

## Uniblitz Shutter Serial Cable

### Abstract

The Uniblitz Shutter can be controlled through a serial port.

### Instructions

#### Cable Information

The cable must have a DB-9 Female to connect to the computer serial port and a DB-9 Male to connect to the Uniblitz serial port. Wire the cable as follows:

| Computer side                                     | Uniblitz side                                     |
|---|---|
| 2   | 3   |
| 3   | 2   |
| 5   | 5   |
| 4 - 6 (short these two together in the connector) | 4 - 6 (short these two together in the connector) |
| 7 - 8 (short these two together in the connector) | 7 - 8 (short these two together in the connector) |

**Document ID**  
T10163

**Product**  
MetaMorph  
MetaFluor  
MetaGFP

**Created**  
08-Oct-2000

### Serial Communications Protocol

No Parity, 8 data bits, 1 stop bit. The BAUD rate can be found by looking in the Vincent Associates Uniblitz manual for your shutter. Old shutters run at 300 baud, while new ones run at 9600 baud.

### Driver

For Meta Imaging Series version 3.5 and beyond, you can use the UniblitzS driver to control the D122 or T122 shutter. For Meta Imaging Series version 4.5.4 and beyond, you can use the UniblitzD3s driver to control the VMM-D3 shutter.

For information on Parallel Port cables, see [T10030](#).

**Keywords:** uniblitz shutter serial cable  
**Issue Type:** Hardware