

# Trademarks labels

ABI Prism® 5700, 7000, 7500, 7700 and 7900 are registered trademarks of The Perkin-Elmer Corp.

Acrydite® is a registered trademark of Mosaic Technologies Corp.

Alexa Fluor® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

AnaSpec® is a registered trademark of AnaSpec, Inc.

AnaTag™ is a trademark of AnaSpec, Inc.

ArrayIt® is a registered trademark of TeleChem International, Inc.

Benzonase® is a registered trademark of Merck KGaA, Darmstadt, Germany.

BHQ-0™, BHQ-1™, BHQ-2™, BHQ-3™ are registered trademarks of Biosearch Technologies, Inc.

BlackBerry® is a registered trademark of Berry & Associates, Inc.

Black Hole Quencher® is a registered trademark of Biosearch Technologies, Inc.

BODIPY® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

CAL Fluor® is a trademark of Biosearch Technologies, Inc.

Cascade Blue® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

CDP Star® is a registered trademark of Applera Corp.

Cherry™ is a trademark of Delphi Genetics S.A.

Cy® is a registered trademark of GE Healthcare Bio-Sciences Corp.

Cy5® Direct is a registered trademark of GE Healthcare Bio-Sciences Corp.

DHL™ is a trademark of AnaSpec, Inc.

Diamond *Taq*® is a registered trademark of Eurogentec S.A.

DNA Engine Opticon® is a registered trademark of Bio-Rad Laboratories, Inc.

Dragonfly Orange™ is a trademark of Eurogentec S.A.

Eclipse® is a registered trademark of Epoch Biosciences, Inc.

Eppendorf® is a registered trademark of Eppendorf AG.

Excel™ is a trademark of Microsoft Corp.

Eurogentec® is a registered trademark of Eurogentec S.A.

Exiqon® is a registered trademark of Exiqon A/S

FLAG® is a registered trademark of Sigma-Aldrich Co

GeneAmp® is a registered trademark of Roche Molecular Systems, Inc.

GenePix® is a registered trademark of Astros Acquisition Sub II.

GenePix 4000® is a registered trademark of Astros Acquisition Sub II.

GetStaby™ is a trademark of Delphi Genetics S.A. Gold'N Oligo® is a registered trademark of Eurogentec S.A.

GO® Peptide is a trademark of AnaSpec, Inc.

GoldStar® is a registered trademark of Eurogentec S.A.

HiLyte™ is a trademark of AnaSpec, Inc.

iCycler iQ® is a registered trademark of Bio-Rad Laboratories, Inc.

iCycler® is a registered trademark of Bio-Rad Laboratories, Inc.

I-MUPID® is a registered trademark of Advance Co., Ltd.

iQ™5 is a trademark of Bio-Rad Laboratories, Inc.

iQ-CHECK® is a registered trademark of Bio-Rad Laboratories, Inc.

LightCycler® is a registered trademark of Roche Diagnostics GmbH.

LNA® is a registered trademark of Exiqon A/S.

Lumavidin® is a registered trademark of Luminex Corp.

Luminex® is a registered trademark of Luminex Corp.

Luminex® 200™ is a trademark of Luminex Corp.

Luminex® XYP™ is a trademark of Luminex Corp.

Luminex® SD™ is a trademark of Luminex Corp.

Luminex® SeroMAP™ is a trademark of Luminex Corp.

MagPlex™ is a trademark of Luminex Corp.

MapMarker® is a registered trademark of BioVentures, Inc.

Marina Blue® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

MasterCycler® is a registered trademark of Eppendorf, Inc

MGB® is a registered trademark of Epoch Biosciences, Inc.

MGB Eclipse® is a registered trademark of Epoch Biosciences, Inc.

MicroPlex™ is a trademark of Luminex Corp.

Mupid® is a registered trademark of Advance Co., Ltd.

Mx3000® is a registered trademark of Stratagene

Mx3000P® is a registered trademark of Stratagene

Mx3005P® is a registered trademark of Stratagene

Mx4000® is a registered trademark of Stratagene

NASBA® is a registered trademark of bioMerieux

Oli&GO 1•2•3™ is a trademark of Eurogentec S.A.

OliGold® is a registered trademark of Eurogentec S.A.

Onco-Mir® is a registered trademark of Ambion Inc.

Opticon® is a registered trademark of MJ Research, Inc.

Oregon Green® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Oyster® is a registered trademark of Denovo Biolabels

Pacific Blue™ is a trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Plexor™ is a trademark of Promega Corporation

Pulsar™ is a trademark of Biosearch Technologies, Inc.

Quantica® is a registered trademark of Techne Ltd and Techne, Inc.

Quasar™ is a trademark of Biosearch Technologies, Inc.

Qiagen® and QiaQuick® are trademarks of Qiagen, Inc.

QXL™ is a trademark of AnaSpec, Inc.

Rhodamine Red™ is a trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Rotor-Gene® is a registered trademark of Corbett Research Pty Ltd

ROX™ is a trademark of Applera Corp.

RP-Cartridge•Gold™ is a trademark of Eurogentec S.A.

SELEX™ is a trademark of Gilead Sciences, Inc.

SensoLyte® is a registered trademark of AnaSpec, Inc.

Smartcycler® is a registered trademark of Cepheid, Inc.

StabyCloning™ is a trademark of Delphi Genetics S.A.

StabyExpress™ is a trademark of Delphi Genetics S.A.

Staby™ Codon is a trademark of Delphi Genetics S.A.

StepOne® is a registered trademark of Applera Corp. or its subsidiaries

SuperScript® is a registered trademark of Invitrogen Corp.

SYBR® Green is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

Taqman® is a registered trademark of Applied Biosystems, Inc.

Texas Red® is a registered trademark of Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corp.)

TRIZol® is a registered trademark of Molecular Research Center Inc.

UltraPureGold™ is a trademark of Eurogentec S.A.

VersArray® is a registered trademark of Bio-Rad Laboratories, Inc.

Vic® is a registered trademark of Applera Corp.

Word™ is a trademark of Microsoft Corp.

xPONENT® is a registered trademark of Luminex Corp.

xMAP® is a registered trademark of Luminex Corp.

Yakima Yellow® is a registered trademark of Epoch Biosciences, Inc.

Z-Fish® is a registered trademark of AnaSpec, Inc.