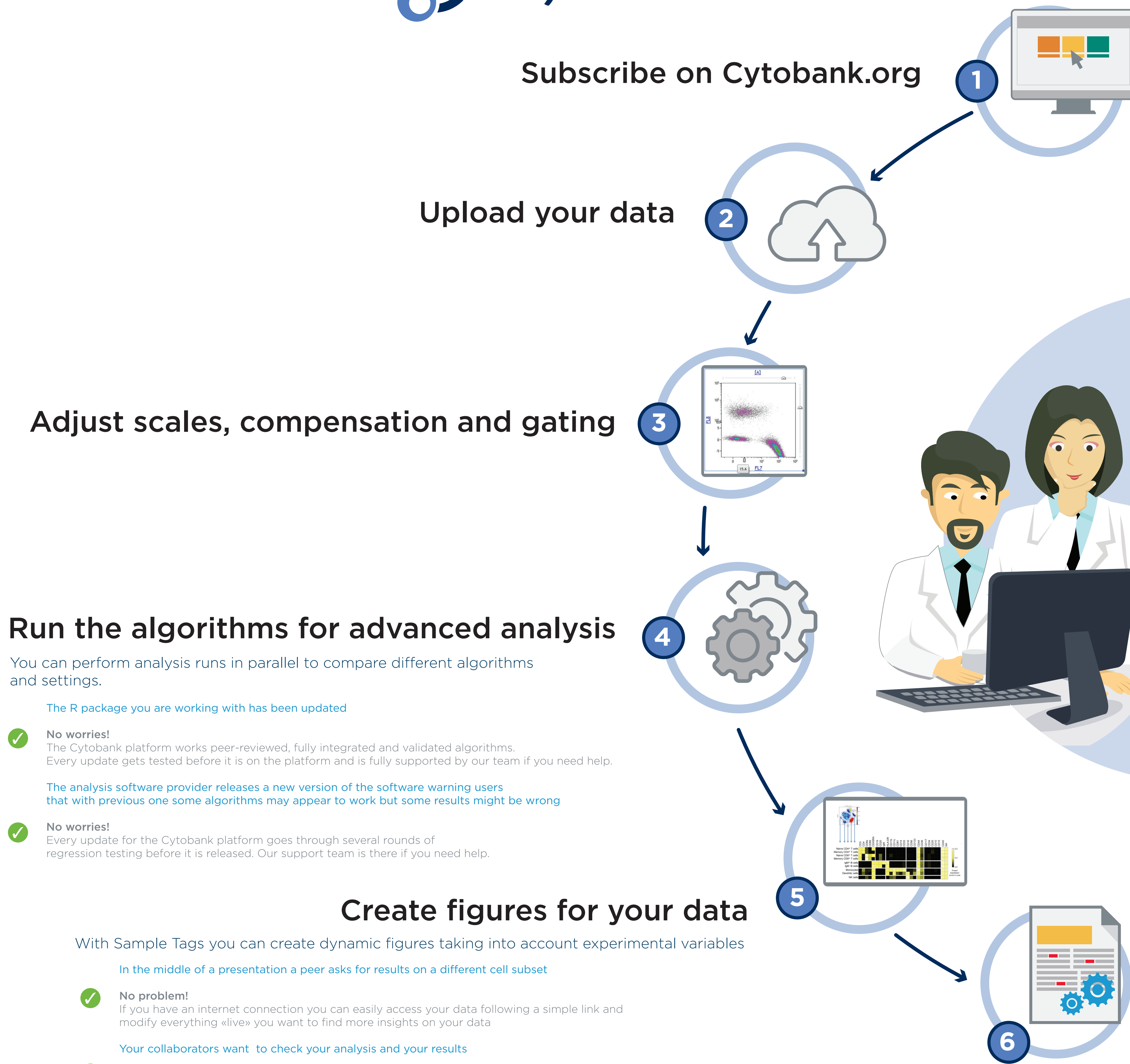


I'm a Biologist, not a Programmer!



Other Solutions



Subscribe on Cytobank.org

Upload your data

Adjust scales, compensation and gating

Run the algorithms for advanced analysis

You can perform analysis runs in parallel to compare different algorithms and settings.

The R package you are working with has been updated

✓ No worries!
The Cytobank platform works peer-reviewed, fully integrated and validated algorithms. Every update gets tested before it is on the platform and is fully supported by our team if you need help.

The analysis software provider releases a new version of the software warning users that with previous one some algorithms may appear to work but some results might be wrong

✓ No worries!
Every update for the Cytobank platform goes through several rounds of regression testing before it is released. Our support team is there if you need help.

Create figures for your data

With Sample Tags you can create dynamic figures taking into account experimental variables

In the middle of a presentation a peer asks for results on a different cell subset

✓ No problem!
If you have an internet connection you can easily access your data following a simple link and modify everything «live» you want to find more insights on your data

Your collaborators want to check your analysis and your results

✓ No worries!
• Create a project, click on «share» on the Experiment Manager and decide what level of access to give your collaborators
• You can collaborate with several users on the same analysis file.
• You get your reviewed analysis and your insights

You can submit your paper

Reviewer #2 asks for the results of a run with different settings

✓ No worries!
In the Cytobank platform everything is archived and settings remain associated with results

The Journal you submitted your paper to requires making your data publicly available

✓ No worries!
In the Cytobank platform you can make your experiment public with just one click



Get your dongle/web-based license

Download the correct plugin for the algorithm

- Add the plugin of interest to the plugins folder on your machine and restart
- Be sure to put the plugin in the correct folder so that it is recognized by your analysis software. If not, specify folder's file path and restart
- If the plugin requires R environment, make sure you have the R package installed and all the algorithm dependencies it requires in R
- If your user accounts does not have Admin rights you might need your IT department help

Adjust scales, compensation and gating

Run the algorithms for advanced analysis

⚠ You chose the wrong settings
Repeat 4

⚠ You chose the wrong algorithm
Repeat 2

The R package you are working with has been updated

✗ There was a bug in a previous version of the package that was later fixed. Now you don't know if your findings are correct and can't easily re-trace your steps. Repeat 2

The analysis software provider releases a new version warning users that with previous one some algorithms may appear to work but some results might be wrong

✗ Again, you don't know if your findings are correct or not. Repeat 2

Create figures for your data

You need to check lab book for notes on experimental variables every time you modify figures

In the middle of a presentation a peer asks for results on a different cell subset

✗ You have an internet-based license for your analysis so you can use it everywhere but you realize you should have copied all your data and analysis and took them with you to be able to change something!

Your collaborators want to check your analysis and your results

✗

- You need to find an empty hard disk to copy your data on
- Data and analysis are big: go and get a coffee while you wait
- Data and analyses are really big: go and get a second coffee while you wait
- Send data to your collaborators
- They change analysis to version _V1
- You make the requested changes and change analysis to version _V2
- They forgot to make a small change and change analysis to version _V2.1
- You get your insights from version _V2.1.2.3

You can submit your paper

Reviewer #2 asks for the results of a run with different settings

✗ Where did I put that analysis?????

The Journal you submitted your paper to requires making your data publicly available

✗ You need to create a new account on a repository and upload all your files