

KENNEN SIE...

ΧΡΩΜΑΤΟΓΡΑΦΙΑ
ΧΡΩΜΑΤΟΓΡΑΦΙΑ

AGILENT CHROMATOGRAPHIE
Seminar Tour 2013

WILLKOMMEN

zum Agilent Chromatographie-Seminar

In BERLIN

27. Februar 2013

AGENDA – 27. Februar 2013 - Berlin

09:00 – 09:15	Begrüßung - Einführung, Agenda und Logistik Joachim Werner, Agilent Technologies
09:15 – 09:45	Neue Agilent LC Produkte für unterschiedliche Einsatzgebiete Sebastian Krahe / Agilent Technologies
09:45 – 10:15	Tücken der UHPLC - Wissenswertes im Bereich < 2 Mikronpartikel Ulrike Jegle / Agilent Technologies
10:15 – 10:45	Säulen- und Phasenauswahl in der (U) HPLC - muss es immer vollporöses Material sein? Ulrike Jegle / Agilent Technologies
10:45 – 11:00	Kaffeepause
11:00 – 11:45	HPLC: Tipps zur Wartung, Fehlervermeidungen, Problemlösung Christoph Menzel / Agilent Technologies
11:45 – 12:30	Vorstellung: neuer GC7890B, neues GC-MSD5977 Claudia Wolf / Agilent Technologies
12:30 – 13:30	Mittagessen – Buffet



AGENDA – 27. Februar 2013 - Berlin

Parallel:	Vorstellungen zur GC (RAUM HUMBOLD)	HPLC WORKSHOP (HIER)
13:30 – 14:30	Neues aus der Gaschromatographie Claudia Wolf / Agilent Technologies	WORKSHOP HPLC – Schwerpunkt "Einführung in tägliche Wartungsarbeiten" Christoph Menzel / Agilent
14:30 – 15:30	CSD: Hocheffiziente und hochsensitive GC Ulrike Jegle / Agilent Technologies	
15:30 – 15:45	Kaffeepause	
15:45 – 16:15	LabAdvisor- ein leistungsfähiges Werkzeug zur Überprüfung Ihrer LC Sebastian Krahe / Agilent Technologies	
16:15 – 16:30	Zusammenfassung	



A Brief History of Agilent

Our journey so far

1939–1998: The Hewlett-Packard years

60-year heritage of leadership and innovation

1999: Agilent is born

Biggest IPO in Silicon Valley: \$2.1 billion

2000–2004: Growing pains

Survived dot-com crash via major restructuring

2005: The new Agilent

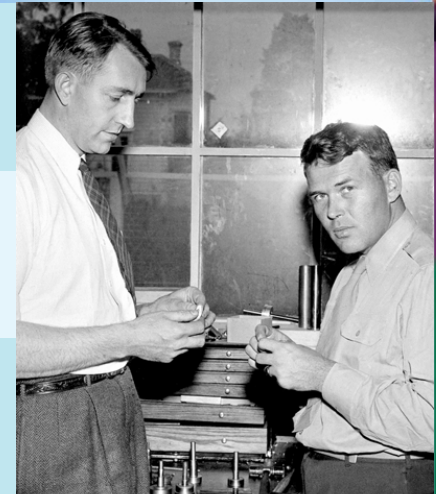
Focus on measurement

2006–2010: Transformation

Acquisitions, divestitures and restructuring

2011: The world's premier measurement company

\$6.6 billion revenue: best operational performance in 11-year history



Agilent at a Glance

REVENUE IN FY11

\$6.6 billion (more than 70% generated outside the U.S.)

