

SARTORIUS

Simplifying Progress

Batch Process Modeling – Step-By-Step Guide

Lennart Eriksson, Ph.D., Assoc. Prof.
Senior Lecturer and Principal Data Scientist



Born in Data Analytics



- Company founded in 1987 by Professor Svante Wold, in Umeå, Sweden
 - Originator of Chemometrics and the SIMCA® Methodology
- Patented technologies in Design of Experiments and Multivariate Data Analysis
- We help our customers bring high-quality products to market faster
- Part of Sartorius Stedim Biotech since April 2017
- Products like MODDE®, SIMCA® and SIMCA®-online
- Global strength with local presence

Business Growth Through the Entire Product Lifecycle

Active Dashboard
Interactive performance
insight

Control Advisor
Avoid problems before
they arise

MODDE®
Get it right from the start

SIMCA®
Turn data into growth

SIMCA®-online
Ensured manufacturing success

Umetrics® Suite

Education/training and consulting
Transferring expertise to you

Upcoming Webinars

(<https://www.sartorius.com/en/company/exhibition-conferences>)

The screenshot displays a web interface for Sartorius webinars. At the top, there are navigation tabs: 'UPCOMING EVENTS' (selected), 'PAST EVENTS', 'CALENDAR' (with a calendar icon), and 'SCHEDULE' (with a plus icon). Below the tabs is a search bar with a magnifying glass icon and the text 'Search'. The main content area is divided into two columns of event cards. Each card features a green square icon with a white checkmark, a 'STANDARD' tag, the event title, and the date and time. A vertical ellipsis menu is located to the right of each card. The events listed are:

Event Title	Date	Time
Design of Experiments (DOE) for the Beginner	TUE, JAN 26, 2021	03:00 PM - 04:00 PM CET
Multivariate Data Analysis (MVDA) for the Beginner	THU, JAN 28, 2021	03:00 PM - 04:00 PM CET
Lean-and-clean DOE using One-click analysis	TUE, FEB 16, 2021	03:00 PM - 04:00 PM CET
OPLS® in process modeling	THU, FEB 18, 2021	03:00 PM - 04:00 PM CET
Robust optimization made easy	TUE, MAR 2, 2021	03:00 PM - 04:00 PM CET
Analyzing batch process data, a step-by-step guide	THU, MAR 4, 2021	03:00 PM - 04:00 PM CET
From Design of Experiments to Design Space Estimation	TUE, MAR 23, 2021	03:00 PM - 04:00 PM CET
Multiblock Orthogonal Component Analysis (MOCA) - A Novel Tool for Data Integration	THU, MAR 25, 2021	03:00 PM - 04:00 PM CET

Contents

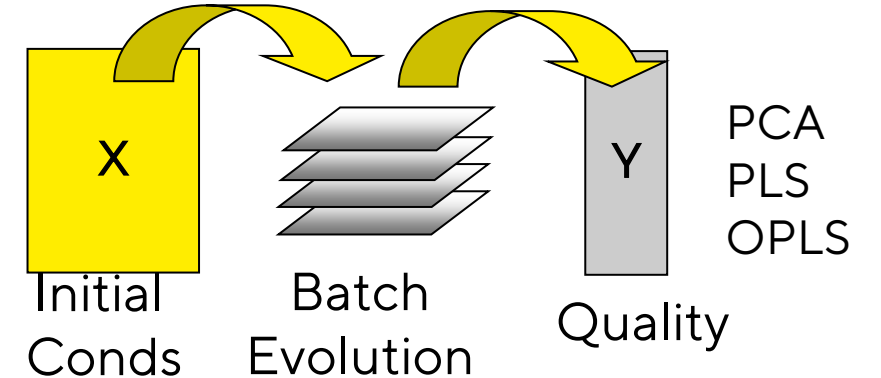
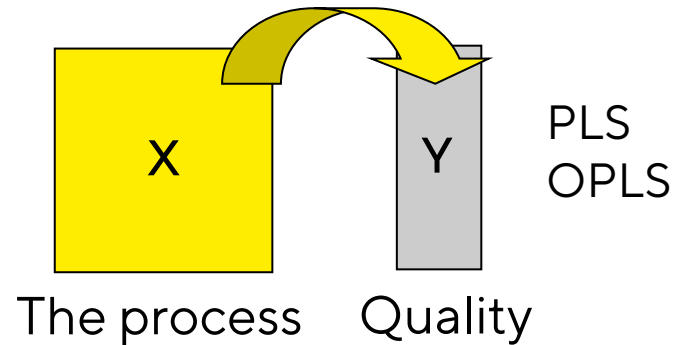
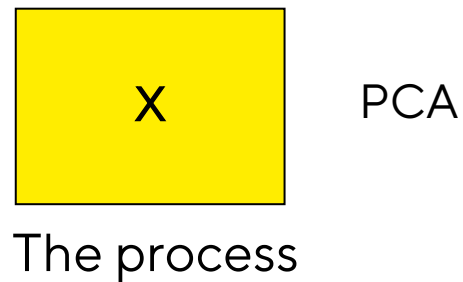
- Multivariate Process data
 - Different blocks of data and methods for their analysis
 - Statistical Process Control (SPC); SPC extensions: MSPC & BSPC
- Data lay-out and data sources
 - Define process and data sources
 - Database import possibilities
 - Process measurements and batch conditions
 - Two model perspectives, BEM & BLM
- Modeling and monitoring a fermentation process producing baker's yeast
 - Batch control charts
 - Early fault detection and interpretation of deviations
 - Model validation
- Modeling a cell culture process
 - Using the BEM to establish NOC
 - Using the BLM to predict final titer
- Demo & Summary

Multivariate Process Data

- Monitoring a process
 - Early warning of disturbances
 - Diagnostics - finding "assignable causes"

