

The Power of Precision

Precision Melt Analysis software is a convenient, easy-to-use application that imports and analyzes data files generated from the CFX96™ or CFX384™ real-time PCR detection 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 SNPs, identifying insertions/deletions or other unknown mutations, and determining the percentage of methylated DNA in unknown samples.

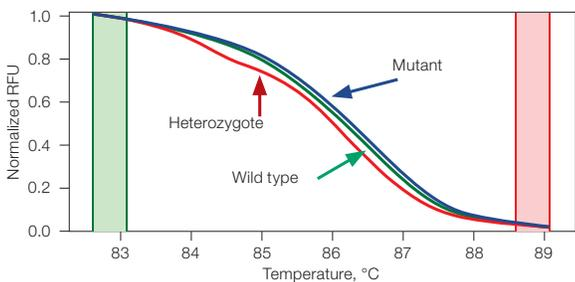
Precision Melt Analysis software makes it easy for you to:

- Streamline your data analysis using the customizable default analysis settings
- Utilize the multiple data view options to manually assign sample genotypes by tailoring the software to the appropriate analysis
- 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

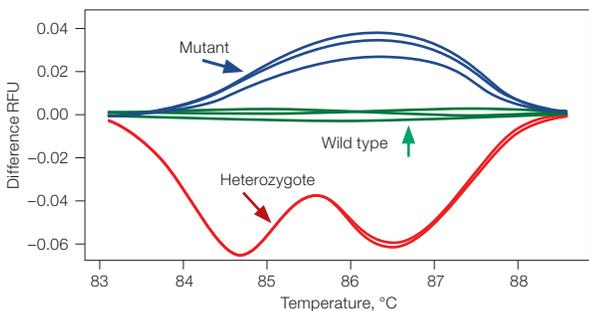
Efficient and Effective Analysis

Precision Melt Analysis software saves analysis time by assigning sample genotypes automatically based on cluster analysis, or manually using 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.

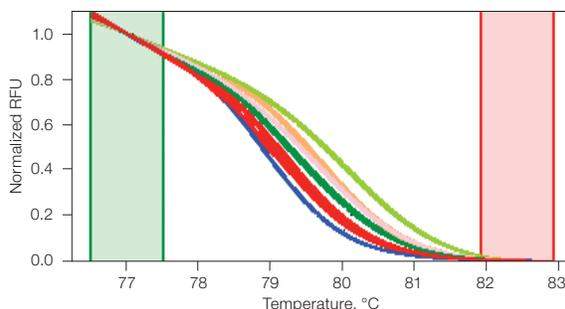
Precision Melt Analysis software enables data comparison between multiple file runs by combining data into a single Melt Study. Develop a standard library of melt curve runs to analyze an unlimited number of melt experiments without having to export data.



SNP genotyping by Precision Melt Analysis software using data generated by the CFX384 real-time PCR detection 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.



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.



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

Ordering Information

Catalog #	Description
184-5025	Precision Melt Analysis Software , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit
185-5096	CFX96 Real-Time PCR Detection System , includes C1000™ thermal cycler chassis, CFX96 optical reaction module, CFX Manager™ software, license for qbase ^{PLUS} software, communication cable, reagents, consumables
185-5384	CFX384 Real-Time PCR Detection System , includes C1000 thermal cycler chassis, CFX384 optical reaction module, CFX Manager software, license for qbase ^{PLUS} software, communication cable, reagents, consumables

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen dye for use in real-time PCR, for research purposes only.

Purchase of SsoFast EvaGreen supermix includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Notice regarding Bio-Rad thermal cyclers and real-time systems:

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under one or more of U.S. Patent Numbers 5,656,493; 5,333,675; 5,475,610 (Claims 1, 44, 158, 160-163, and 167 only); and 6,703,236 (Claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems Corporation. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied Biosystems' United States Patent Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000