

---

## Linkam Temperature Controlled XY Stage

### Abstract

MetaMorph controls the Linkam MDS 600 motorized stage, using the Linkam TMS93 temperature controller. This is a heating and cooling stage.

### Instructions

#### Manufacturer Information

Information on the MDS 600 stage can be obtained by clicking on <http://www.linkam.co.uk/mds.html>.

Information on the TMS93 temperature controller can be obtained by clicking on <http://www.linkam.co.uk/programmers.html>.

**Document ID**  
T10254

**Product**  
MetaMorph

**Created**  
08-Jul-2001

**Keywords:** metamorph, linkam, hardware  
**Issue Type:** Hardware