

The Format of IMG Files

Abstract

The Format of IMG Files.

Instructions

MetaMorph-family software, Image-1/AT and Image-1/FL and can each load and store images in IMG files. MetaMorph calls these UIC format images. Image-1/AT refers to these as Image-1 format images. All Image-1/FL image files have this format. IMG files come in two varieties. One stores eight-bit images, and the other stores 16-bit images. Eight-bit IMG files have a 6-byte header, organized as three 16-bit unsigned integers. The first integer has the value 16. The second is the image width. The third is the image height. The image data follows the header. The image data and the integers in the header are stored less significant byte-first.

Document ID

T10128

Product

MetaMorph

MetaFluor

Image-1

Created

17-Jul-1998

Keywords: metamorph, metafluor, image-1, image-1/fl, image files

Issue Type: Application