>
- Swedish Energy Agency:
<https://report.whistleb.com/sv/Energimyndigheten>
- Swedish Board of Agriculture:
<https://jordbruksverket.se/e-tjanster-databaser-och-appar/ovriga-e-tjanster-och-databaser/anmal-offentliga-kontroller-visselblasarfunktion>
- Swedish Board for Accreditation and Conformity Assessment:
<https://www.swedac.se/visselblasning/>
- Swedish Radiation Safety Authority:
<https://www.stralsakerhetsmyndigheten.se/kontakt/visselblasarfunktion/>
- Swedish Transport Agency:
<https://www.transportstyrelsen.se/sv/Om-transportstyrelsen/visselblasning/>

For Sweden please note:

An external reporting channel is a channel established by an authority for handling reporting on breaches that are within the authority's field of responsibility. As with internal reporting, the protection of the Swedish Whistleblower Act applies if the reporting is made to the relevant competent authority via the external reporting channel.

In addition to the Swedish Whistleblower Act please consider the following: The Freedom of the Press Act (SFS 1949:105) and the Fundamental Law on Freedom of Expression (SFS 1991:1469) provide for the right to submit information on any subject for publication in certain media and the right to acquire information on any subject for the purpose of communicating it for publication in certain media.