



Zürich, March 11th 2014



Full Year Results 2013

Analyst & Media Conference



Speakers



Dr. David Martyr, CEO

Dr. Rudolf Eugster, CFO

Agenda

- Introduction Dr. David Martyr
- Financial Results Dr. Rudolf Eugster
- Priorities Dr. David Martyr
- Outlook & Summary Dr. David Martyr
- Questions & Answers

Financial Performance 2013

- Financial objectives set at start of the year only partly met
- Sales growth of +0.1% in local currencies below original expectations
 - One clear cause: lower Life Sciences Business instrument sales in established markets in tough market environment
 - Continued solid growth in Partnering Business, despite faster than anticipated phase out of products
 - Double-digit growth in both segments in China and in consumables business
- Profitability target achieved
 - EBIT margin improved by 60 basis points
 - Net profit margin increased by 100 basis points
- Improved cash flow from operating activities



Operational Highlights 2013

- Important progress in the implementation of strategic priorities
- Dako Omnis development (P16) concluded
 - Instrument successfully launched by partner Dako in September 2013
- Good progress in other development programs with launch of ORTHO Vision™ (P14) in Partnering Business and next generation liquid handling platform in Life Sciences Business expected in 2014
- Strengthening of management team with three new members of the Management Board and additional targeted appointments at different levels of organization and in regions

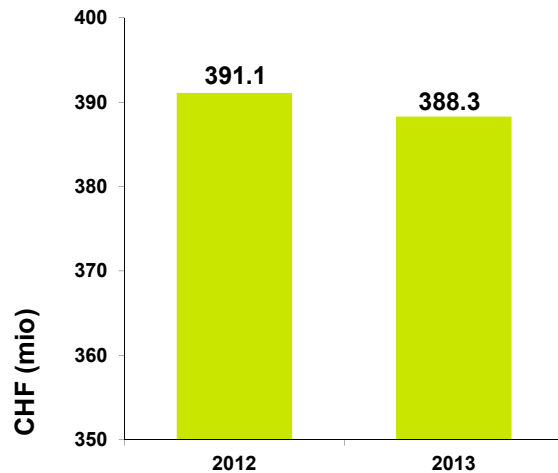


Agenda

- Introduction Dr. David Martyr
- Financial Results Dr. Rudolf Eugster
- Priorities Dr. David Martyr
- Outlook & Summary Dr. David Martyr
- Questions & Answers

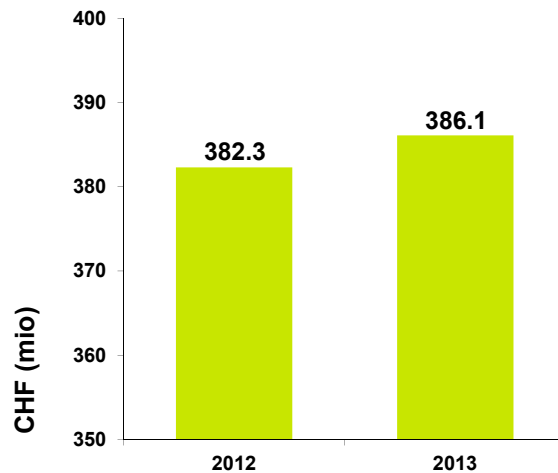
2013 Sales Performance and Order Entry

Sales FY 2013



- Sales FY 2013 vs. FY 2012
 - -0.7% in Swiss Francs
 - +0.1% in local currencies
- Sales H2 2013 vs. H2 2012
 - -1.1% in Swiss Francs
 - +0.8% in local currencies

