

**applied biosystems**

by Thermo Fisher Scientific

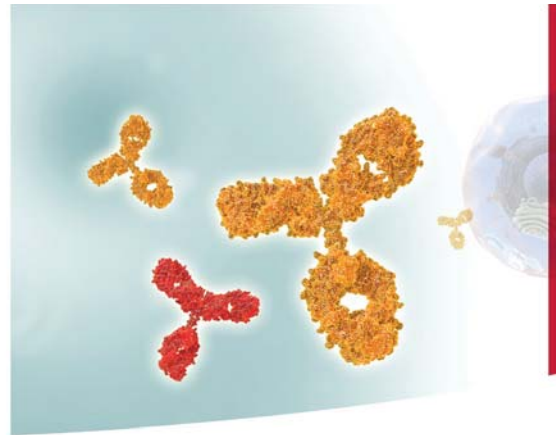
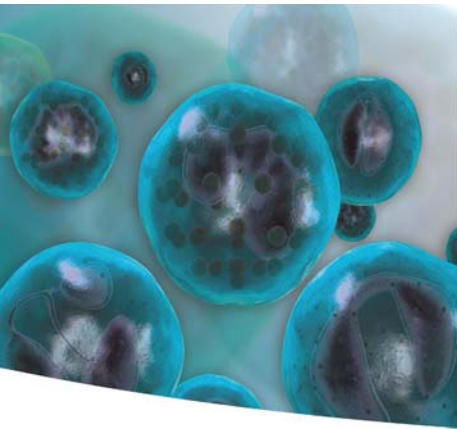
Now a part of our Microarray solutions portfolio

Visit us at [thermofisher.com/microarray](http://thermofisher.com/microarray)

# PharmacoScan™ Assay 24-Array Format

## Manual Protocol

## SitePreparationGuide



eBioscience

GeneChip

USB

## Trademarks

Affymetrix®, Axiom®, GeneChip®, Command Console®, CytoScan®, DMET™, Eureka™, Eureka Genomics®, Eureka Genotyping™, Expression Console™, GeneAtlas®, GeneChip-compatible™, GeneTitan®, Genotyping Console™, myDesign™, MyGeneChip™, NetAffx®, OncoScan®, PharmacoScan™, Powered by Affymetrix™, PrimeView®, and ViewRNA® are trademarks or registered trademarks of Affymetrix, Inc. Please see [www.affymetrix.com/trademarks](http://www.affymetrix.com/trademarks) for a complete list of Affymetrix trademarks. All other trademarks are the property of their respective owners.

## Limited License

Affymetrix hereby grants to buyer a non-exclusive, non-transferable, non-sublicensable license to Affymetrix' Core Product IP to use the product(s), but only in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix. "Core Product IP" is the intellectual property owned or controlled by Affymetrix as of the shipment date of a product that covers one or more features of the product that are applicable in all applications of the product that are in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix. The license granted herein to buyer to the Core Product IP expressly excludes any use that: (i) is not in accordance with the product labels, inserts, manuals and written instructions provided by Affymetrix, (ii) requires a license to intellectual property that covers one or more features of a product that are only applicable within particular fields of use or specific applications, (iii) involves reverse engineering, disassembly, or unauthorized analysis of the product and/or its methods of use, or (iv) involves the re-use of a consumable product. Buyer understands and agrees that except as expressly set forth, no right or license to any patent or other intellectual property owned or controlled by Affymetrix is granted upon purchase of any product, whether by implication, estoppel or otherwise. In particular, no right or license is conveyed or implied to use any product provided hereunder in combination with a product or service not provided, licensed or specifically recommended by Affymetrix for such use. Furthermore, buyer understands and agrees that buyer is solely responsible for determining whether buyer possesses all intellectual property rights that may be necessary for buyer's specific use of the product, including any rights from third parties.

## Copyright

© 2016 Affymetrix Inc. All rights reserved.