- Modelling a process output
 - Monitor Quality of final product

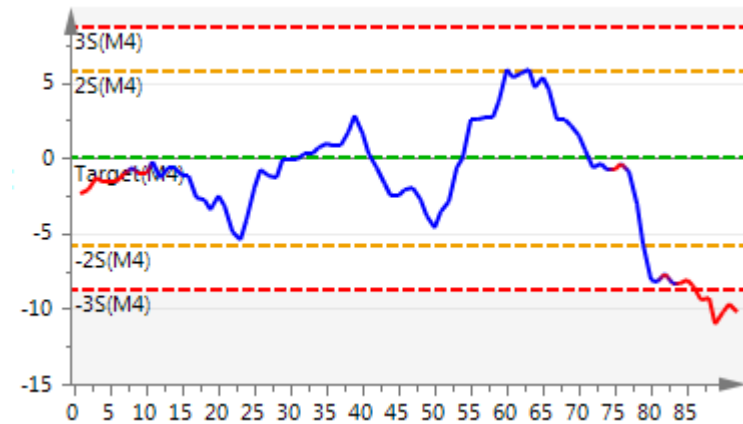
- Modelling Batch Processes
 - Majority of industrial processes
 - More complex analysis



Statistical Process Control, SPC

Walter A. Shewhart ~1930

- Assumption
 - A process is in a state of “statistical control” unless a special event occurs
- Approach
 - Devise a test to detect the occurrence of any special event
- Response
 - Look for “assignable cause” for the special event. Correct the process back to target
- Result
 - Robust process performance, long term process improvement

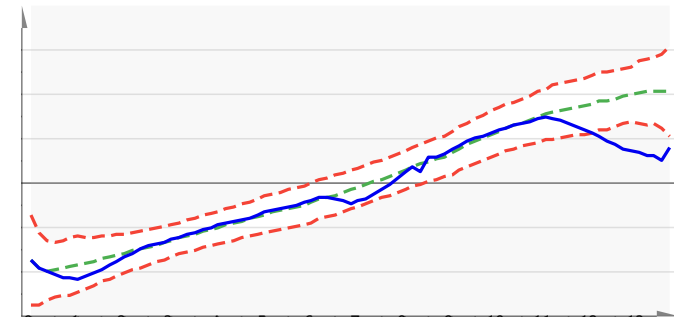


Using data analytics, the transition from SPC to multivariate SPC (MSPC) is obvious

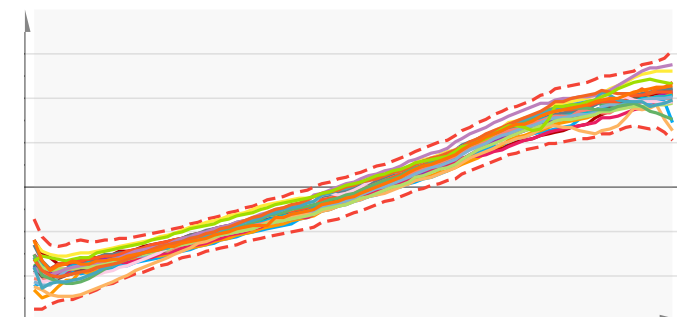
Batch statistical process control, BSPC

- The same philosophy is the basis for Batch Multivariate SPC
- For applications in Batch processes the evolution of the batch must be taken into account
- A process “road” for normal evolution is constructed
 - A reliable monitoring model should be able to detect when the process does not evolve in a normal way
- Any deviations are considered as abnormal process events
 - To be analyzed and acted upon

BakersYeast.M3, PS-Complement Batches, Model 3
Predicted Scores [Comp. 1]



BakersYeast.M3
Scores [Comp. 1]



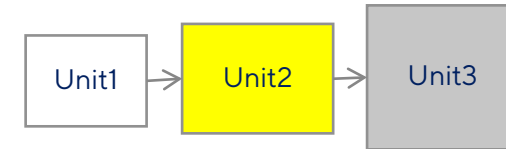
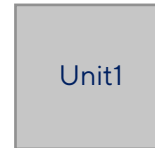


Data Lay-Out and Data Sources

CONNECTION
ANALYSIS
DATA
SEARCHING
VERIFICATION

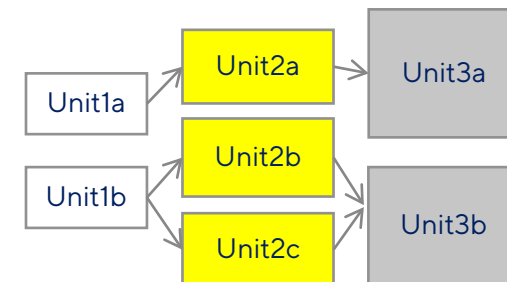
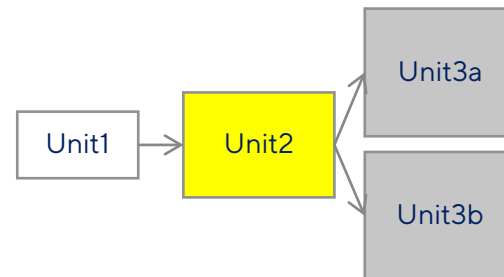
Define Process

- What does the process look like?
 - How many phases?
 - Process steps/ units
 - Split/ Merge batches



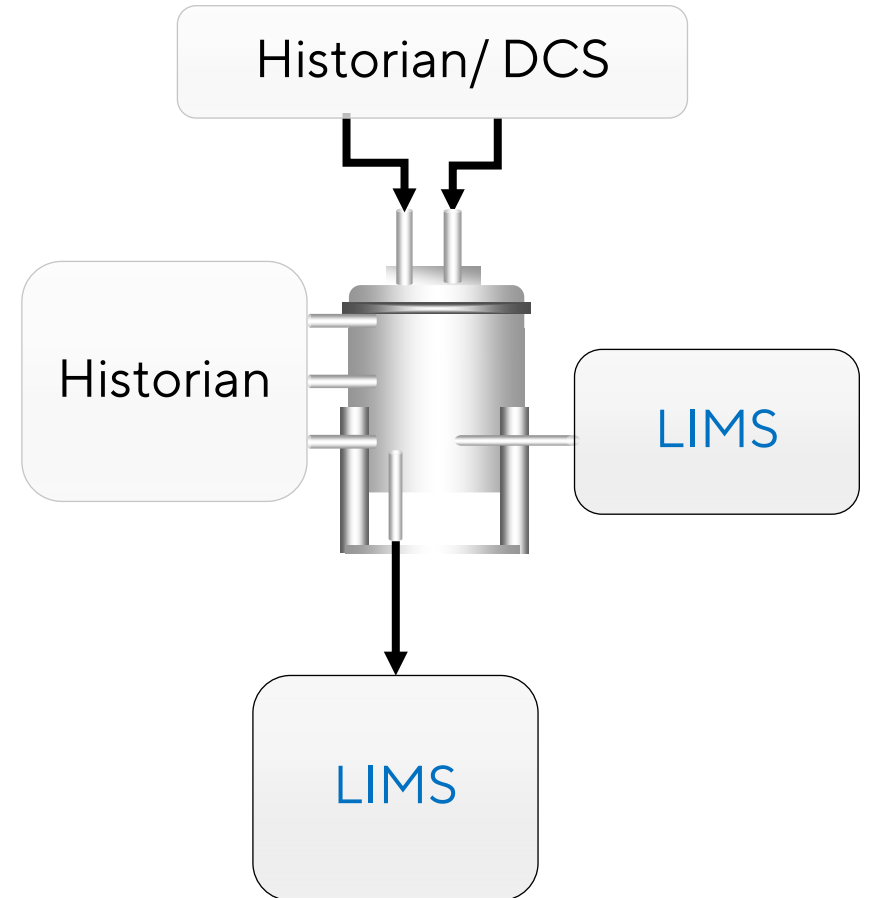
- Which parts of the process should be modeled?