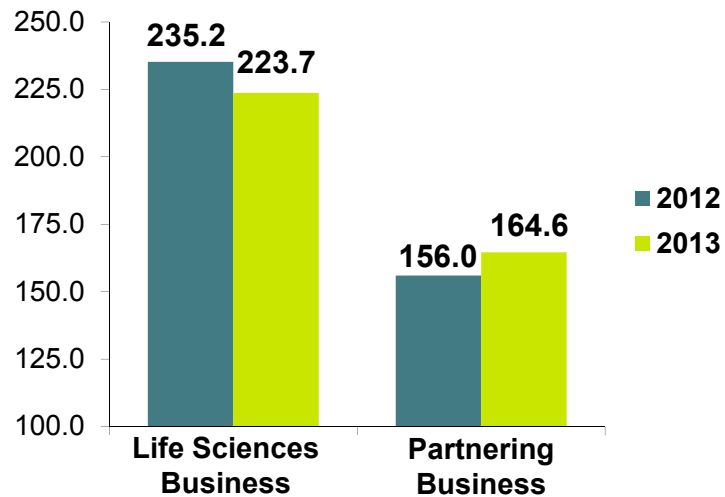
Order Entry FY 2013



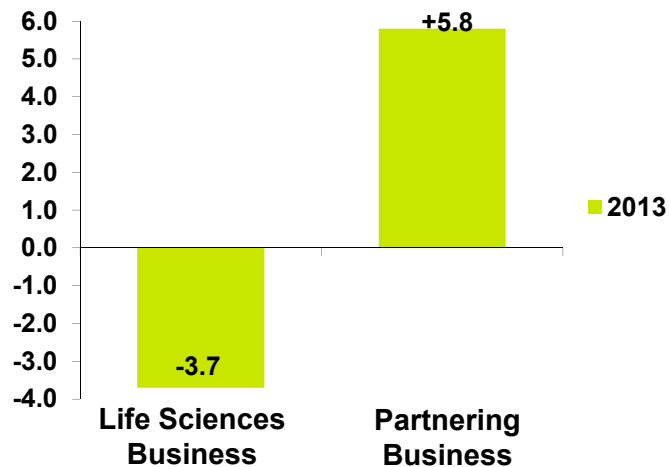
- Order Entry FY 2013 vs. FY 2012
 - +1.0% in Swiss Francs
 - +1.9% in local currencies
- Order Entry H2 2013 vs. H2 2012
 - -2.9% in Swiss Francs
 - -1.0% in local currencies

2013 Segment Sales

Sales (in CHF millions)



Sales in local currency (change in %)



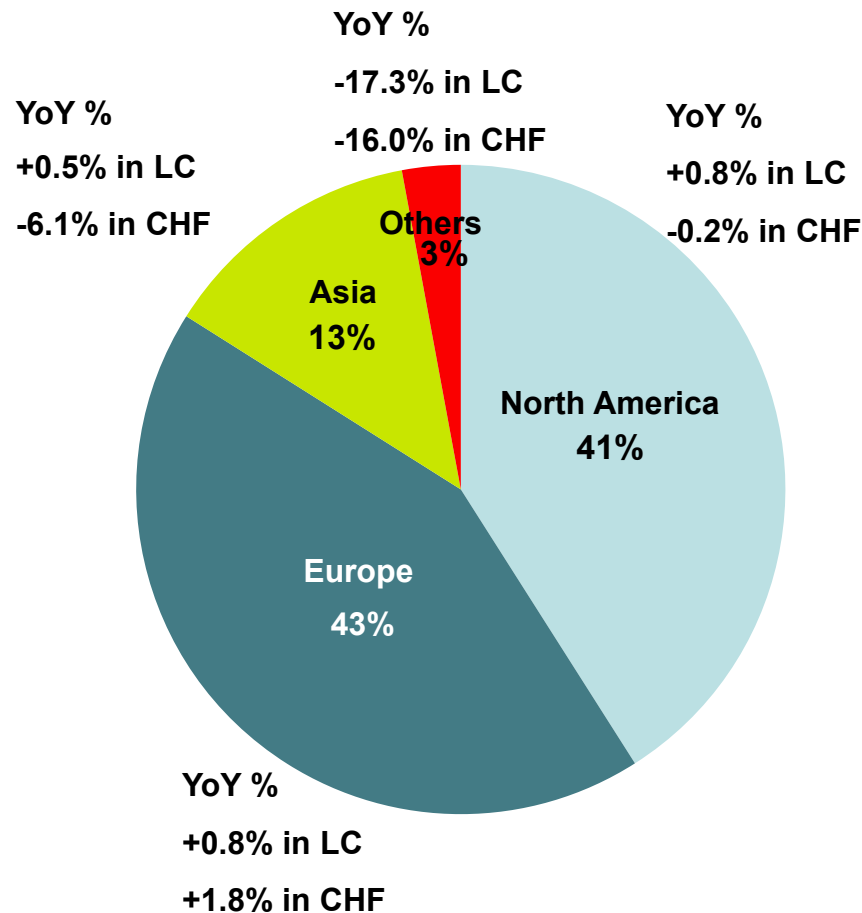
• Life Sciences Business:

- FY 2013: Sales 4.9% below prior-year in CHF and 3.7% lower in local currencies
- Decline largely due to fewer instruments sold in established markets
- Sales in China and of consumables recorded strong growth
- Order Entry also down, but exceeded sales
- H2 2013: Sales down by 6.1% in CHF and 3.8% lower in local currencies

• Partnering Business:

- FY 2013: Sales up by 5.5% in CHF and 5.8% in local currencies
- Strong growth in Services & Consumables, good performance in Components, instruments only slightly down
- Order Entry also saw good growth
- H2 2013: Sales accelerated further and increased by 7.8% in CHF and 8.9% in local currencies

