



The sky's the limit with Tecan's Te-Cloud™

Take your Te-Care Service Contract to new heights with Te-Cloud

Tecan is going above and beyond to bring you our newest service offering, Te-Cloud, a state of the art Freedom EVOware® back-up and computer protection service.

Add Te-Cloud to your Te-Care Complete Service Contract today to give your team peace of mind with secure and simple file recovery.

Te-Cloud service includes daily automatic on-line backups of all your Freedom EVOware configuration files with three months of saved history. This backup schedule is fully managed by Tecan with file encryption for security and confidentiality. Computer replacement is also included in the event of a hardware failure.

With Te-Cloud you will:

- Save time recreating your scripts and custom files with lightning fast restore
- Receive a replacement computer in the event of a hardware failure
- Ensure maximum safety with automatic back-ups and file encryption
- Free up space on your internal servers with cloud-based storage
- Avoid unexpected expenses when failures occur



Take advantage of free installation by upgrading your current Te-Care Complete service contract. Te-Cloud services are available as a service contract upgrade or a stand-alone offering.

Services are currently available for our US and Canada customers. For more details contact us at InsideSalesUS@tecan.com today!

www.tecan.com

Call: +1 919 361 5200 Email: InsideSalesUS@tecan.com