- Which data is relevant?
 - Process tags
 - Quality information



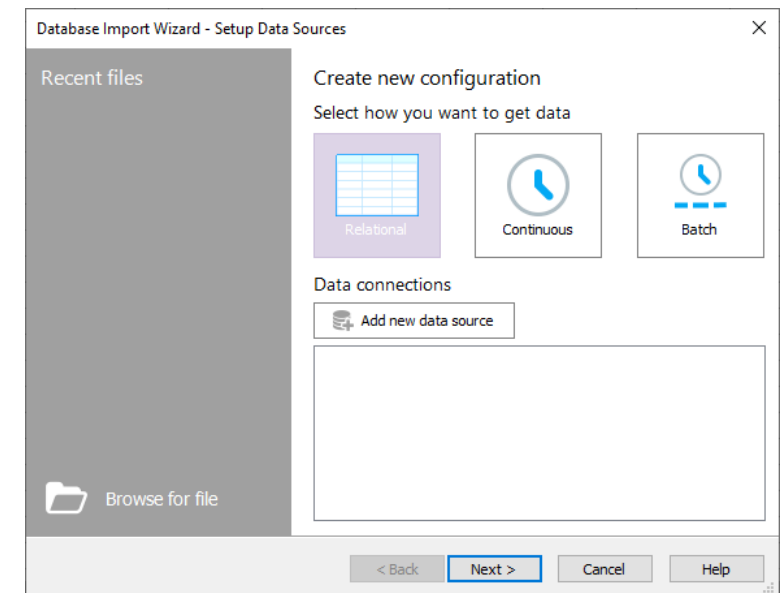
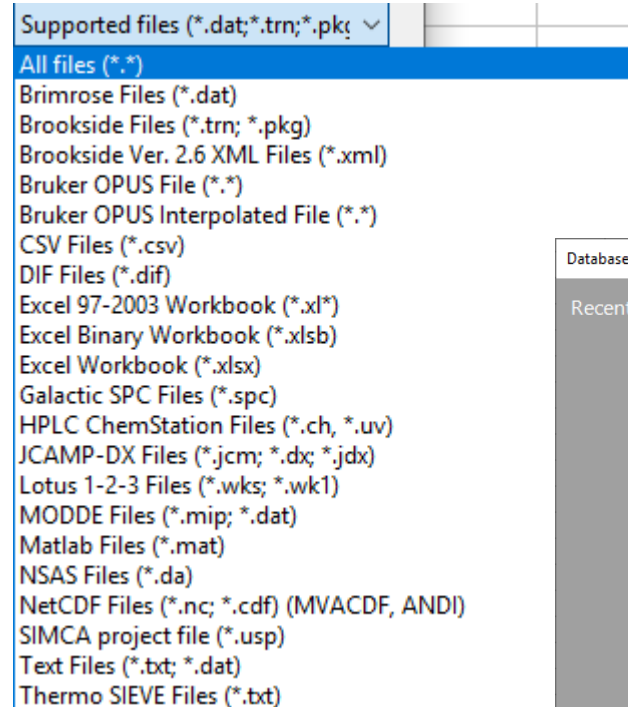
Available Data Sources

- Different types of data
 - Start-point information
 - Set-points pH, temp
 - Process parameters
 - Controlled parameters
 - Controlling parameters (actuators)
 - IPC / At-line data (ex. daily measurements)
 - Quality attributes
- Data stored in different places
 - LIMS, Historian...
- Depending on objective different data is used
- For analysis they need to be correctly synchronized



Process Database Data Extraction

- Export from DB to SIMCA supported file format
 - Most databases support data extraction to Excel
 - Data must be organized correctly
 - Data may be distributed over several files or excel sheets
- SIMCA can connect directly to a process data base
 - SIMCA import makes sure data is presented as required by SIMCA
 - Requires SimApi compatible with DB



SimApi

	Aspe n IP21	CSV	Disc over ant	June 5	MFC S	ODB C	OPC HDA	OPC UA	OSIs oftPI	SIPA T	SynT Q	Won derw are
Current	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>
Historical	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>
Discrete			<input type="radio"/>			<input type="radio"/>						
Batch	<input type="radio"/>		<input type="radio"/>		<input type="radio"/>	<input type="radio"/>			<input type="radio"/>	<input type="radio"/>		
Write back	<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>
Write back discrete						<input type="radio"/>						
Write back batch						<input type="radio"/>			<input type="radio"/>			
Node hierarchy	<input type="radio"/>						<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			
Array tag expansion	<input type="radio"/>						<input type="radio"/>					
Multiple data sources	<input type="radio"/>			<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>			
Connection resiliency		<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>			
3 rd party developed				<input type="radio"/>	<input type="radio"/>			<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	

