

Transmit Your Screen Contents View Online – Worldwide

... With Security

Factsheet
Security & Technology

With the use of pcvisit, you and your customers can utilize unique communication advantages.

To ensure your data to be secure on their way through the Internet, we have equipped pcvisit with many security mechanisms.

The communication via pcvisit is not only comfortable, but is secure too.

Individual visitor number provides undisturbed communication

The pcvisit session starts after the guest and the host have telephonically communicated an eight-digit connection number created by a random number generator. Not before the correct connection number has been input, the connection between guest and host is established. The individual connection number for each pcvisit session ensures that the two call partners will have a communication completely undisturbed by thirds.

128-bit encryption prevents bugging operations

Symmetrical encryption methods (128-bit Blowfish algorithms) as well as a secure data transmission via multilevel security layers offer a "bugging-proof line". Moreover, during the entire connection time, pcvisit actively prevents the data from being manipulated by thirds. Thus, the authenticity of both sides is fully guaranteed. With this high security standard, a pcvisit visit becomes so secure as online banking does.

View direction change with permission only

At the beginning of the pcvisit session, the guest initially sees her/his host's screen. If the view direction shall be reversed, the reversal can be performed not before the guest has agreed by a mouse click. In this way – with acknowledgment – there can be switched over as often as desired. Moreover, at the same time, there is ensured that either can only see the information destined for her/him.

Remote control rights with permission only

Each pcvisit session starts with showing the host's screen contents. Neither the host nor the guest have access rights to the call partner's computer.

By a mouse click, the remote control right is released. Not before the person opposite has acknowledged, the remote PC can be controlled remotely. When this is done, each activity can be traced via the moved mouse pointer. Here, invisible manipulations are being excluded.

The ongoing remote control is shown via an activated button. By a mouse click, the remote access or even the entire pcvisit session can be aborted immediately.

Comfortable technology in the background ... with security

With http, pcvisit utilizes the conventional Internet protocol that is accepted by any firewall and any proxy server. Thus, in 99% of all cases, a secure data connection can also be established behind firewalls of sealed-off company networks.

A pcvisit visit is completely browser-independent. To set up a pcvisit session, a Windows PC with Internet connection is sufficient. This provides additional security reserves in the cooperation with various call partners and computer configurations. After the pcvisit visit, the guest's computer is in its original state. On the guest computer, no changes such as registry entries are made. No installation is performed. In this way, pcvisit combines simple utilization and data security.

Verisign certificate confirms program code authenticity

The internationally accepted certification authority Verisign has given pcvisit a code signing certificate. Thus, there is ensured that each pcvisit module downloaded in the Internet actually represents the original pcvisit software. On each download, there is checked whether the program code corresponds to the original software and is transmitted without any changes.

A manipulation of the transmitted software by thirds is excluded. With the code signing certificate, you thus get an additional security from a neutral authority whenever a pcvisit session takes place.