EMPLOYEES

20,000

PRESIDENT AND CEO

William P. (Bill) Sullivan

CUSTOMER LOCATIONS

100+ countries

MANUFACTURING AND
R&D LOCATIONS

U.S., Europe, Asia Pacific

NYSE

A



Agilent Technologies

Addressing critical measurement challenges

FY11 Revenue: \$6.6B



Electronic Measurement Group

- Wireless technologies
- Mobile phone R&D and manufacturing
- Aerospace/defense systems and maintenance
- Computer & industrial research & manufacturing

**FY11 Revenue
\$3.3B**



Chemical Analysis Group

- Food safety, quality
- Energy research, production
- Quality of air, water, soil
- Forensics, drugs of abuse

**FY11 Revenue
\$1.5B**



Life Sciences Group

- Pharmaceutical research and manufacturing
- Academic and government research

**FY11 Revenue*
\$1.5B**



Diagnostics and Genomics Group

- Cancer diagnostics
- Cytogenetics research
- Genomics tools for molecular analysis

**FY11 Revenue*
\$277M**

Agilent Research Laboratories
Enabling technology breakthroughs across Agilent

**DGG segment created
in Q312; FY11 revenue
restated*



Agilent Research Laboratories

Instrumental to our future

Applied research

Breakthrough technologies

Measurement synergies

World-class research teams, in core and emerging markets



Agilent's Core Technologies

Communications/Electronics

Life Sciences / Chemical Analysis

- Measurement science
- Electronic circuit and systems design
- Applications software and solutions integration

- Radio frequency/microwave/wireless simulation and test
- Digital design validation and test
- General-purpose instrumentation
- Parametric test solutions
- Manufacturing test solutions

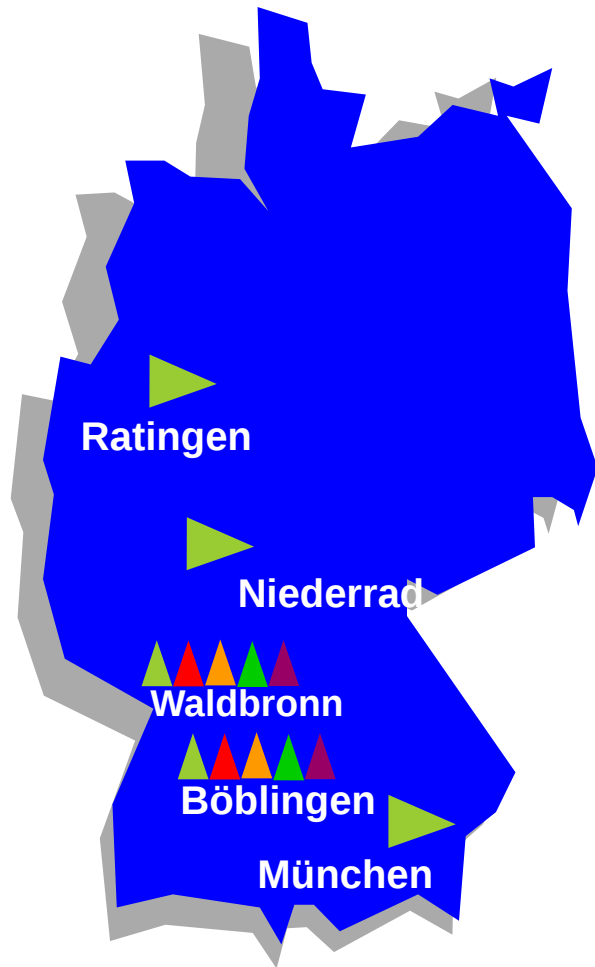
- Liquid and gas chromatography
- Mass spectrometry
- Microfluidics
- Microarrays
- Reagents chemistry
- LC and GC columns
- Software and informatics
- Lab automation
- Nuclear magnetic resonance
- Vacuum technology
- X-ray
- Atomic spectroscopy

Agilent Research Laboratories



Agilent Technologies in Germany

Locations



Employees: 1,400



Waldbronn:

- Pharma & Biotech Solutions
- Liquid Phase Separations
- LSCA Order Fulfillment WAD
- EMEA Field Support Ctr.

