

Using Visual Basic with MetaMorph and MetaXpress

OVERVIEW

MetaMorph's extended set of Microsoft Visual Basic (VB) programming functions enable you to create Visual Basic programs that can be run in MetaMorph or MetaXpress using the Run User Program drop-in command. Creating VB programs for use in MetaMorph makes it possible for you to process and analyze images within the application using the power and flexibility of a programming language.

For more information on using Visual Basic with MetaMorph and MetaXpress, refer to the MetaMorph Visual Basic Reference Guide, available on the [Molecular Devices MetaMorph support page](#).

Note: The following versions of Microsoft Visual Basic can be used to create programs to run in MetaMorph 7.0:

- Microsoft Visual Basic .NET 2003 Professional or Enterprise editions
- Microsoft Visual Basic .NET 2002 Professional or Enterprise editions
- Microsoft Visual Basic 6.0 Professional or Enterprise editions
- Microsoft Visual Basic 5.0 Professional or Enterprise editions

NOTES:

- Programs created using the Standard and Express editions of the above versions are NOT supported.
- The version of the *MetaMorph Visual Basic Reference Guide* included on the MetaMorph 7.0 CD erroneously states that Visual Basic .NET is not supported in MetaMorph 7.0.

ARTICLE #
T20051

PRODUCTS
MetaMorph® 7.0 and
above
MetaXpress 1.7 and
above

CREATED
03-March-2006

LAST UPDATED
07-March-2006