# Contents

---

<b>Chapter 1</b>	<b>Equipment and Supplies for the PharmacoScan™ Assay 24-Array Format Manual Protocol</b> . . . . .	<b>4</b>
	Affymetrix Equipment, Software, Reagents and Arrays Required . . . . .	4
	Related Documentation . . . . .	5
	Equipment Required for Manual Target Prep . . . . .	5
	Pre-Amplification Area/Amplification Staging Area . . . . .	6
	Oven Requirements . . . . .	6
	Spectrophotometer . . . . .	6
	Thermal Cycler Recommendations and Protocols . . . . .	7
	Thermal Cycler Consumables . . . . .	8
	Plate Centrifuge . . . . .	8
	Sample Plates . . . . .	9
	Pipettes and Tips . . . . .	9
	Other Labware Required . . . . .	10
	Other Reagents and Gels Required . . . . .	10
	Other Lab Equipment Required . . . . .	11

---

<b>Appendix A</b>	<b>Supplier Contact List.</b> . . . . .	<b>12</b>
-------------------	---	-----------



**Table 1.1** Affymetrix Equipment, Software, Reagents and Arrays Required (Continued)

✓	Item	Part Number
<b>PharmacoScan™ Kits</b>		
☐	PharmacoScan™ Training 24F Kit for Manual Protocol Each Kit Includes: <ul style="list-style-type: none"> <li>■ PharmacoScan™ Reagent Kit 4x24 Reactions</li> <li>■ Four PharmacoScan™ 24F Array Plates</li> <li>■ Four Axiom® GeneTitan® Consumables Kits</li> <li>■ Two PharmacoScan™ DNA Training Plates, for 24 format Assay</li> </ul>	903011
☐	PharmacoScan™ Assay 24F Kit for Manual Protocol Each Kit Includes: <ul style="list-style-type: none"> <li>■ PharmacoScan™ Reagent Kit 4x24 Reactions</li> <li>■ Four PharmacoScan™ 24F Array Plates</li> <li>■ Four Axiom® GeneTitan® Consumables Kits</li> </ul>	903010

## Related Documentation

- *PharmacoScan™ Assay 24-Array Format Manual Protocol User Guide*, P/N 703286
- *PharmacoScan™ Assay 24-Array Format Manual Protocol QRC*, P/N 703288
- *Axiom™ Analysis Suite User Guide*, P/N 703307
- *Axiom® Genotyping Solution Data Analysis Guide*, P/N 702961
- *GeneTitan® MC Protocol for Axiom 2.0 Array Plate Processing QRC*, P/N 702988
- *Affymetrix® GeneChip® Command Console® 4.0 User Guide*, P/N 702569
- *Affymetrix® Genotyping Console™ 4.2 User Guide*, P/N 702982
- *GeneTitan® Multi-Channel Instrument User Guide*, P/N 08-0308
- *GeneTitan® Multi-Channel Instrument Site Preparation Guide*, P/N 08-0305

## Equipment Required for Manual Target Prep

**Table 1.2** Additional Instruments Required for the PharmacoScan™ Assay 24-Array Format Manual Protocol

✓	Equipment	Manufacturer/ Distributor	Part Number
☐	Two ovens (see <a href="#">Oven Requirements on page 6</a> )		
☐	Fume Hood (Strongly recommended: see the <i>PharmacoScan™ Assay 24-Array Format Manual Protocol User Guide</i> for more information)		
☐	Thermal Cycler: <ul style="list-style-type: none"> <li>■ ABI 9700</li> <li>■ ABI Veriti™</li> <li>■ ABI ProFlex™</li> <li>■ Eppendorf Mastercycler™ pro S</li> </ul> (see <a href="#">Thermal Cycler Recommendations and Protocols on page 7</a> )	ABI ABI ABI Eppendorf	
☐	One of the following shakers: <ul style="list-style-type: none"> <li>■ Compact Digital Microplate Shaker</li> <li>■ Jitterbug™</li> </ul>	Thermo Scientific Boekel Scientific	88880023 Model 130 000

## Pre-Amplification Area/Amplification Staging Area

Precautions are required when manipulating genomic DNA or setting up amplification reactions to avoid contamination with foreign DNA amplified in other reactions and procedures. It is recommended that genomic DNA manipulations and amplification reaction set up are performed in a dedicated rooms or areas separate from the main laboratory.

These areas should have dedicated sets of pipettes and plasticware. If no dedicated area is available, use of a dedicated bench or a dedicated biosafety hood and dedicated pipettes is suggested. If no dedicated bench or biosafety hood is available, a set of dedicated pipettes is recommended.

Ideally, the pre-amplification and amplification staging areas would be separate; however these areas may be combined due to space and equipment limitations.