- Sales
- Research & Development
- Manufacturing
- Marketing
- Services

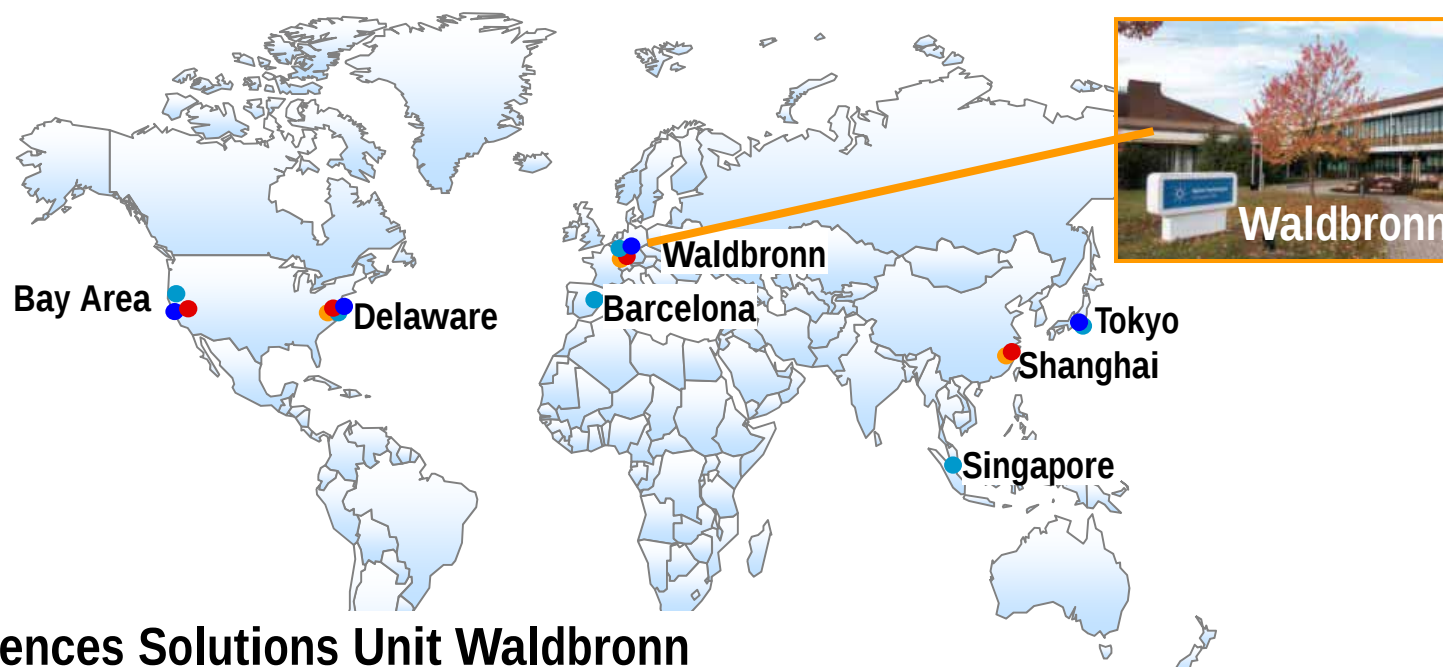
Boeblingen:

- Digital Verification Solutions
- Photonics & Network Test
- GIO Infrastructure Services
- EMG Field Germany



Life Sciences and Chemical Analysis

Waldbronn Site



Life Sciences Solutions Unit Waldbronn

- Liquid Phase Separations Division
- Pharma & Biotech Solutions

Waldbronn Order Fulfillment

- Liquid Phase Separations Platforms
- Pharma & Biotech Solutions
- LSCA Logistics Center Europe

