

## Hamamatsu Orca Cable Information

### Abstract

Definitions of the cables used by Hamamatsu Orca series cameras

### Instructions

Hamamatsu Orca-series cameras (including the Orca 100, Orca ER, Orca 2, C7300-51, C4742-95, C4742-98, and C4880 series) using the RS-422 (non-firewire) interface, require two cables to connect to the computer.

The first cable is the serial cable. This has a DB-9 Female connector on each end. The serial cable is straight-through, meaning pin 1 connects to pin 1, 2 to 2, 3 to 3, and so on, for all 9 pins.

The second cable is the data cable. The data cable has a connector of the same appearance on each end. However, the cable is direction-specific. The end of the cable with RED TAPE must connect to the Hamamatsu camera controller. The end of the cable with BLUE TAPE must connect to the MV-1500 board in the computer.

**Document ID**  
T20003

**Created**  
10-Jan-2002

**Last Reviewed**  
1/10/2002

**Keywords:** hamamatsu, orca, mv-1500, cables  
**Issue Type:** Hardware