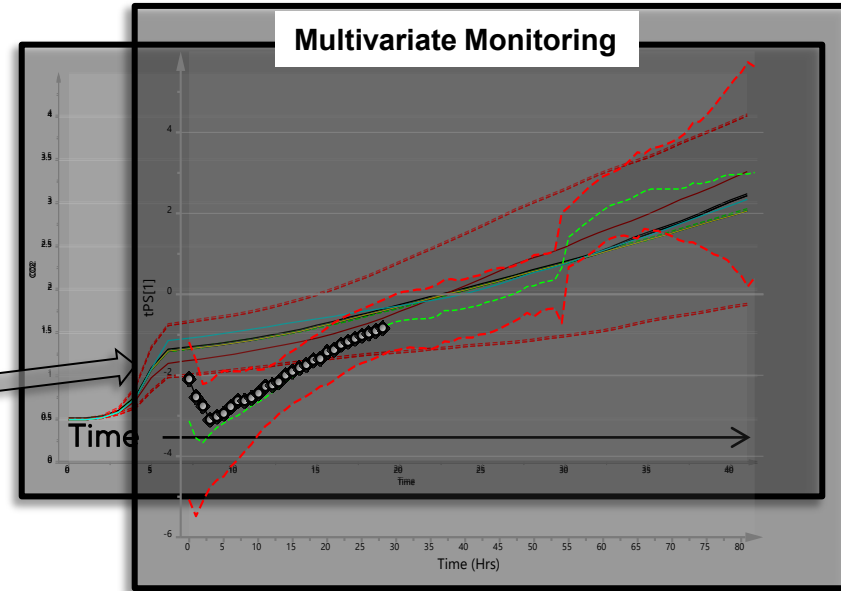
Batch Data

Process Data

Time ↓

Variables →

BatchID	TimeHrs	CL	V	CO2	T	F	Fc	Fb	fg	Ssp	Tsp	pH	
1	0	1.16	100	0.5	297	0	10	0	2	0.02	320	5.1	
1	2	0	1.16	100	0.5	297	0	10	0	0.02	320	5.1	
1	3	0	1.16	100	0.5	297	0	10	0	0.026	330	5.1	
1	4	0	1.16	100	0.5	297	0	10	0	0.02	320	5.1	
1	5	0.08333	1.15695	99.995	0.50013	296.998	0	10	0	2	0.02	310	5.1
1	10	0.16667	1.15172	99.837	0.52552	296.954	0	2	0.076967	1.86708	0.02	310	4.99307
1	20	0.33333	1.13769	100.56	0.59000	297.024	0	2	0.201543	1.75939	0.02	310	4.98243
1	30	0.5	1.09362	103.30	0.74793	297.381	0	2	0.506962	1.71858	0.02	310	4.95818
1	40	0.83333	1.03992	110.61	1.08958	298.377	0	2	1.13705	1.81531	0.02	310	4.92515
1	50	1.16667	1.11777	119.21	1.41539	300.353	0.032947	2	0.002955	1.88646	0.02	310	5.12493
1	60	1.5	1.11907	118.99	1.43866	302.636	0.032809	2	0.047028	1.83693	0.02	310	4.99949
1	70	1.83333	1.11749	118.71	1.4618	304.982	0.033705	2	5.00E-02	1.79198	0.02	310	4.9997
1	80	2.16667	1.11582	118.31	1.48773	307.398	0.034648	2	0.05304	1.75115	0.02	310	4.99969
1	90	2.5	1.11405	117.75	1.51579	309.894	0.035639	4.19453	0.056201	1.71465	0.02	310	4.99968
1	100	2.83333	1.11224	117.09	1.54988	311.286	0.036476	554.565	0.057968	1.68254	0.02	310	4.99996
1	110	3.16667	1.11088	116.31	1.57986	310.745	0.036956	1203.54	0.056805	1.65427	0.02	310	5.00019
1	120	3.5	1.10958	115.67	1.60909	309.532	0.037288	1213.61	0.054588	1.62994	0.02	310	5.00021
1	130	3.83333	1.10827	115.11	1.63833	309.03	0.037722	689.387	0.053342	1.60766	0.02	310	5.00005
1	140	4.16667	1.10683	114.54	1.66773	309.857	0.038383	347.719	0.053973	1.58946	0.02	310	4.99988
1	150	4.5	1.10537	113.91	1.69863	310.815	0.039089	646.424	0.054926	1.57495	0.02	310	4.99996
1	160	4.83333	1.10406	113.25	1.73001	310.684	0.039561	1143.03	0.054368	1.56384	0.02	310	5.00011
1	170	5.16667	1.10292	112.64	1.76105	309.889	0.040002	1278.96	0.052841	1.5557	0.02	310	5.00016
1	180	5.5	1.10182	112.10	1.79088	309.372	0.040425	992.11	0.051669	1.55031	0.02	310	5.00008
1	190	5.83333	1.10069	111.57	1.82053	309.696	0.040982	689.535	0.051634	1.54772	0.02	310	4.99996
1	200	6.16667	1.09953	111.01	1.85108	310.412	0.041612	769.918	0.052185	1.54803	0.02	310	4.99996
1	210	6.5	1.09954	110.14	1.89053	313.938	0.040985	2	0.039404	1.54806	0.014	325	4.99976
1	220	6.83333	1.09819	108.90	1.91548	317.738	0.041787	2	0.042336	1.55154	0.014	325	4.99972
1	230	7.16667	1.09675	107.23	1.93932	321.692	0.042335	2	0.04464	1.55856	0.014	325	4.99982
1	240	7.5	1.09532	105.01	2.01079	325.734	0.043048	100.68	0.044996	1.5692	0.014	325	5.0001
1	250	7.83333	1.09397	102.32	2.07223	327.348	0.043261	1293.94	0.04297	1.58325	0.014	325	5.00017
1	260	8.16667	1.09258	99.694	2.13524	326.148	0.043594	2373.11	0.042058	1.60076	0.014	325	5.00007
1	270	8.5	1.09142	97.459	2.19423	324.076	0.043885	2308.41	0.040861	1.62139	0.014	325	5.00015
1	280	8.83333	1.09044	95.529	2.24977	323.26	0.044152	1354.8	0.039957	1.64456	0.014	325	5.0001
1	290	9.16667	1.08944	93.600	2.30299	324.689	0.044466	725.846	0.038866	1.67018	0.014	325	5.00007
1	300	9.5	1.08843	91.438	2.37079	326.462	0.04474	1237.02	0.037657	1.69834	0.014	325	5.00015
1	310	9.83333	1.0874	89.191	2.43878	326.386	0.044984	2187.11	0.036467	1.72907	0.014	325	5.00009
1	320	10.16667	1.08647	87.145	2.50508	324.973	0.04525	2541.49	0.035535	1.76223	0.014	325	5.0001
1	330	10.5	1.0857	85.356	2.56777	323.883	0.045487	2077.8	0.03442	1.79742	0.014	325	5.00011
1	340	10.83333	1.08498	83.663	2.63099	324.312	0.045751	1478.67	0.0336	1.83442	0.014	325	5.00007
1	350	11.16667	1.08423	81.975	2.69393	325.672	0.046001	1551.6	0.032784	1.87246	0.014	325	5.0001
1	360	11.5	1.08346	79.985	2.75893	326.198	0.046213	2214.82	0.031715	1.91403	0.014	325	5.0001
1	370	11.83333	1.08272	78.179	2.81448	325.446	0.046439	2711.22	0.030852	1.95669	0.014	325	5.00008
1	380	12.16667	1.08209	76.549	2.87124	324.402	0.046653	2598.47	0.029954	2.00101	0.014	325	5.00009
1	390	12.5	1.08154	75.039	2.92752	324.351	0.046873	2170.21	0.02915	2.04674	0.014	325	5.00007
1	400	12.83333	1.08107	73.515	3.00058	325.208	0.047093	2062.9	0.028439	2.09389	0.014	325	5.00008