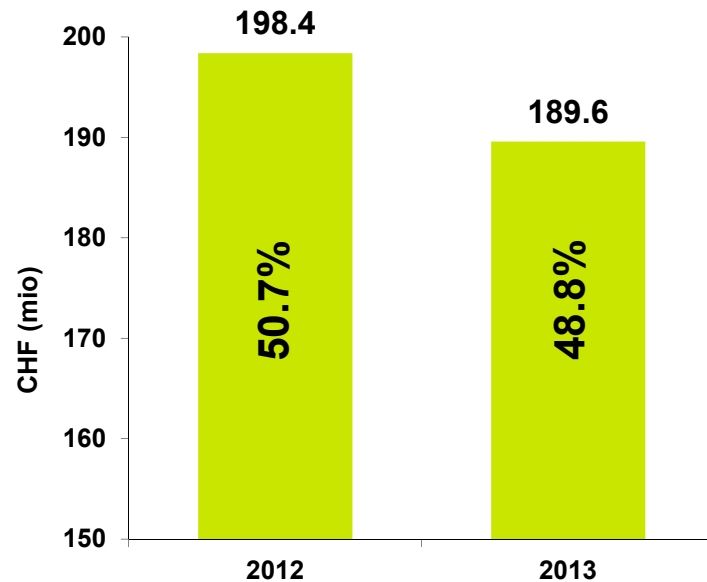
2013 Regional Sales Development



- Europe: growth in the Partnering Business accelerated further in the second half, driven by instrument deliveries to Dako; lower Life Sciences Business sales due to continuing challenging economic situation
- North America: Partnering Business increased, driven by the components business; Life Sciences Business sales below the prior-year as public budget cuts and the government shutdown in October had a negative impact
- Asia: both segments recorded clearly double-digit sales growth in China; declining Life Sciences Business sales in Japan and Asia-Pacific

FY 2013 Gross Profit

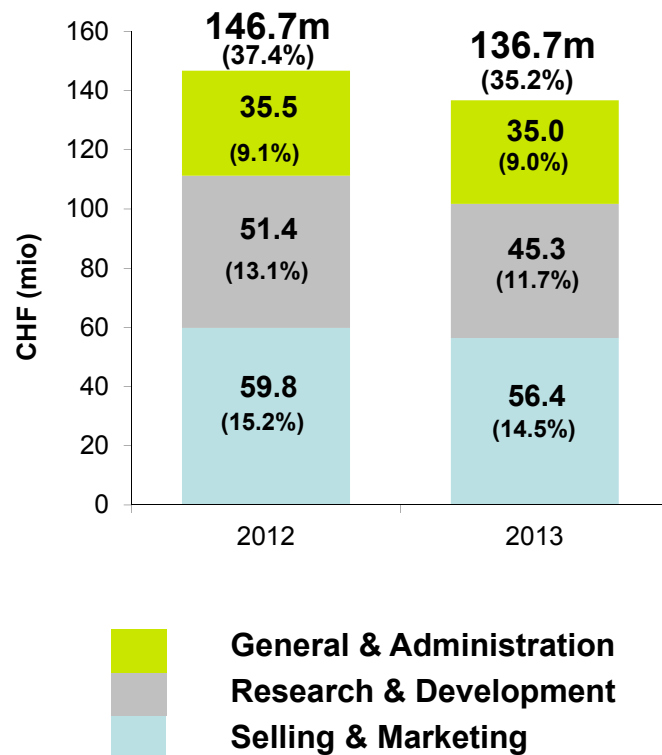
Gross Profit (% = as a % of sales)



- Gross profit decreased to CHF 189.6m
 - CHF 8.8m or 4.5% below 2012
- Gross profit margin at 48.8%
 - Product mix impact: more sales from Partnering Business and Services & Consumables with lower GP margins and less sales from higher margin Life Sciences Business instruments
 - Slight price increases
 - Less costs of OEM development programs offset by medical device tax and higher non-standard cost of sales

FY 2013 Cost Structure

Operating Expenses (% = as a % of sales)

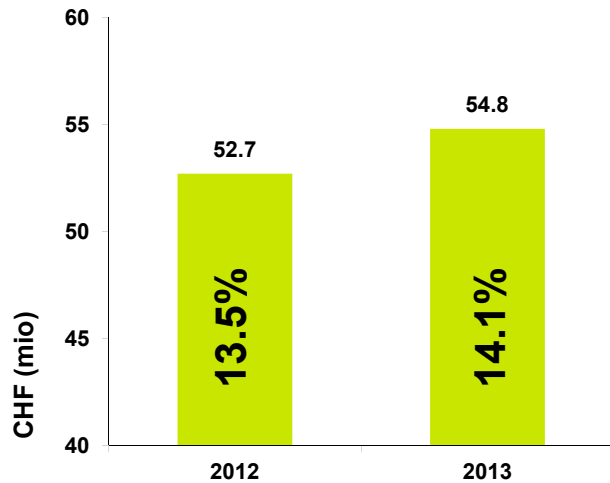


- Costs decreased by CHF 10.0m (-6.8%)