EMEA Field Operations

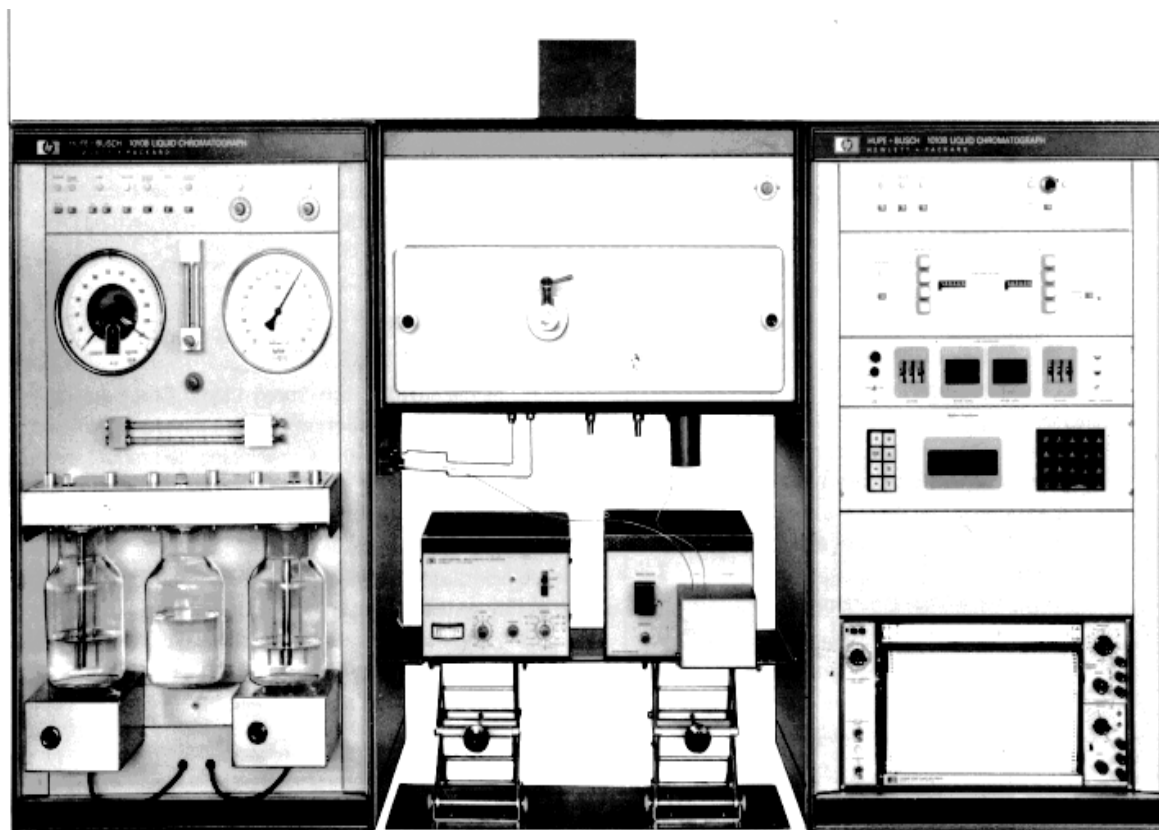
- Field Support Center
- Customer Care & Training Centers
- Repair Center
- German Field Operations





