

TIME MATTERS: VISUALIZE YOUR RESULTS, FASTER

New, faster algorithms for dimensionality reduction

Join the Cytobank platform and access the computational power of the cloud; use faster implementations of dimensionality reduction algorithms with no need for powerful desktop hardware or programming skills.

U UMAP Widely believed to preserve global structure better than other solutions. Join the team and analyze up to 10M events using GPU-UMAP.

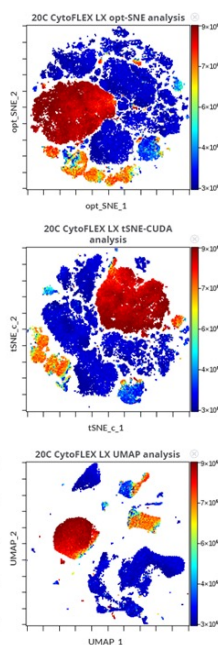
tSNE-CUDA Run up to 10M events in less than an hour* and get the SNE maps you're familiar with. Capture even subtle differences within cellular subsets.

opt-SNE Enable optimal data resolution in t-SNE space and precise data interpretation by automating the selection of settings. Start your run with up to 3M events.

* Time might vary depending on dataset

Accelerate the time to results with the "Compare" function

Want to explore all the available option on your data or compare results of different algorithms? Easy, just run UMAP, tSNE-CUDA or opt-SNE on the same data set. Visualize and study results directly on the Analysis Setup page.



The Cytobank platform is a cloud-based single-cell analysis solution for high-dimensional data, visual data management and controlled sharing.



Get Personalized Support

Our Application Specialists at Cytobank_Support@beckman.com
Your local sales representative at Cytobank_Sales@beckman.com



Access online training resources

Support Articles at support.cytobank.org
Videos and more in the Cytobank Learning Center on beckman.com



Access Machine Learning-assisted analysis with no need to code



Upload any type of data in a tabular format



Securely Store & Easily Share Data



Organize Data & Create Auditable Workflows