

## Privacy Policy

**Effective Date: June 9th 2017**

This privacy statement describes how SignalChem and its affiliates (“SignalChem Lifesciences and SignalChem Pharmaceuticals Inc.”) collects and uses the personal information you provide on our website: [www.signalchem.com](http://www.signalchem.com). It also describes the choices available to you regarding our use of your personal information and how you can access and update this information.

### Collection and Use of Personal Information.

We collect the following personal information from you:

- Contact Information such as name, email address, mailing address, phone number
- Billing Information such as credit card number and billing address
- Unique Identifiers such as user name, account number, password

We also collect the following information from you:

- Information about your business such as company name and business type
- Demographic information such as education, interests and postal/zip code

As is true of most websites, we automatically gather information about your computer such as your IP address, browser type, referring/exit pages, operating system and store it in log files.

We may combine this automatically collected log information with other information we collect about you. We do this to improve services we offer you.

We use this information to:

- Send general communications
- Send you newsletters
- Send you promotional communications
- Process and track transactions
- Provide products and services to you
- Recommend products and series that SIGNALCHEM believes will be of interest and provide value to you

### Choice/Opt-Out

You may choose to stop receiving our newsletter or promotional communication by following the unsubscribe instructions included in these emails or you can contact us at [info@signalchem.com](mailto:info@signalchem.com).



#110-13120 Vanier Place  
Richmond, BC V6V2J2  
CANADA  
1-866-9 KINASE  
Tel: 1-(604)-232-4600  
Fax: 1-(604)-232-4601  
[www.signalchem.com](http://www.signalchem.com)

### **Refer-A-Friend**

If you choose to tell a friend about our site, we will ask you for your friend's name and email address. We will automatically send your friend a one-time email inviting him or her to visit the site. SignalChem stores this information for the sole purpose of sending this one-time email and tracking the success of our referral program.

Your friend may contact us at [info@signalchem.com](mailto:info@signalchem.com) to request that we remove this information from our database.

### **Information Sharing**

We will share your personal information including your email address with third parties only in the ways that are described in this privacy statement. We do not sell your personal information to third parties. We may provide your personal information including your email address to companies that provide services to help us with our business activities such as shipping your order or offering customer service. These companies are authorized to use your personal information only as necessary to provide these services to us.

In certain situations, SignalChem may be required to disclose personal data in response to lawful requests by public authorities, including to meet national security or law enforcement requirements.

Also, we may also disclose your personal information as required by law such as to comply with a subpoena, or similar legal process, when we believe in good faith that disclosure is necessary to protect our rights, protect your safety or the safety of others, investigate fraud, or respond to a government request, if SignalChem is involved in a merger, acquisition, or sale of all or a portion of its assets, you will be notified via email and/or a prominent notice on our website of any change in ownership or uses of your personal information, as well as any choices you may have regarding your personal information, to any other third party with your prior consent to do so.

### **Tracking Technologies**

Technologies such as: cookies, beacons, tags and scripts are used by SignalChem and our online support providers. These technologies are used in analyzing trends, administering the site, tracking user's movements around the site and to gather demographic information about our user base as a whole. We may receive reports based on the use of these technologies by these companies on an individual as well as aggregated basis.

We use cookies for our shopping cart, to remember users' settings (e.g. language preference) and for authentication. Users can control the use of cookies at the individual browser level. If you reject cookies, you may still use our site, but your ability to use some features or areas of our site may be limited.

### **Links to Other Web Sites**

Our website includes links to other websites whose privacy practices may differ from those of SignalChem. If you submit personal information to any of those sites, your information is governed by their privacy statements. We encourage you to carefully read the privacy statement of any website you visit.

### **Social Media Buttons and Widgets**

Our website includes Social Media Features, such as the Facebook, LinkedIn and Twitter buttons and widgets, such interactive mini-programs that run on our site. These features may collect your IP address, which page you are visiting on our site, and may set a cookie to enable the feature to function properly. Social Media Features and widgets are either hosted by a third party or hosted directly on our site. Your interactions with these features are governed by the privacy policy of the company providing it.

### **Surveys or Contests**

From time-to-time we may provide you the opportunity to participate in surveys on our site. If you participate, we will request certain personal information from you. Participation in these surveys is completely voluntary and you therefore have a choice whether or not to disclose this information. The requested information typically includes contact information (such as name and shipping address), and demographic information (such as zip code).

We use this information to notify contest winners and award prizes, to send participants an email, or to collect customer testimonials.

### **Security**

The security of your personal information is important to us. We follow generally accepted industry standards to protect the personal information submitted to us, both during transmission and once we receive it. No method of transmission over the Internet, or method of electronic storage, is 100% secure, however. Therefore, we cannot guarantee its absolute security. If you have any questions about security on our website, you can contact us at [info@signalchem.com](mailto:info@signalchem.com). When you enter sensitive information (such as a credit card number) on our order forms, we encrypt the transmission of that information using secure socket layer technology (SSL).

### **PCI Compliance**

We are PCI compliant because nothing is more important than keeping our customer's payment data secure. The PCI Data Security Standard represents a common set of industry tools and measurements to help ensure the safe handling of sensitive information. The standard provides an actionable framework for developing a robust account data security process - including preventing, detecting and reacting to security incidents.



#110-13120 Vanier Place  
Richmond, BC V6V2J2  
CANADA  
1-866-9 KINASE  
Tel: 1-(604)-232-4600  
Fax: 1-(604)-232-4601  
[www.signalchem.com](http://www.signalchem.com)

### **Customer Testimonials**

We post customer testimonials on our website which may contain personal information. We do obtain the customer's consent prior to posting the testimonial to post their name along with their testimonial. To request removal of your personal information from testimonials please contact us at [info@signalchem.com](mailto:info@signalchem.com).

### **Updating/Accessing/Correcting Personally Identifiable Information**

SignalChem will provide you with information about whether we hold any of your personal information. If you're personally identifiable information or your email address changes, or if you no longer desire our service, you may correct, update, amend, delete/remove or deactivate it by making the change on our member information page or by emailing our Customer Support at [info@signalchem.com](mailto:info@signalchem.com) or by contacting us by telephone or postal mail at the contact information listed below. We will respond to your request to access within 30 days.

### **Data Retention**

We will retain your information for as long as your account is active or as needed to provide you services. If you wish to cancel your account or request that we no longer use your information to provide you services contact us at [info@signalchem.com](mailto:info@signalchem.com). We will retain and use your information as necessary to comply with our legal obligations, resolve disputes, and enforce our agreements.

### **Notification of Privacy Policy Changes**

We may update this privacy policy to reflect changes to our information practices. If we make any change in how we use your personal information including your email address we will notify you by email (sent to the e-mail address specified in your account) or by means of a notice on this site prior to the change becoming effective. We encourage you to periodically review this page for the latest information on our privacy practices.

SignalChem  
110-13120 Vanier Place Richmond,  
BC, Canada, V6V 2J2.  
[info@signalchem.com](mailto:info@signalchem.com)  
Tel: 1.604.232.4600.  
Toll Free Tel: 1.866.9.KINASE