## Oven Requirements

Two ovens are required for PharmacoScan manual target preparation.

**Table 1.3** Ovens Required for the PharmacoScan Assay 24-Array Format Manual Protocol

✓	Equipment	Manufacturer	Part Number
☐	Two ovens, any combination of the following types:		
	<ul style="list-style-type: none"> <li>■ ED 56 Drying and Heating Chamber*</li> </ul>	BINDER	ED056UL-120V Voltage: 120 V 1~60 Hz P/N 9010-0334
			ED056-230V Voltage: 230 V 1~50/60 Hz P/N 9010-0333
	<ul style="list-style-type: none"> <li>■ Affymetrix® GeneChip® Hyb Oven 645†</li> </ul>	Affymetrix	00-0331

\* Replaces BINDER Model ED 53.

† The GeneChip® Hybridization Oven 640 is currently not supported with the PharmacoScan Assay 24-Array Format Manual Protocol.

## Spectrophotometer

We recommend using one of the spectrophotometers listed in [Table 1.4](#).

**Table 1.4** Spectrophotometers

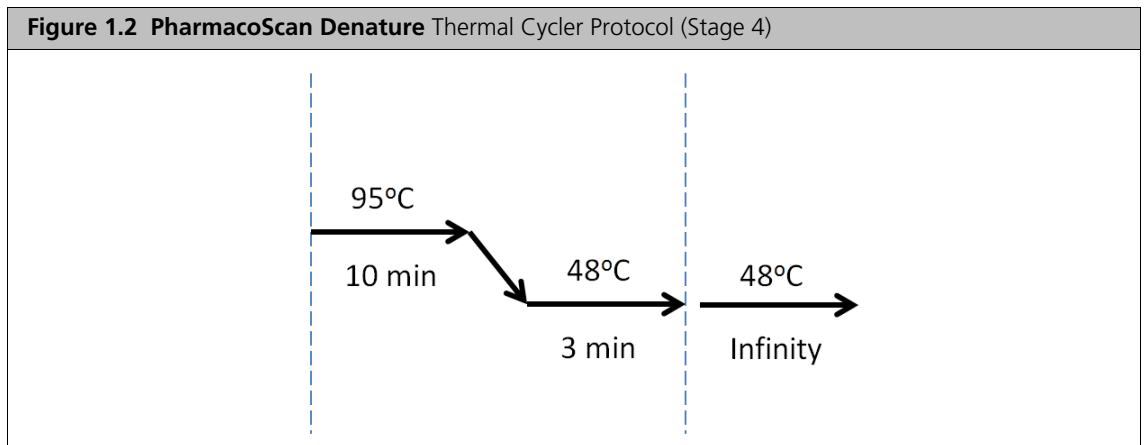
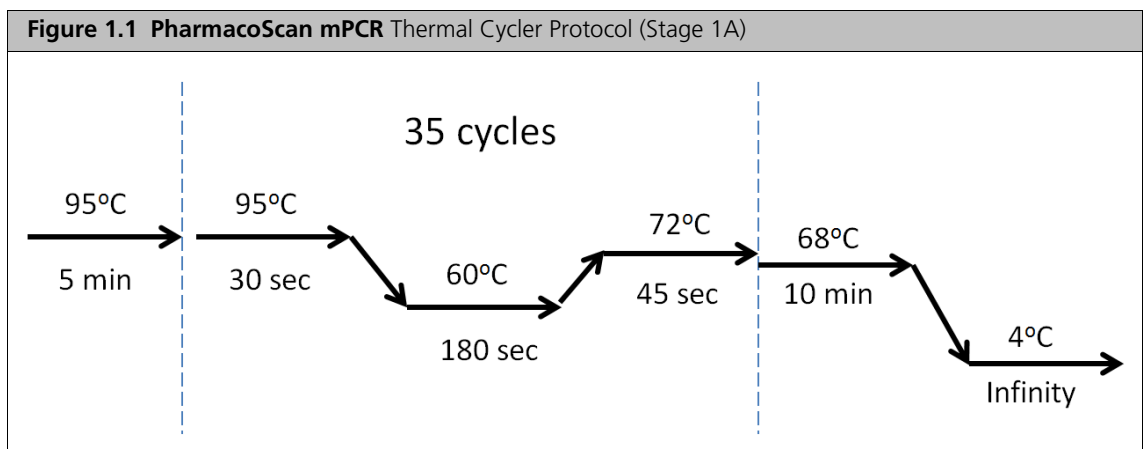
✓	Item	Supplier	Part Number
☐	One of the following spectrophotometers:		
	<ul style="list-style-type: none"> <li>■ DTX 880 Multimode Detector, with:               <ul style="list-style-type: none"> <li>☐ Genomic Filter Slide</li> </ul> </li> </ul>	Beckman Coulter	987921 – detector
	or	Beckman Coulter	A30184 – filter slide
	<ul style="list-style-type: none"> <li>■ SpectraMax High throughput Microplate Spectrophotometer</li> </ul>	Molecular Devices	Plus384

