

The Power of Precision



Precision Melt Analysis Software is a convenient, easy-to-use application that imports and analyzes data files generated from the CFX Opus 96, CFX96 Touch, CFX96, CFX Opus Deepwell, CFX96 Touch Deep Well, CFX Opus 384, CFX384 Touch, CFX384, CFX Duet, or CFX Connect Real-Time PCR System to genotype samples based on the thermal denaturation properties of double-stranded DNA. The software can be used for a variety of genotyping applications, including scanning for new gene variants, screening DNA samples for single nucleotide polymorphisms (SNPs), identifying insertions/deletions or other unknown mutations, and determining the percentage of methylated DNA in unknown samples.

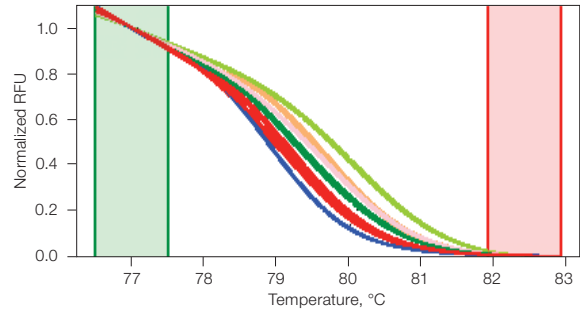
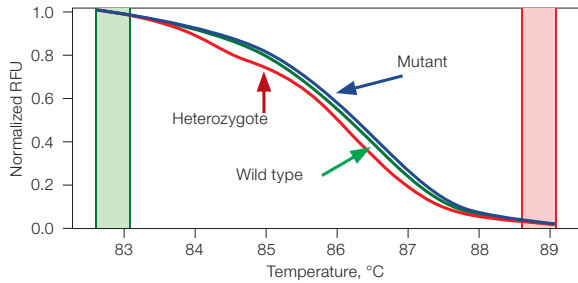
Precision Melt Analysis Software makes it easy for you to:

- Streamline your data analysis using the customizable default analysis settings
- Tailor the software to the appropriate analysis by using multiple data view options to manually assign sample genotypes
- Examine results from a number of melt files, without having to export data, using the Melt Study module
- Analyze multiple experiments from a single plate using the Well Groups feature
- Publish your data in several formats by easily exporting data and images

Efficient and Effective Analysis

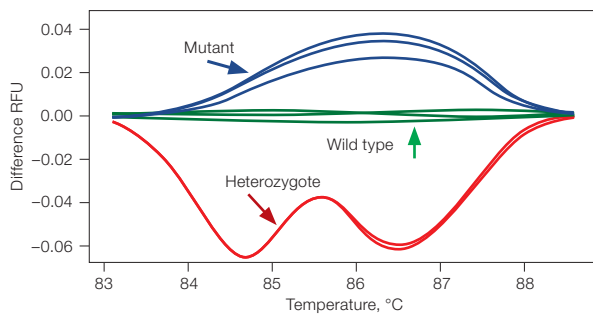
Precision Melt Analysis Software saves analysis time by assigning sample genotypes automatically based on cluster analysis, or allowing manual use of multiple data view options to tailor the software to the appropriate analysis. Use the normalized melt curves plot feature to generate a basic representation of the different clusters based on curve shifting (for homozygotes) and curve shape change (for heterozygotes). Difference curve plots of a sample fluorescence versus a selected control at each temperature transition provide a convenient visual aid to interpret the data.

Precision Melt Analysis Software enables data comparison between multiple file runs by combining data into a single Melt Study. Develop a standard library of melt curve runs to analyze an unlimited number of melt experiments without having to export data.



SNP genotyping by Precision Melt Analysis Software using data generated by the CFX384 Real-Time PCR System. Discrimination of human hemochromatosis S65C SNP genotypes (A to T substitution). Data from homozygous wild-type (■), mutant (■), and heterozygote (■) samples are shown on a normalized melt curve plot. RFU, relative fluorescence units.

Multifile melt study analysis. Compare melt curves from different files using the Melt Study module in Precision Melt Analysis Software. RFU, relative fluorescence units.



A representative difference curve plot generated from data normalized to a mutant sample. Data from homozygous wild-type (■), mutant (■), and heterozygote (■) samples are shown. RFU, relative fluorescence units.

Ordering Information

Catalog #	Description
1845025	Precision Melt Analysis Software , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit
12011319	CFX Opus 96 Real-Time PCR System , includes CFX Opus 96 System and communication cables
12011452	CFX Opus 384 Real-Time PCR System , includes CFX Opus 384 System and communication cables
12016658	CFX Opus Deepwell Real-Time PCR System , includes CFX Opus Deepwell System and communication cables
12016265	CFX Duet Real-Time PCR System , includes CFX Duet System, power cord, USB cable

Visit bio-rad.com/PrecisionMeltAnalysisSoftware for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

