

## Discovery-1 Software Parallel Port Pin Numbering

### Abstract

Getting the Discovery-1 software numbers parallel port pins to match the numbers on the connector.

### Instructions

Getting the Discovery-1 software numbers parallel port pins to match the numbers on the connector.

The following pins are inputs, and can carry a signal from the outside to your imaging software: 10, 11, 12, 13, and 15.

The following pins are outputs, and can carry a signal from your imaging software to external devices: 2, 3, 4, 5, 6, 7, 8, 9, 14 and 16. Pins 14 and 16 are commonly reserved for shutters; use them only if you need more than eight outputs and will attach no shutters to the associated parallel port.

Parallel Port Pins 1 and 17 should never be used.

Parallel Port Pins 18-25 are grounds.

**Document ID**  
D10124

**Product**  
Discovery-1 software

**Created**  
24-Jun-1998

**Last Reviewed**  
14-Jul-2003

**Keywords:** discovery-1 software parallel port pin numbering  
**Issue Type:** geninfo