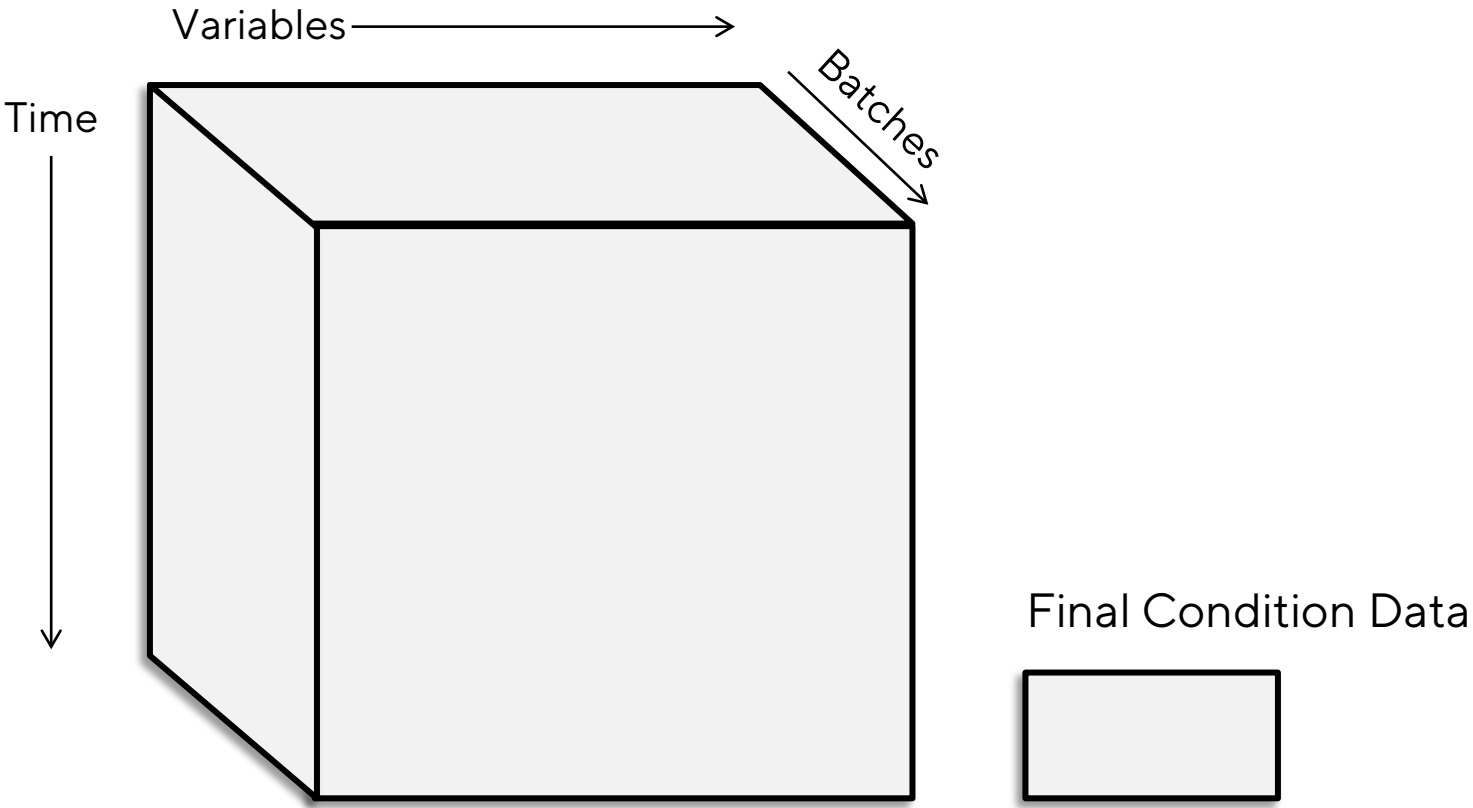


Final Condition Data

Batch ID	[Pen]	Amount	Volume
1	2.305952	164.535	71.3523
2	2.140447	164.238	76.7307
3	2.397035	146.426	61.0863
4	0.467192	76.7307	164.238
5	0.418107	73.515	175.828

Batch Data

- The challenge is to model the relationship between process variation and quality (final conditions)



