

Using 12 bit images with Application Modules

Abstract:

This document describes how to use 12 bit images with application modules, which require 16 bit images.

Instructions:

MetaMorph® software works with and stores 5 types of images: binary (0, 1), 8 bit (0-255), 16 bit (0 - 65535), 24 bit (0-255, 0-255, 0-255), and 48 bit (0-65535, 0-65535, 0-65535).

When a camera is called 12 bit it is a 12 bit dynamic range (grey values 0 – 4095). Images from a 10, 12, or 14 bit cameras are stored within MetaMorph® software as 16 bit images that you can use with any of the application modules.

ARTICLE #T20130

PRODUCTS
MetaFluor®
MetaMorph®

CREATED
8-July-2009

LAST UPDATED
31-July-2009