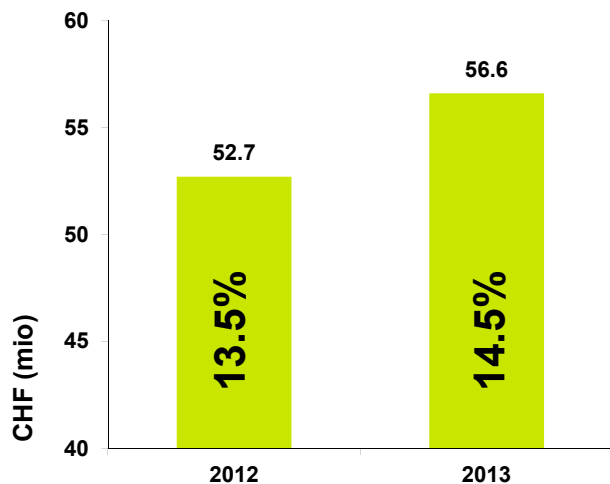
- Selling & Marketing
 - Less costs, despite increased investments in sales organization (especially in China)
- Research & Development
 - Net R&D decreased by 11.8% to CHF 45.3m
 - Net R&D expenses down to 11.7% of sales
 - Gross R&D investments decreased by 9.2% to CHF 104.1m (2012: CHF 114.6m),
 - Gross R&D includes CHF 51.2m for OEM partners and CHF 10.2m capitalized R&D costs (2012: CHF 4.0m)
- General & Administration
 - Decreased by 1.6% with less cost on corporate level

FY 2013 EBIT

EBIT (% = as a % of sales)



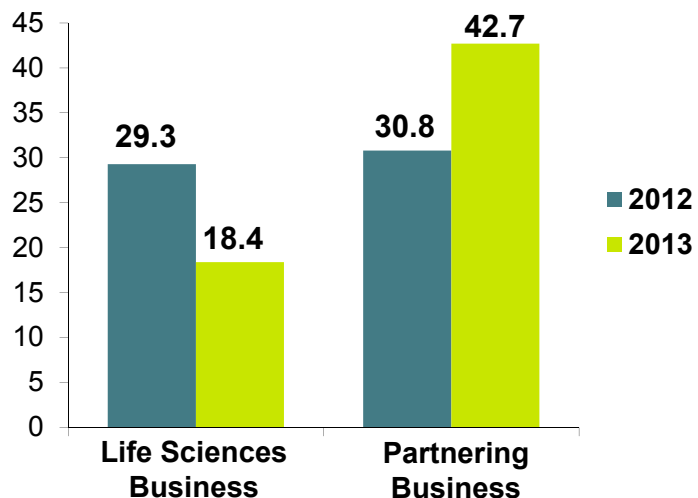
EBIT 2013 at 2012 constant currencies



- EBIT increased to CHF 54.8m
 - CHF 2.1m or 4.0% above 2012
- EBIT margin at 14.1%
 - 60 basis points higher, fulfilling target of improvement of “around 50 basis points”
 - Main effects impacting the EBIT margin:
 - (-) GP margin 190 bps lower
 - (+) Expenses 250 bps better
 - At 2012 constant exchange rates EBIT margin reached 14.5% of sales

FY 2013 Segment Profitability

EBIT (in CHF millions)



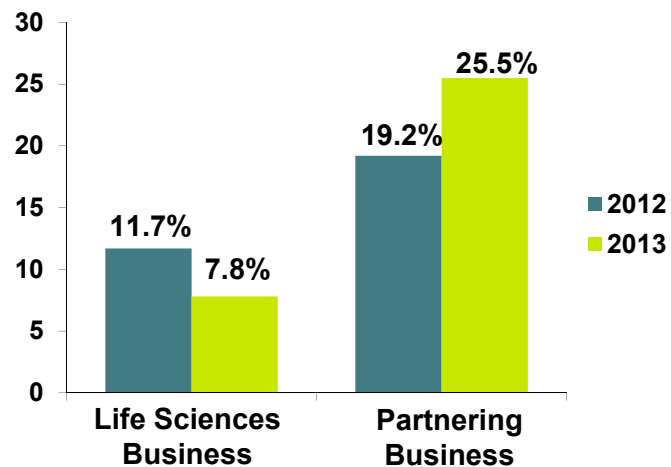
• **Life Sciences Business:**

- EBIT down by 37% to CHF 18.4m
- EBIT margin decreased to 7.8%
 - Significant negative volume and mix effect
 - Reduced R&D expenses, but negative FX and Medical Device Tax effect