Process Data and Batch Conditions

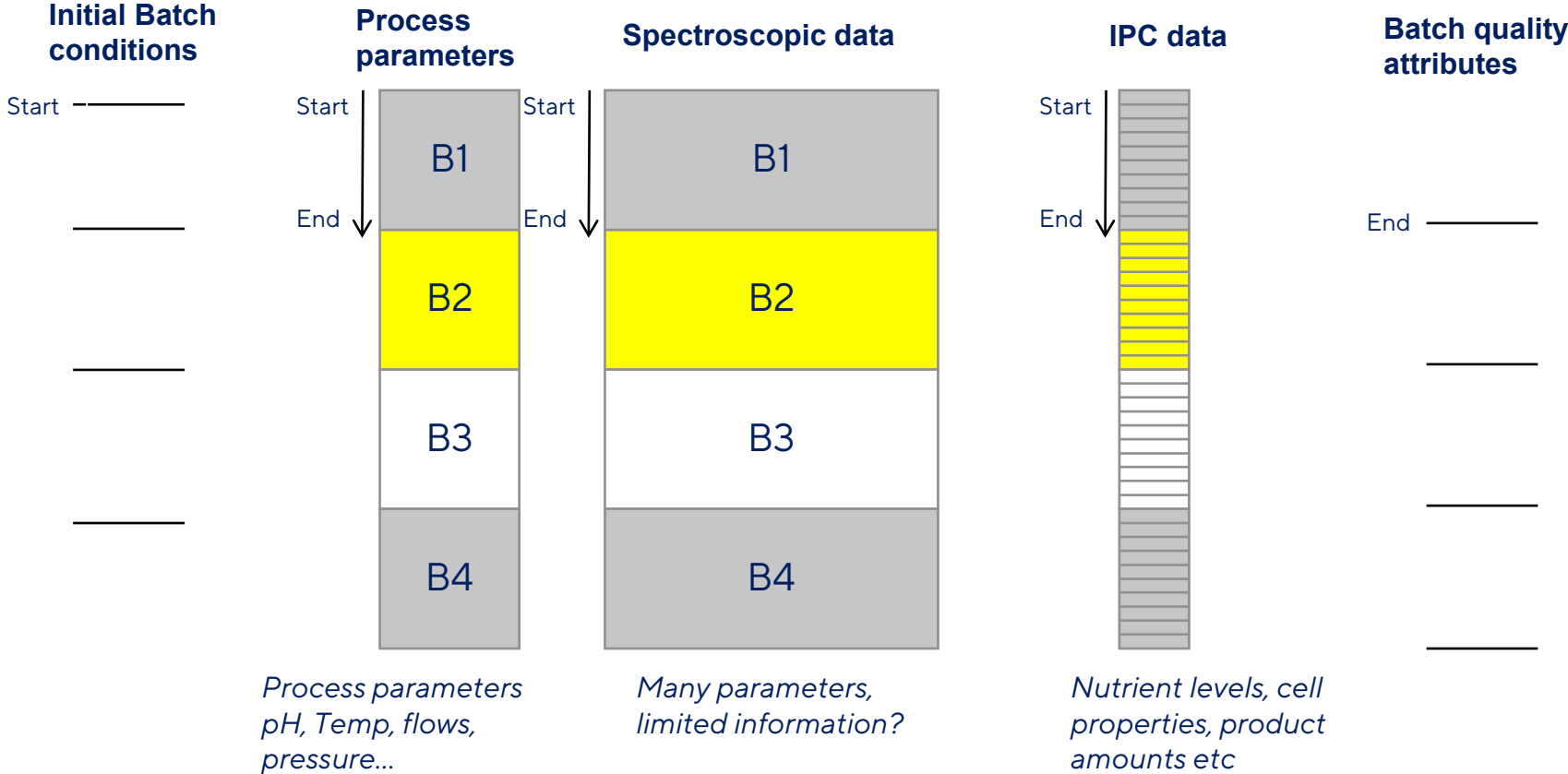
- Process data describes the process evolution
 - Timely measurements
 - All parameters are extracted at the same time frequency

- Batch conditions does not have an evolution
 - The same value true for the entire batch
 - Starting conditions, intermediate quality, final batch properties

	2	3	4	5	6	7	8	9	10
1	SBatchID	Batch Age	Ethanol	Temp	Molasses	NH3	Air	Level	pH
153	Ba	11,1667	0,0577	30,28	3252,92	98,5	6704	54,00	5,379
154	Ba	11,3333	0,0528	30,28	3263,22	98,5	6693	54,24	5,380
155	Ba	11,5	0,0497	30,31	3268,86	98,3	6718	54,53	5,376
156	Ba	11,6667	0,0469	30,35	3273,44	98,2	6704	54,89	5,369
157	Ba	11,8333	0,0444	30,38	3279,19	98,0	6689	55,13	5,363
158	Ba	12	0,0419	30,32	3279,62	10,0	6700	55,47	5,376
159	Ba	12,1667	0,035	30,29	3270,92	0,10	6700	55,72	5,445
160	Ba	12,3333	0,0303	30,23	3250,12	0,12	6709	55,91	5,541
161	Ba	12,5	0,0270	30,21	3223,85	0,10	6695	56,21	5,650
162	Ba	12,6667	0,0244	30,21	3202,48	0,11	6701	56,45	5,764
163	Ba	12,8333	0,0223	30,25	3174,19	0,10	6697	56,72	5,880
164	Ba	13	0,0203	30,28	3122,73	0,12	6702	57,06	5,994
165	Ba	13,1667	0,0184	30,28	2679,03	0,11	6692	57,25	6,123
166	Ba	13,3333	0,0165	30,14	2014,34	0,09	6701	57,41	6,277
167	Ba	13,5	0,0148	30,06	1389,3	0,10	6706	57,70	6,447
168	Ba	13,6667	0,0134	29,90	859,207	0,12	6718	57,58	6,591
169	Ca	0	0,0662	29,82	667,187	47,4	2306	41,23	3,913
170	Ca	0,166667	0,1449	29,69	997,483	73,4	2496	41,30	4,024
171	Ca	0,333333	0,3917	29,61	1165,18	86,2	2672	41,39	4,466
172	Ca	0,5	0,7348	29,55	1087,61	80,4	2857	41,48	4,806
173	Ca	0,666667	1,0436	29,46	995,638	73,8	3033	41,56	5,001
174	Ca	0,833333	1,2598	29,52	929,693	68,3	3204	41,63	5,075
175	Ca	1	1,3986	29,53	892,962	66,2	3366	41,71	4,979
176	Ca	1,16667	1,5052	29,44	871,689	63,8	3533	41,78	5,078
177	Ca	1,33333	1,5369	29,53	883,85	65,6	3700	41,85	4,989
178	Ca	1,5	1,5637	29,55	905,478	65,8	3870	41,93	5,026
179	Ca	1,66667	1,5785	29,42	942,776	72,3	4029	42,00	5,072
180	Ca	1,83333	1,5812	29,49	974,933	74,9	4195	42,08	5,080
181	Ca	2	1,5930	29,57	1006,21	78,5	4362	42,17	5,035
182	Ca	2,16667	1,5873	29,51	1051,41	82,8	4527	42,24	5,091
183	Ca	2,33333	1,5810	29,50	1088,24	84,3	4690	42,34	5,066
184	Ca	2,5	1,5668	29,45	1141	88,9	4857	42,42	5,08
185	Ca	2,66667	1,5518	29,45	1192,38	93,0	5026	42,53	5,091
186	Ca	2,83333	1,5296	29,50	1263,22	99,2	5195	42,63	5,060
187	Ca	3	1,5114	29,55	1324,52	103,	5359	42,73	5,109

