



Short instruction pcvisit "Guest - Module"

Please enter the number given to you by your session partner (host module) in the field „session no.“ .



If you have no dedicated line to the internet, make a manual connection to the internet. Enter the number and then press the receiver symbol.



If the connection is successful, the pcvisit window looks like this:



Now you see the desktop of your session partner.

Enjoy the presentation!