• **Partnering Business:**

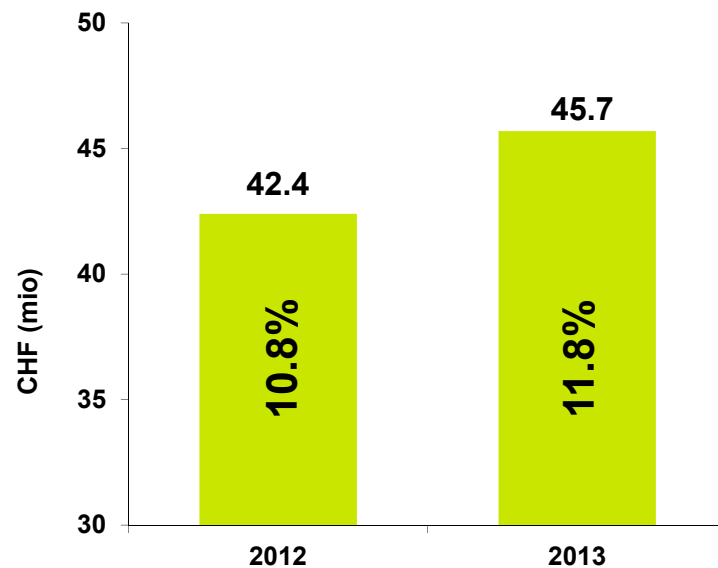
- EBIT increased by 39% to CHF 42.7m
- EBIT margin increased to 25.5% due to
 - Higher volumes
 - Lower engineering costs from development program and reversal of bad debt accrual

EBIT Margin as a % of sales



FY 2013 Net Profit

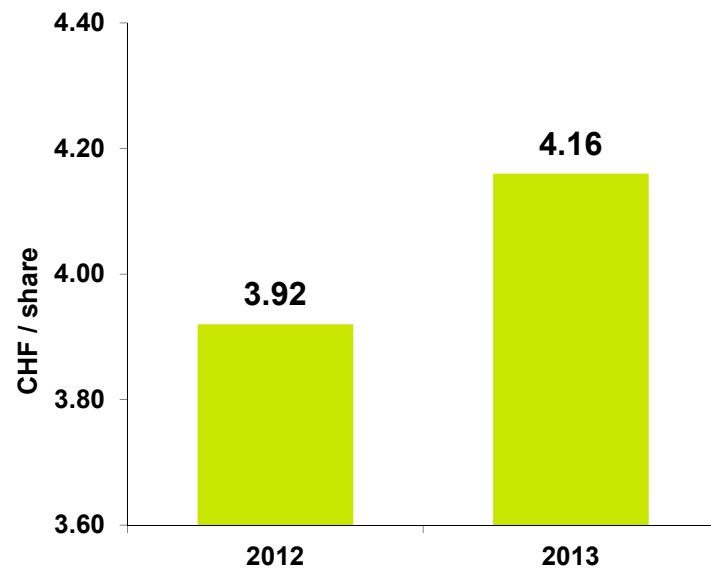
Net Profit (% = as a % of sales)



- Net profit of CHF 45.7m
 - CHF 3.3m or 7.8% above 2012
 - Positive effects below EBIT line: improved financial result and lower tax rate
- Net profit margin reached 11.8%
- Tax rate at 17.7% (2012: 19.6%)