Acquisition of
Hupe & Busch
by Hewlett-Packard

1973



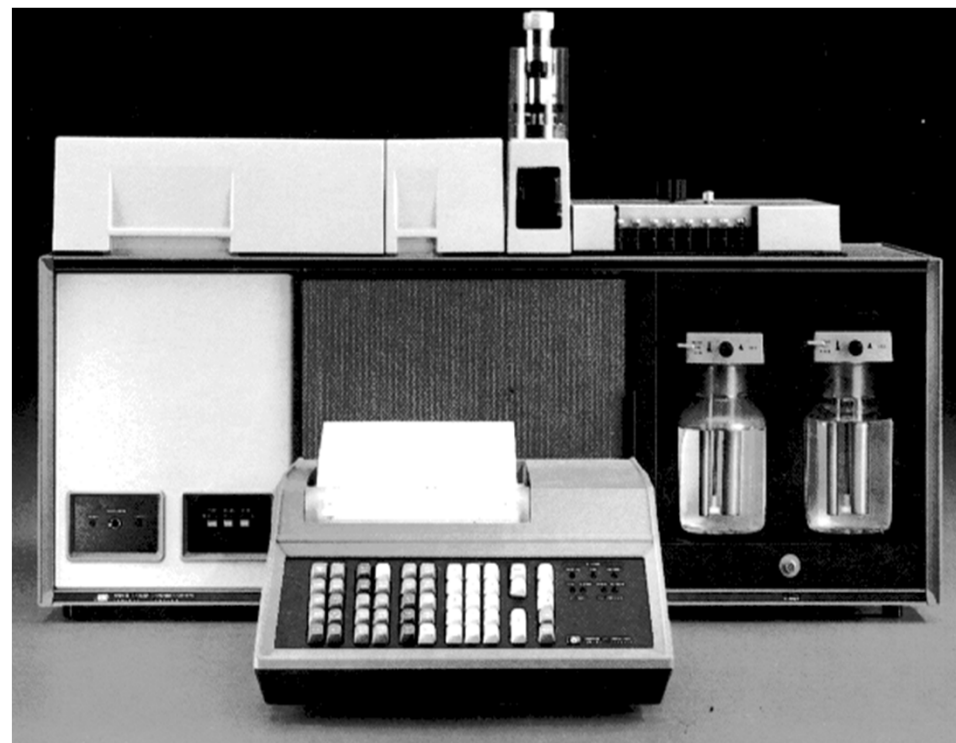
HP 1010B Liquid Chromatograph





Move to Waldbronn Building 1

1978



**HP 1084B HPLC System
with microprocessor control**



**New Sales Office
Now: EMEA Field Support Ctr.**



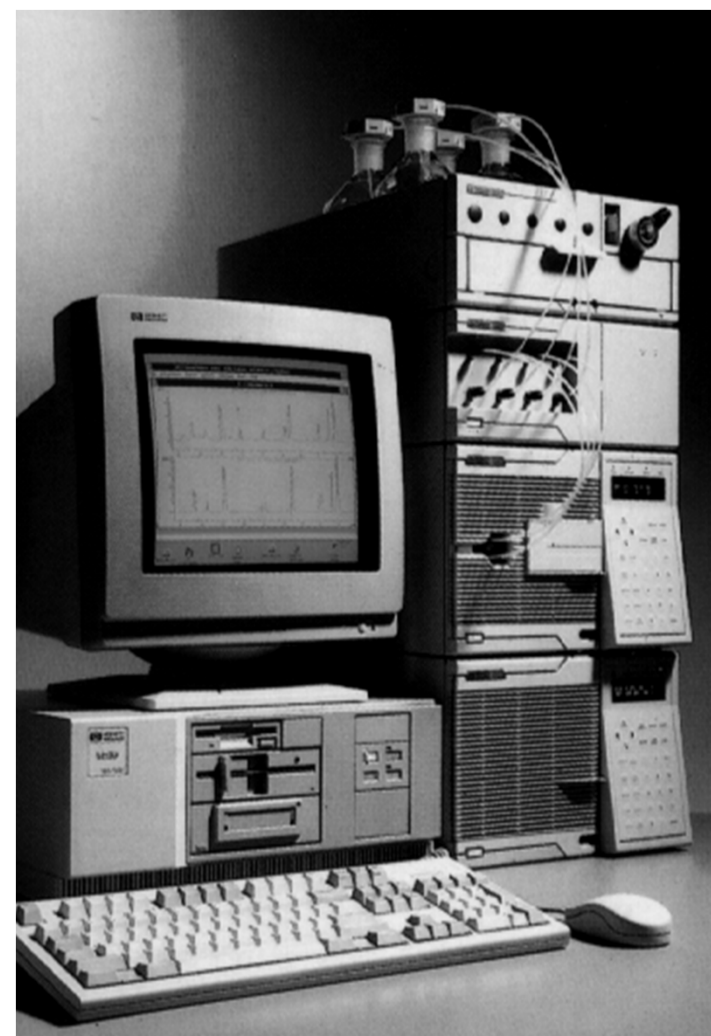
**HP 1090
High Performance HPLC System**