## Thermal Cycler Recommendations and Protocols

The following thermal cycler has been validated for the PharmacoScan Assay 24-Array Format Manual Protocol:

- ABI 9700 (with gold-plated or silver block)
- ABI Veriti
- ABI ProFlex
- Eppendorf Mastercycler pro S

**!** **IMPORTANT:** Always use the heated lid option when programming protocols. The PharmacoScan mPCR protocol was validated using the “9600 mode” on the ABI 9700, ABI Veriti, and ABI ProFlex thermal cyclers. The “Safe” mode was used for the Eppendorf Mastercycler pro S. Refer to the manufacturer’s user guide for programming information.



**▲** **WARNING:** Evaporation during denaturation can negatively impact assay performance. Use the recommended thermal cycler consumables and sealing film to eliminate condensation and evaporation.

PharmacoScan Assay 24-Array Format Manual Protocol has been validated with the ABI 9700 (with gold-plated or silver block) ABI Veriti, ABI ProFlex, and Eppendorf Mastercycler pro S. Use of other thermal cyclers for this stage may result in assay failure and may violate the array and reagent replacement policy.

## Thermal Cycler Consumables

Table 1.5 provides details into the consumables to be used with each thermal cycler.

**Table 1.5** Thermal Cycler Consumables for the PharmacoScan Assay 24-Array Format Manual Protocol

Thermal Cycler Model	PCR Plate Type	Seal*
ABI 9700	■ BioRad Hard-Shell Full-Height 96-Well Semi-Skirted PCR Plate (P/N HSS-9641)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
ABI Veriti	■ BioRad Hard-Shell Full-Height 96-Well Semi-Skirted PCR Plate (P/N HSS-9641)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
ABI ProFlex	■ BioRad Hard-Shell Full-Height 96-Well Semi-Skirted PCR Plate (P/N HSS-9641)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)
Eppendorf Mastercycler pro S	■ BioRad Hard-Shell Low Profile 96-well Full-Skirt PCR Plate (P/N HSP-9631)	MicroAmp Clear Adhesive Film from Applied Biosystems (P/N 4306311)

\*Microseal 'B' film from BioRad (P/N MSB-1001) may be used in place of MicroAmp Clear Adhesive Film for the ABI thermal cycler.

## Plate Centrifuge

One plate centrifuge is required for the PharmacoScan Assay 24-Array Format Manual Protocol. We recommend the plate centrifuges listed in Table 1.6. When centrifuging and drying pellets, the centrifuge must be able to spin down plates at:

- rcf: 3200 xg with an appropriate rotor - bucket combination (4000 RPM for the Eppendorf 5810R configuration described below)
- Temperature: 4°C

In addition, the bottom of the rotor buckets should be soft rubber to ensure that the deep well plates do not crack. Do not use buckets where the plates sit directly on a metal or hard plastic bottom. For the Eppendorf 5810R, do not use the A-4-62 rotor with a WO-15 plate carrier (hard bottom).