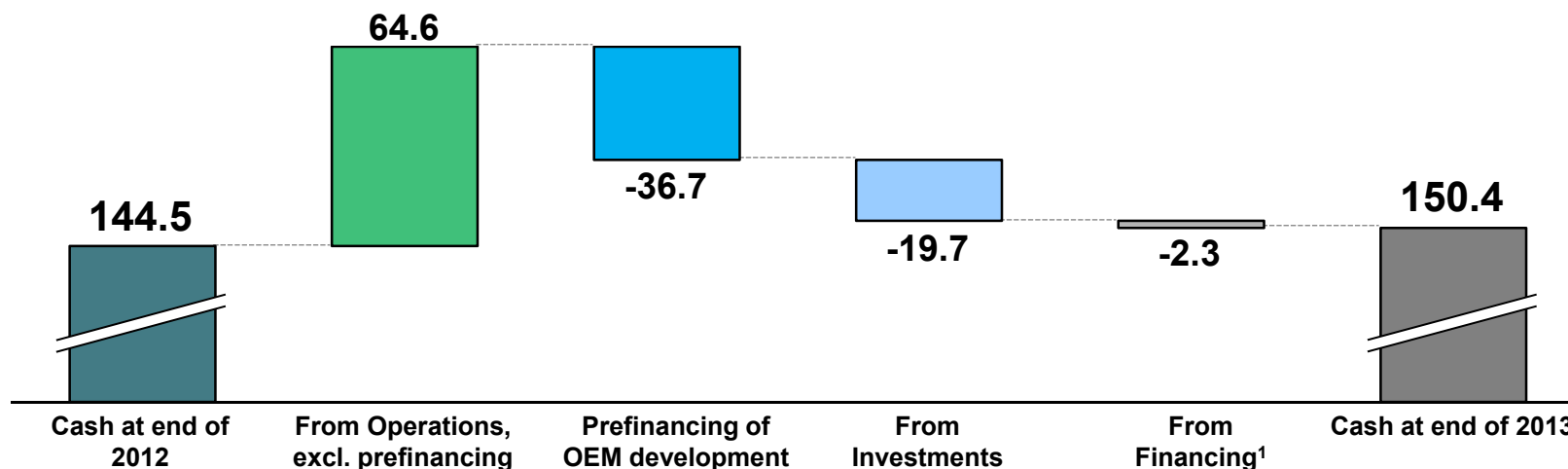
FY 2013 Basic Earnings per Share

Basic Earnings per Share



- Earnings per share increased by 6.1% to CHF 4.16
- Average number of shares at 11.0m (2012: 10.8m)
- Unchanged dividend of CHF 1.50 per share proposed

FY 2013 Cash Flow



- Cash flow from operations increased to CHF 27.9m (2012: CHF 2.4m)
 - Excluding prefinancing of CHF 36.7m for OEM development project, cash flow from operations reached CHF 64.6m
- Net working capital excluding pre-financing of OEM development stable
 - Days Sales Outstanding at 51 days (2012: 56)
- Investments of CHF 19.7m (2012: CHF 13.3m), exceed CHF 10.3m for amortization & depreciation (2012: CHF 10.3m)
- Cash flow from financing activities includes dividend payments of CHF 16.5m and proceeds of CHF 10.8m from sales of treasury shares (mainly due to exercise of options)
- Net liquidity² increased to CHF 143.4m (Dec. 31, 2012: CHF 141.3m)

¹ Includes translation differences of CHF -0.1m

² Net Liquidity = cash and cash equivalents minus bank liabilities and loans

FY 2013 Key Figures

CHF m	2012	2013	Δ 12/13
Sales	391.1	388.3	-0.7%
Sales in local currencies	387.8	388.3	+0.1%
Gross Profit	198.4	189.6	-4.5%
in % of sales	50.7%	48.8%	
R&D	51.4	45.3	-11.8%
in % of sales	13.1%	11.7%	
OPEX	146.7	136.7	-6.8%
in % of sales	37.4%	35.2%	
Operating profit / EBIT	52.7	54.8	+4.0%
in % of sales	13.5%	14.1%	
EBIT at 2012 constant currencies	52.7	56.6	+7.4%
In % of sales	13.5%	14.5%	
Net profit	42.4	45.7	+7.8%
in % of sales	10.8%	11.8%	
EPS	3.92	4.16	+6.1%
Return on net assets (RONA)	42%	32%	-23.8%
Net Added Value Index ⁽¹⁾	1.38	1.38	0.0%
Net liquidity December 31 ⁽²⁾	141.3	143.4	+1.5%
Working capital ⁽³⁾	122.9	159.5	+29.8%
Equity	293.6	336.2	+11.0%
Cash Flow (operating)	2.4	27.9	+1062.5%

(1) Net Added Value Index (NAVI) = (EBIT + personnel expenses)/personnel expenses

(2) Net Liquidity = cash and cash equivalents minus bank liabilities and loans

(3) Includes customer-specific development costs capitalized of CHF 110.4m as of Dec, 31, 2013 (Dec, 31, 2012: CHF 73.7m)

Agenda

- Introduction Dr. David Martyr
- Financial Results Dr. Rudolf Eugster
- **Priorities** Dr. David Martyr
- Outlook & Summary Dr. David Martyr
- Questions & Answers

Review of Priorities 2013

	Commence instrument shipments for Dako Omnis (P16)	<ul style="list-style-type: none"> • Development successfully completed • Instrument launched in September; very positive end-user feedback
	Commence instrument shipments for Ortho Vision (P14)	<ul style="list-style-type: none"> • Large batch of instruments for validation was supplied to partner in October • Preparations for series production and market launch started
	Increase sales momentum in China	<ul style="list-style-type: none"> • Sales again increased with a clear double-digit growth rate • Sales exceeded CHF 25m (2012: >20m)
	Improving the innovation process	<ul style="list-style-type: none"> • Ongoing changes to processes and adjustments within R&D organization • On track to launch next generation liquid handling platform in 2014
	Continue to drive components business	<ul style="list-style-type: none"> • Posted solid sales growth; refined strategy • Strong growth in Next Gen Sequencing and increased footprint in China