1983



1988

Expansion – Building 2



**HP 1050
Modular HPLC System**



1985



**HP ChemStation
PASCAL**

1993



**Agilent ChemStation
MS WINDOWS**

2002



Cerity Pharma QA / QC



1993

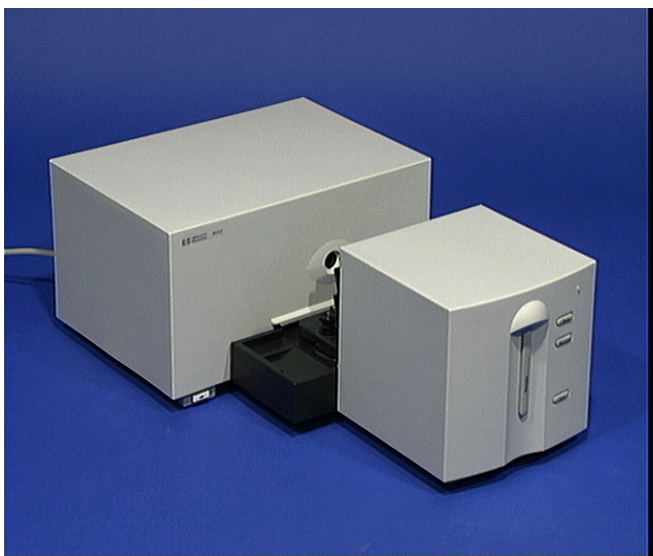


Agilent 3D Capillary Electrophoresis System



1995

Agilent 8453 UV / Vis System



Agilent 1100 HPLC Series



1999



Agilent 2100 Bioanalyzer

2005



Agilent LC-Chip / MS System



Expansion - Building #4



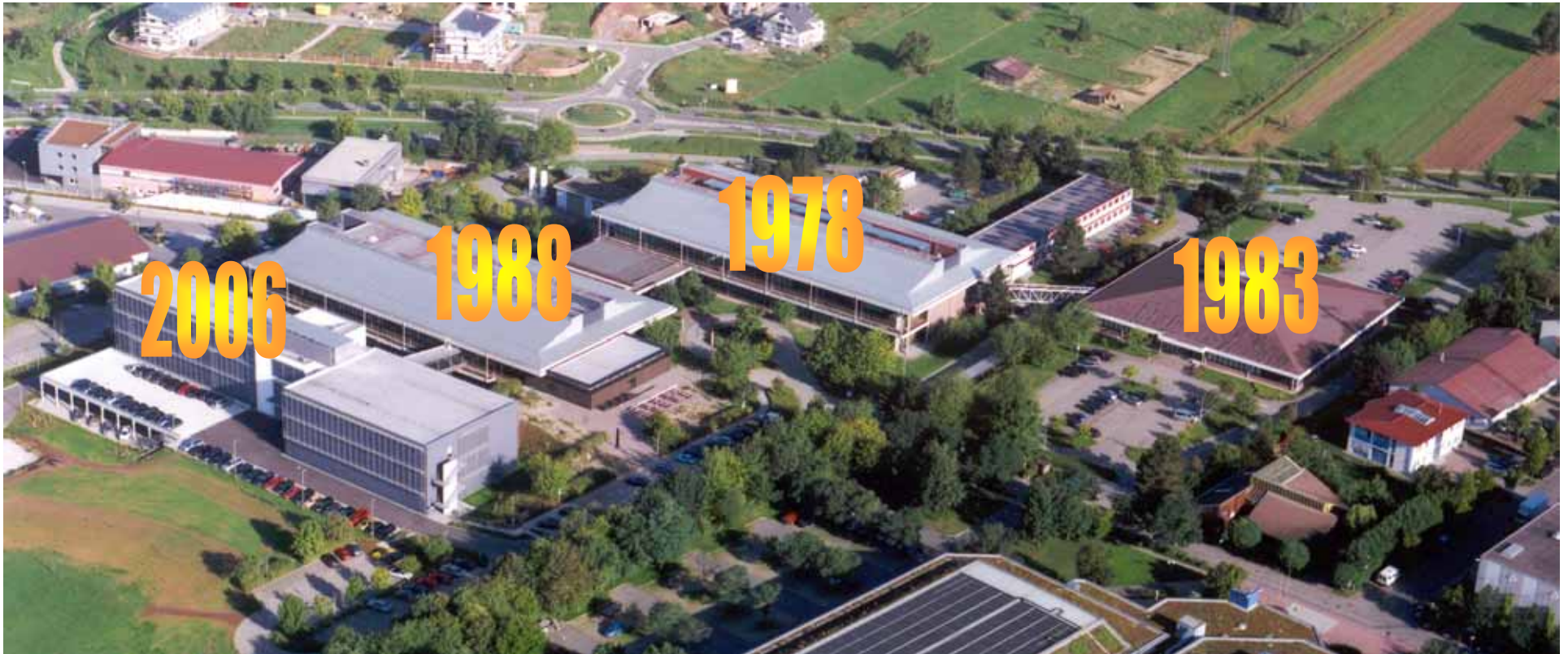
2006



**Agilent 1200 LC Series
Incl. Rapid Resolution LC**



LSCA Waldbronn Site: A History of Growth . . .



Agilent Acquisitions

Extensions of our measurement solutions

Chemical Analysis

Yokogawa Analytical Systems

Chemical analysis
Sulfur/nitrogen detectors

GE Chemiluminescence

Particle size and zeta potential

Colloidal Dynamics

GC & GC/MS workflow

RVM Scientific

Particle characterization

Particle Sizing Systems

Life Sciences

Silicon Genetics

Comparative genomic hybridization

Computational Biology

Location analysis

Scientific Software

Lab informatics software

SynPro Corp

Stratagene

Synthetic nucleic acid

Kalabie

Life sciences and diagnostics

Electronic Laboratory Notebook

Velocity11

Life science workflow solutions

Varian

Spectroscopy, NMR, x-Ray, Vacuum

Dako

Clinical Diagnostics



Agilent Portfolio

Life Sciences and Chemical Analysis



