

## EFOS/EXFO X-Cite

### Abstract

The EFOS/EXFO X-Cite is an illumination source which can be shuttered by computer control.

### Instructions

The EFOS X-Cite is an illumination source that consists of a metal-halide lamp, liquid light guide, and mechanical shutter. Metamorph controls the shutter as a standard shutter through an available parallel port. The shutter is not fast.

In order to control the device, a custom cable must be configured.

Parallel Port (25pin Dsub Male)	Aux. Connector on X- Cite Controller (9pin Dsub Male)
Pin 15	Pin 2*
Pin 18	Pin 3*, Pin4
Pin 3	Pin 5
Pin 4	Pin 7

Note: Pin3 and Pin4 on the X-Cite controller must be connected together.

\* Note: You must also connect a 1K-Ohm resistor from Pin2 to Pin 3 on the X-Cite controller.

**Document ID**  
T20008

**Created**  
13-Jun-2001

**Last Reviewed**  
20-Apr-2002

**Keywords:** efos exfo x-cite shutter illumination  
**Issue Type:** Hardware