Priorities for 2014

Focus on driving growth in Life Sciences Business, especially in Europe and North America

Supporting Partnering Business customers with delivery ramp-up (Dako Omnis and Ortho Vision)

Continue to focus on building out China (Life Sciences Business and Partnering Business)

Improving operational excellence and reducing manufacturing costs (COGS)

Driving appropriate, thoughtful M&A as catalyst to core business and to support evolution into solutions

Driving Growth in the Life Sciences Business

- Improved economic and market environment in Europe and North America
- Recent and upcoming product introductions
 - Several new products with focus on improved user-friendliness launched early 2014
 - High-throughput ELISA workstation
 - AC Extraction plate: fast, convenient, pipette and shake sample preparation for Mass Spectrometry
 - Launch of next generation liquid handling platform focused on specific applications starting in 2014
- Strengthening the organization
 - Key regional hirings with extensive industry-specific experience
 - Expanding market coverage by rebalancing ratio of back-office / field functions



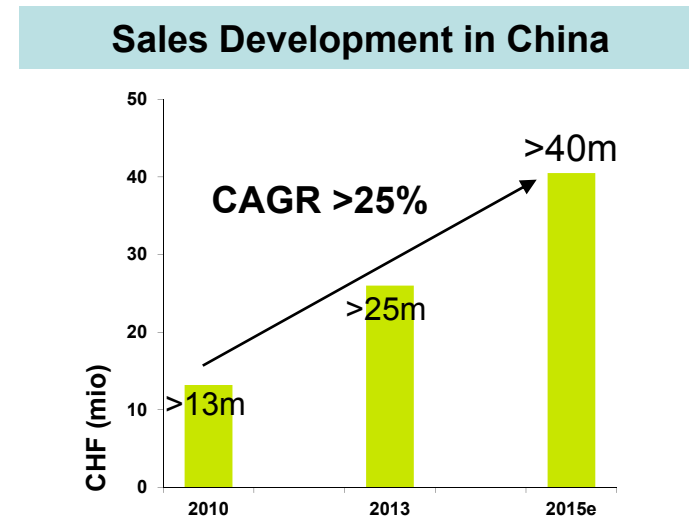
Partnering Business Delivery Ramp-up

- Partner Dako (an Agilent Technologies Company) successfully launched the Dako Omnis in 09/2013
 - Considerably higher production volumes of instrument expected in 2014 compared to 2013
- ORTHO Vision™ instrument will be launched by partner Ortho Clinical Diagnostics into several regional markets during 2014
 - Continued development of platform extension ORTHO Vision™ Max
- Tecan to ramp up serial production in support of phased regional product launches
- Close interaction between partners' support & complaint handling and Tecan expertline
- Continued close cooperation between R&D teams



Continue to Focus on Building out China

- Sales expected to continue to grow with a CAGR of around 25% to reach over CHF 40m by 2015
- Strengthened leadership and continuing to add “feet on the street”
- Expanding regional market coverage, dealer channels and service capabilities
- Added focus on Partnering Business sales
- High-throughput ELISA workstation, well suited to Chinese market



Improving Operational Excellence

- A multi-year project to reduce manufacturing costs (COGS) was launched at the start of 2014
- The biggest potential is lowering material costs
- Improved supply chain management and sourcing are key factors to develop profitability further
- Increased sourcing from around the world through supplier relocation and consolidation
- Freight cost optimization
- Additional longer-term benefits to be realized with the start of new product development programs
 - Building modular, flexible platforms
 - Increased re-use of common modules



Driving Appropriate, Thoughtful M&A

- M&A is a key element of the corporate strategy
- Team and approach
 - New experienced Corporate Development team with range of skills (scientific background, deep industry knowledge, finance expertise, project management)
 - Corporate strategy to define areas of focus
 - Active identification of targets, build-up of target funnel and cultivation of targets
 - Disciplined approach on valuation
 - Early involvement of key functions to ensure successful integration
- Areas of focus
 - Catalyst to core business and to support evolution into solutions
 - Ideally high consumables/reagents content, focusing on niches where Tecan is advantaged owner
 - Also interested in bolt-ons to expand addressable market



Agenda

- Introduction Dr. David Martyr
- Financial Results Dr. Rudolf Eugster
- Priorities Dr. David Martyr
- Outlook & Summary Dr. David Martyr
- Questions & Answers

