

Nikon Remote Focus Accessory - Configuration Information

Abstract

This document contains the DIP switch configuration information for the Nikon Remote Focus Accessory.

Instructions

Configuration Information

The Nikon Remote Focus Accessory has a 4-position DIP switch visible on the rear of the unit. The switches are labelled S1, S2, S3, S4. Each switch can be either Open or Closed. The default MetaMorph configuration is as follows:

S1: Open

S2: Closed

S3: Closed

S4: Open

This configures the device for 1200 baud, 8 data bits, 1 stop bit, no parity, and for use as Unit #0.

Document ID
T10222

Product
MetaMorph
MetaVue

Created
13-Aug-1999

Keywords: nikon remote focus accessory

Issue Type: Hardware