**Table 1.6** Plate Centrifuges Recommended for the PharmacoScan Assay 24-Array Format Manual Protocol

✓	Item	Supplier	Part Number
□	One of the following centrifuges:		
■	Allegra® 25R Refrigerated Benchtop Centrifuge	Beckman Coulter	369434 (60 Hz, 280 V) 369435 (50/60 Hz, 200 V) 369436 (50 Hz, 230 V)  368954 S5700 Microplate rotor
■	Centrifuge 5810R (refrigerated)	Eppendorf	022625501 (60 Hz, 120 V) 022625101 (50 Hz, 120 V, 20 AMP version) 022625551 (50 Hz, 230 V)
□	A-4-81 MTP/Flex swinging bucket rotor with 4 microtest-plate buckets		022638807
■	Sorvall® Legend® XTR (refrigerated), with:	Thermo Scientific	75004521 (60 Hz, 120 V) 75004520 (50 Hz, 230 V) 75004523 (50/60 Hz, 230 V USA and Canada)
□	TX-750 high visibility swinging rotor bucket		75003607
□	Carrier for microplates (including plate trays and neoprene pads)		One of either: ■ Set of two carriers (75003795) ■ Set of four carriers (75003617)



Relative centrifugal force (*rcf*) can be calculated as follows:

$$rcf = (1.118 \times 10^{-5}) R S^2$$

where R is the radius of the rotor in centimeters, and S is the speed of the centrifuge in revolutions per minute.

## Sample Plates

**Table 1.7** Sample Plates Required for PharmacoScan Assay 24-Array Format Manual Protocol

✓	Plate Description	Manufacturer/ Distributor	Part Number
☐	One of the following PCR Plates: <ul style="list-style-type: none"> <li>■ BioRad Hard-Shell Full-Height 96-Well Semi-Skirted PCR Plate (if using ABI 9700, ABI Veriti, ABI ProFlex thermal cyclers)</li> <li>■ BioRad Hard-Shell Low Profile 96-well Full-Skirt PCR Plate (if using Eppendorf Master pro S thermal cyclers)</li> </ul>	Bio-Rad	HSS-9641
☐	ABgene® 96 Square Well Storage Plate Mark II, 2.2 mL	Thermo Scientific	AB-0932
☐	96-well UV Star Plates, 370 µL/well	E&K Scientific VWR, Sigma-Aldrich	25801 655801

## Pipettes and Tips

Pipettes and tips recommended for PharmacoScan target preparation are listed in [Table 1.8](#).

**Table 1.8** Recommended Pipettes and Tips for PharmacoScan Assay 24-Array Format Manual Protocol

✓	Equipment	Manufacturer/ Distributor	Part Number
☐	Pipet-Lite™, Magnetic-Assist single channel P20	Rainin	L-20
☐	Pipet-Lite™, Magnetic-Assist single channel P200	Rainin	L-200
☐	Pipet-Lite™, Magnetic-Assist single channel P1000	Rainin	L-1000
☐	Pipette, 12-channel P20	Rainin	L12-20
☐	Pipette, 12-channel P50 (optional)	Rainin	L12-50
☐	Pipette, 12-channel P200	Rainin	L12-200
☐	Pipette, 12-channel P1200	Rainin	L12-1200
☐	Pipette tips GP = refill	Rainin	GP-L10F
☐	Pipette tips GP = refill	Rainin	GP-L200F
☐	Pipette tips GP = refill	Rainin	GP-L1000F
☐	Pipette tips RT = with rack	Rainin	RT-L10F
☐	Pipette tips RT = with rack	Rainin	RT-L200F
☐	Pipette tips RT = with rack	Rainin	RT-L1000F

## Other Labware Required

**Table 1.9** Other Labware Required for the PharmacoScan Assay 24-Array Format Manual Protocol

✓	Item	Manufacturer/ Distributor	Part Number
☐	Microcentrifuge tubes and tube holder	various	various
☐	8 well strip tubes with caps	various	various
☐	50 mL conical centrifuge tube, polypropylene, sterile	various	various
☐	15 mL conical centrifuge tube, polypropylene, sterile	various	various
☐	Tube Holder, 15 and 50 mL	various	various
☐	Matrix™ 25 mL divided reservoir	Thermo Fisher Scientific	8095
☐	96 well plate metal chamber	Diversified Biotech	CHAM-1000
☐	Serological Pipettes (10 mL)	VWR	89130-898
☐	Serological Pipettes (5 mL)	VWR	89130-896
☐	Adhesive film for 96-well plates – use one of the following: <ul style="list-style-type: none"> <li>■ MicroAmp® Clear Adhesive Film</li> <li>■ Microseal 'B' Film</li> </ul>	Applied Biosystems Bio-Rad	4306311 MSB1001
☐	Kimwipes®	Your choice	—