	2	3	4	5	6	7
1	SBatchID	Innoc	QP1	QP2	Amount	Yield
2	bb	914	93	77	5365	0,482205
3	Ga	932	86	77	6089	0,48
4	Ha	940	81	71	5904	0,46
5	hb	943	88	82	5977	0,470986
6	gb	950	89	78	5982	0,47437
7	cb	952	91	83	5875	0,500618
8	db	952	91	82	5835	0,505364
9	eb	952	90	80	5973	0,4816
10	Za	952,32	86	77	5541	0,45427
11	fb	960	81	72	5402	0,46254
12	Va	964	70	63	4689	0,45
13	ab	967	85	81	5749	0,492269
14	ib	973	91	80	5977	0,497475
15	jb	973	91	79	6147	0,480445
16	kb	980	88	82	6154	0,485785
17	lb	981	93	83	5919	0,481485
18	Xa	993,28	84	68	6597	0,504218
19	Ua	1027	94	91	6427	0,46
20	Aa	1049	92	93	5715	0,42
21	Ta	1049,04			6960	0,505668
22	Ma	1050,56	89	82	6658	0,479405
23	Pa	1050,56	89			0,556606
24	La	1051	83	76	6515	0,55
25	Oa	1051	94	89	4442	0,44

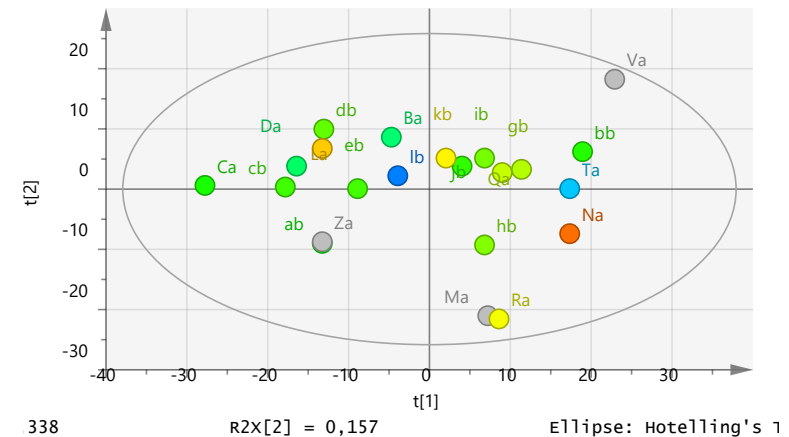
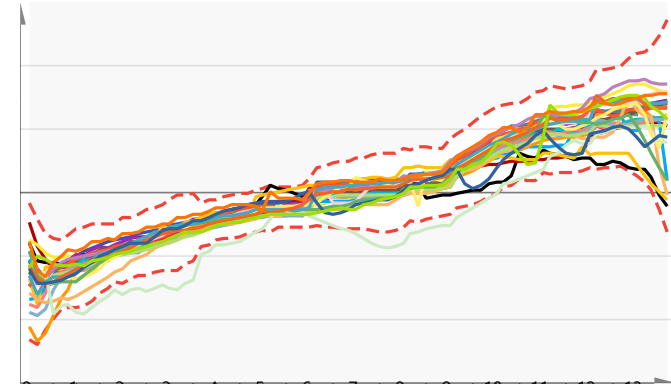
Data Overview



BSPC - From Two Perspectives

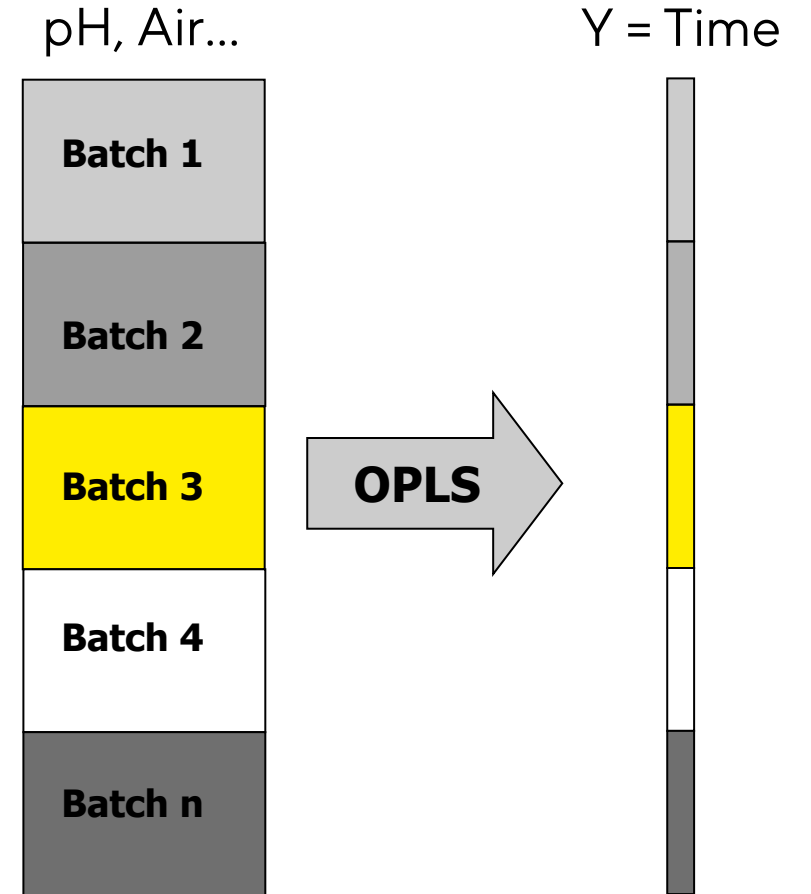
- Batch processes evolve over time
 - Many measurements over time
 - Batch Evolution Model (BEM)