MR Imaging



Prep LC



NMR



Bravo Automated
Liquid Handling
Platform



Microarray



Capillary Electrophoresis,
Bioanalyzer portfolio



Dissolution



Ion Trap



X-Ray



LC Q-TOF



LC-Technologies



LC-QQQ



Columns



UV/VIS



ICP OES



GC - Technologies



ICP MS QQQ



GC QQQ



GC qTOF



Vacuum Technologies



Agilent Technologies

KENNEN SIE...

ΧΡΩΜΑΤΟΓΡΑΦΙΑ
ΧΡΩΜΑΤΟΓΡΑΦΙΑ

AGILENT CHROMATOGRAPHIE
Seminar Tour 2013

Bitte ...

Geben Sie am Ende der Veranstaltung
einen Feedback-Bogen ab. Danke!

AGENDA – 27. Februar 2013 - Berlin

- | | |
|---------------|---|
| 09:00 – 09:15 | Begrüßung - Einführung, Agenda und Logistik
Joachim Werner, Agilent Technologies |
| 09:15 – 09:45 | Neue Agilent LC Produkte für unterschiedliche Einsatzgebiete
Sebastian Krahe / Agilent Technologies |
| 09:45 – 10:15 | Tücken der UHPLC - Wissenswertes im Bereich < 2 Mikronpartikel
Ulrike Jegle / Agilent Technologies |
| 10:15 – 10:45 | Säulen- und Phasenauswahl in der (U) HPLC - muss es immer vollporöses Material sein? Ulrike Jegle / Agilent Technologies |
| 10:45 – 11:00 | Kaffeepause |
| 11:00 – 11:45 | HPLC: Tipps zur Wartung, Fehlervermeidungen, Problemlösung
Christoph Menzel / Agilent Technologies |
| 11:45 – 12:30 | Vorstellung: neuer GC7890B, neues GC-MSD5977
Claudia Wolf / Agilent Technologies |
| 12:30 – 13:30 | Mittagessen – Buffet |

