



Dear Valued Customer,

Molecular Devices is writing to inform you of a significant development regarding its partner, Core Life Analytics (CLA). Due to ongoing financial challenges and a sharp decline in revenue caused by the current economic climate, CLA has been unable to secure new funding. As a result, CLA no longer has the resources or team capacity to continue their operations.

This difficult decision means that access to StratoMineR cannot be guaranteed beyond **Friday, Feb 7th, 2025** and work on collaborative projects will unfortunately cease. Please rest assured that CLA will maintain the security of all data in their systems until it can be safely and permanently deleted.

If you have critical data stored with CLA that is not already backed up locally, we strongly recommend downloading it before the above date. Should you require assistance with this process, please reach out to CLA directly at support@corelifeanalytics.com, and they will provide the necessary support. Please reach out to me if you have any questions regarding cease of operations of CLA.

Please direct your questions to any of the following methods:

Molecular Devices North America Customer Support: 1.800.635.5577 or email support@moldev.com

Core Life Analytics Support: Please email support@corelifeanalytics.com

As always, we remain committed to supporting your research and providing you with the highest quality of products and services. Thank you for the opportunity to continue to support and help advance your research and discovery.

Please let us know if we can answer any questions.

Thank you!