

DATA ONLINE

A powerful web-based, secure system to view your study information in real-time.

■ DOWNLOAD YOUR DATA AND REPORTS DIRECTLY

Cerep's Data Online service can be accessed using Internet Explorer 5+, Netscape Navigator 6+ and Mozilla.

To experience the power and convenience of Data Online, go to <https://www.cerep.com/secure>, and use "Demo" as login and password.

■ TWO OPTIONS

- The realtime data option enables the sponsor to view study data as they are produced, and also to convert them into MSExcel.
- The final report option enables the sponsor to download complete study reports and other related documents as they are generated.

■ SECURITY

- To deliver the world's highest level of trust, a company that specializes in providing High-Assurance SSL Certificates authenticates our organization, enabling end users to verify our site and communicate via state-of-the-art SSL encryption.
- Our server supports all browser SSL encryption types (40 bit, 56 bit, and 128 bit strong encryption).
- Passwords are not stored in any database; they are owned and encrypted by the operating system security mechanism. Only the sponsor has knowledge of the password.
- You can safely enter your password and download files from the secure server.
- Our network topology to access the Data Online web server is also secured via state-of-the-art network security solutions.
- Server operating systems and software are updated immediately upon receipt of the provider's security bulletin.

To experience the power and convenience of Data Online, please go to www.cerep.com **DATA ONLINE**

■ FAQs

Here are answers to the most frequently asked questions about Data Online.

For further information, please contact customer.support@cerep.fr.

- **In the Final Report section, I click on MSWord or MSExcel document, Data OnLine asks me for a username and password. What's wrong?**
This is a known bug (see Microsoft reference) within version 2000 of MSOffice. Ask your IT team to update your version to Service Release 2 (SR-2).
- **I made a shortcut to Data Online on my computer but when I use it I am redirected to Data Online homepage.**
For security reasons you cannot access Data Online without entering your personal and confidential information.
- **When I click on the logout button and then access Data Online again, I am not asked for my login/password.**
This is a bug in Microsoft Internet Explorer version 4. When a secure site is accessed and the user logs out without shutting down the browser and then accesses the site a second time, the system will not ask for identification. To avoid this, logout from Data Online and close your browser. You can also update your system to Internet Explorer 5+.
- **Cerep has informed me by e-mail that my final study report is available but I cannot see any new document in the Final report section?**
The service that is indexing the documents is working asynchronously so when your document is published to Data Online it might take a little time to appear. To see all your documents, switch to explorer view.

Ask your Cerep representative: sales@cerep.com
customer.support@cerep.fr for access information and technical assistance.



FRANCE

Le Bois l'Evêque
86600 CELLE L'EVESCAULT
tel. +33 (0)5 49 89 30 00

(Headquarters)

155 boulevard Haussmann
75008 PARIS
tel. +33 (0)1 45 64 44 60

USA

15318 N.E. 95th Street
REDMOND, WA 98052
tel. +1 (425) 895 8666

JAPAN

Namiki Shoji Co., Ltd.
Kenseishinjuku Bldg. 5-5-3
Shinjuku, Shinjuku-ku
TOKYO, 160-0022
tel. +81 (0)3 3354 4026
fax +81 (0)3 3352 2196

CHINA

Ai Di Sheng (Edison) Road 326,
302-1 room
Zhangjiang High-Tech Park
SHANGHAI
tel. +86 18702160370