

PIFOC E-662 RS-232 Serial Cable

Abstract

This technical note describes the connectors and pinouts for use with the Physik Instrumente PIFOC E-662 controller, so that the device can be communicated with by computer over the RS-232 Serial Port

Instructions

The Physik Instrumente PIFOC E-662 Controller can be controlled by either analog voltage or serial communications. This document describes the cable connectors and pinouts for the serial communications cable.

Each connector on the cable must be a DB-9 female connector. The pinout diagram is as follows:

PIFOC	Computer
1	1
2	3
3	2
4	6
5	5
6	4
7	8
8	7
9	9

Document ID
T20027

Product
MetaMorph
MetaFluor
MetaVue

Platform
Windows 95/98
Windows NT
Windows 2000

Created
25-Jun-2004

Last Reviewed
25-Jun-2004

Keywords: pifoc e-662 e662 rs232 rs-232 serial cable
Issue Type: Hardware