## Other Reagents and Gels Required

**Table 1.10** Other Reagents Required

✓	Item	Supplier	Part Number
☐	QIAGEN Multiplex PCR <i>Plus</i> Kit (100)	QIAGEN	206152
☐	Quant-iT PicoGreen dsDNA Assay Kit (recommended for DNA quantitation)	Thermo Fisher Scientific	P7589
☐	Reduced EDTA TE Buffer (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA)	Affymetrix	75793
☐	2-Propanol, anhydrous, 99.5% (Isopropanol)	Sigma-Aldrich	278475

**Table 1.11** Reagents and Gels Required to Run QC Steps

✓	Item	Supplier	Part Number
☐	Mother E-Base™ Device	Thermo Fisher Scientific (formerly Life Technologies™ /Invitrogen)	EB-M03
☐	Daughter E-Base™ Device		EB-D03
☐	E-Gel® 48 4% agarose gels (for Sample QC)		G8008-04
☐	TrackIt™ 25 bp DNA Ladder (for Sample QC)		10488-022
☐	TrackIt™ Cyan/Orange Loading Buffer (for Sample QC)		10482-028
☐	E-Gel® 48 1% agarose gels (for gDNA QC)		G8008-01
☐	RediLoad™ (for gDNA QC)		750026
☐	E-Gel® 96 High Range DNA Marker (for gDNA QC)		12352-019
☐	E-Gel® 48 2% agarose gels (optional mPCR QC)		G8008-02
☐	Water, nuclease-free, ultrapure MB grade		Affymetrix
☐	50 bp DNA ladder (optional mPCR QC)	New England Biolabs Inc.	N3236S

## Other Lab Equipment Required

The remaining equipment required is listed in [Table 1.12](#).

**Table 1.12** Other Equipment Required

✓	Item	Supplier	Part Number
☐	Freezer, -20°C	Any vendor	—
☐	Refrigerator, 2 to 8°C	Any vendor	—
☐	Vortex-Genie® (for plates and microtubes)*	Scientific Industries	SI-0236 (120 V/60 Hz) SI-0246 (230 V/50 Hz)
☐	Mini Microcentrifuge, for 2 mL tubes and PCR strip tube adapter*	VWR	93000-196 (120V) 93000-196 (230V)
☐	Bel-Art Cryo-Safe Mini Cooler, -15°C*	VWR	47751-730
☐	Ice bucket, 4 to 9 liters	Any vendor	—
☐	Markers, permanent, fine point	Your choice	—
☐	Pipet aid*	VWR	53106-220
☐	Zerostat Anti-static Gun	Affymetrix	74-0014

\* Equivalent items from other manufacturers are acceptable.

# Appendix A

## Supplier Contact List

**Table A.1** Supplier Contact List

<b>Supplier</b>	<b>Website Address</b>
Affymetrix	<a href="http://www.affymetrix.com">www.affymetrix.com</a>
Applied Biosystems	<a href="http://www.appliedbiosystems.com">www.appliedbiosystems.com</a>
Beckman Coulter	<a href="http://www.beckmancoulter.com">www.beckmancoulter.com</a>
Bio-Rad	<a href="http://www.bio-rad.com">www.bio-rad.com</a>
E&K Scientific	<a href="http://www.eandkscientific.com">www.eandkscientific.com</a>
Eppendorf	<a href="http://www.eppendorf.com">www.eppendorf.com</a>
New England BioLabs Inc.	<a href="http://www.neb.com">www.neb.com</a>
QIAGEN	<a href="http://www.qiagen.com">www.qiagen.com</a>
Rainin	<a href="http://www.shoprainin.com">www.shoprainin.com</a>
Scientific Industries	<a href="http://www.scientificindustries.com">www.scientificindustries.com</a>
Sigma-Aldrich	<a href="http://www.sigmaaldrich.com">www.sigmaaldrich.com</a>
Thermo Fisher Scientific (Thermo Scientific, Applied Biosystems, Invitrogen, Fisher Scientific)	<a href="http://www.thermofisher.com">www.thermofisher.com</a>
USB Corporation (Affymetrix)	<a href="http://www.usb.affymetrix.com">www.usb.affymetrix.com</a>
VWR	<a href="https://us.vwr.com/store/">https://us.vwr.com/store/</a>
Biometra	<a href="http://www.biometra.de">www.biometra.de</a>