Outlook for 2014

Sales

At least mid single-digit sales growth in local currencies

Profitability

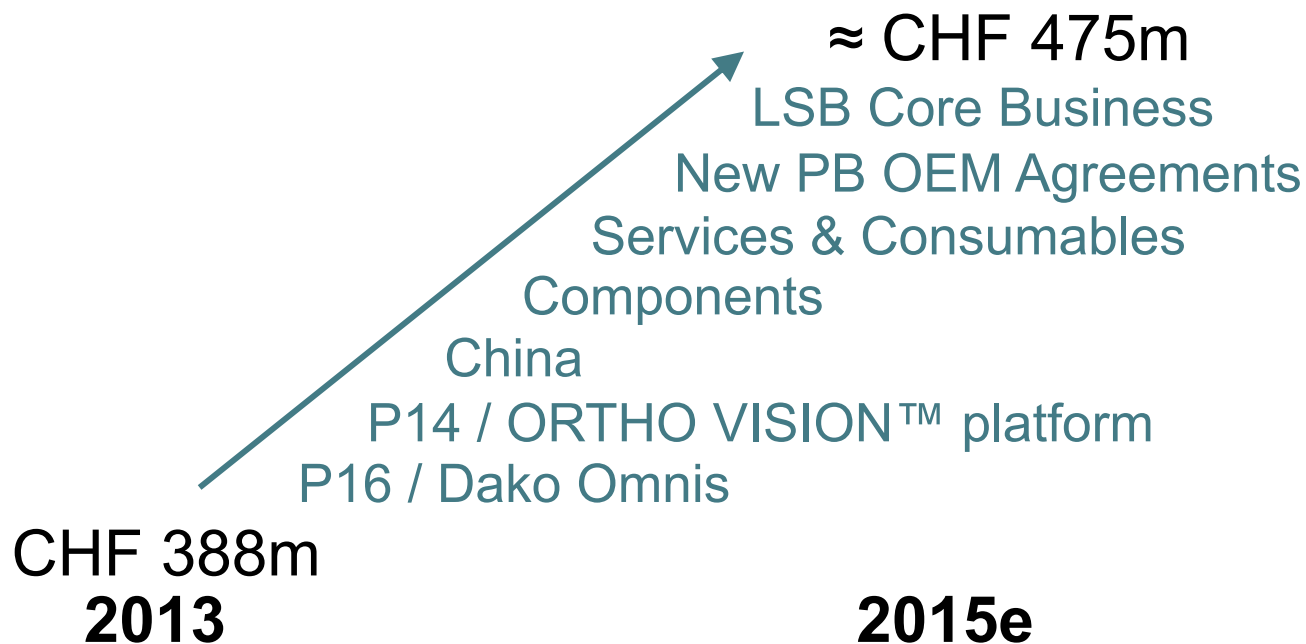
Further increase of EBIT Margin by around 50 bps
Based on average FX rates of: 1.21 EUR/CHF and 0.92 USD/CHF



Medium-Term Targets for 2015

Sales of around CHF 475 million at current currencies and increased profitability

- Updated objective reflecting negative FX impact since March 2013 (estimated CHF 15 million) and 2013 business results
- Actual rates in 2015 may differ and therefore can have a negative or positive effect
- Major growth drivers are intact, significant growth step expected in 2015 (P14)



Summary and Conclusion

- Financial objectives for 2013 only partly met
- Topline below original expectations, profitability target achieved
- Good progress in development programs
- Strengthened organization
- Refined strategy in place, including focus on M&A
- Major organic growth drivers for medium-term targets are intact
- Strong balance sheet and key financial ratios
- Unchanged dividend proposed



Agenda

- Introduction Dr. David Martyr
- Financial Results Dr. Rudolf Eugster
- Priorities Dr. David Martyr
- Outlook & Summary Dr. David Martyr
- Questions & Answers

Next Events

2014:

March 25: Publication of full Annual Report 2013

April 14: Annual Shareholder Meeting

August 13: Half Year Results 2014

Contact

Martin Braendle

Head of Corporate Communications & Investor Relations

Phone: +41 (0) 44 922 84 30

investor@tecan.com

www.tecan.com



Safe Harbor Statement



All statements in this presentation not referring to historical facts are predictions of the future and constitute no guarantee whatsoever of future performances. They are subject to risks and uncertainties including, but not limited to, future global economic conditions, exchange rates, legal regulations, market conditions, activities of competitors and other factors outside the Company's control.

Tecan Group AG is making all efforts to include accurate and up-to-date information into this presentation. Yet, it cannot be ruled out that omissions or errors might have occurred. Therefore, Tecan Group AG cannot make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this presentation. Changes in this presentation can be made any time without notice. All mentioned trademarks are protected by law.

© Tecan Group AG, Tecan Trading AG, Switzerland, all rights reserved.