- Quality, yield etc are summarized per batch
 - One measurement per batch
 - Batch Level Model (BLM)



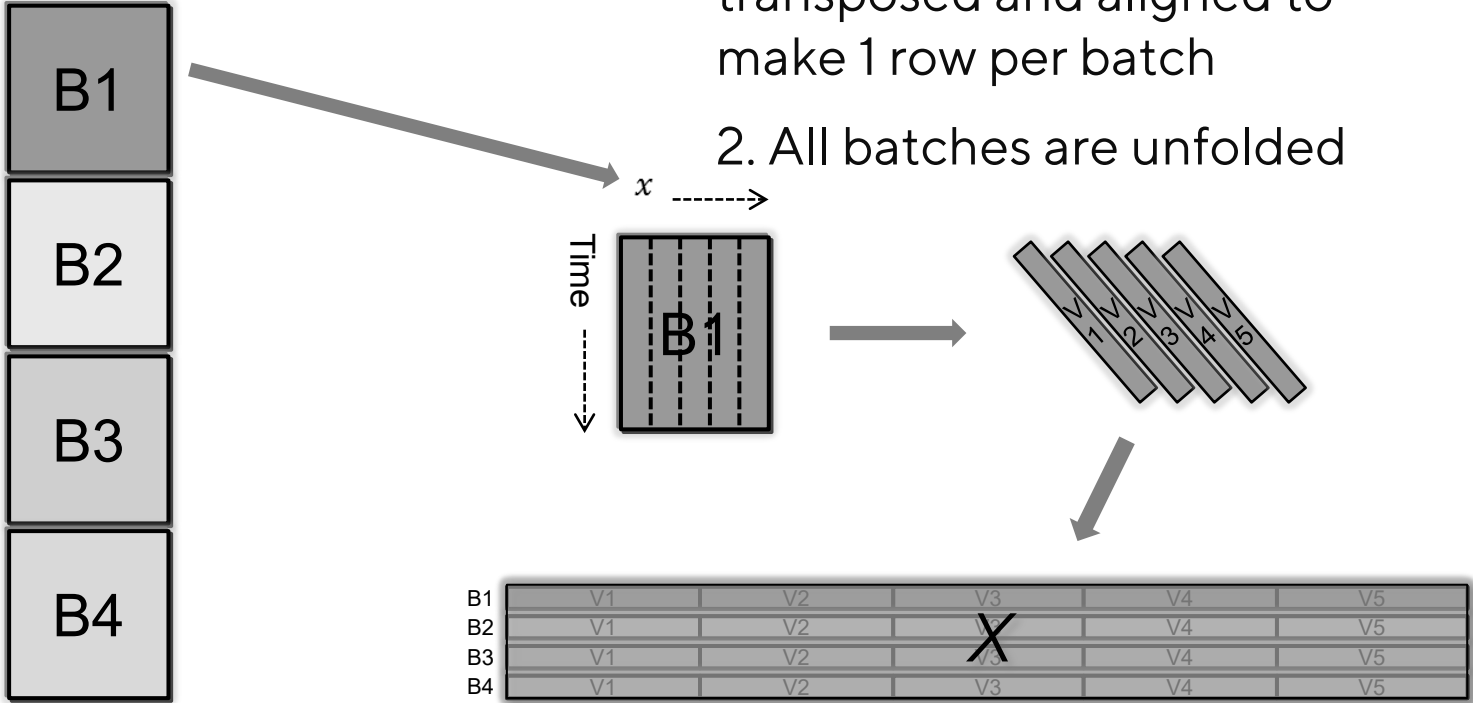
The Batch Evolution Model (BEM)

- Time (or maturity) is used as a Y-variable to give the model a direction
- Maturity need not be time. It could be say, for example, be Ethanol in beer brewing



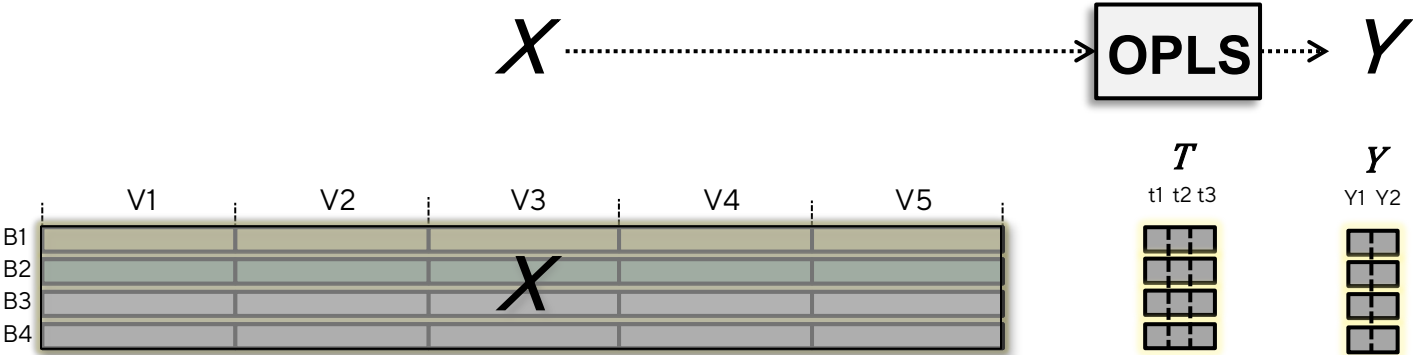
Batch Data; BEM and BLM

- Batch data has time dependency
 - A table of data is generated for each batch
 - Variables measured over time



The Batch Level Model (BLM)

- OPLS can handle the highly correlated batch unfolded dataset



OPLS is used to map the correlation of the process trajectory (X) to final batch conditions (Y)



Example: Baker's Yeast (BEM)

Modelling of a Batch Process

- Example: Baker's yeast production
 - Dataset from Jästbolaget AB in Sweden
- 33 batches represented by 7 process variables and 5 batch conditions
 - Batch conditions: X parameter Inoculum and Y parameters QP1, QP2, amount and yield
- Objective: Establish model for normal process evolution
 - Same approach as for a continuous process
 - Model is built on well performing and behaving batches

Data Structure

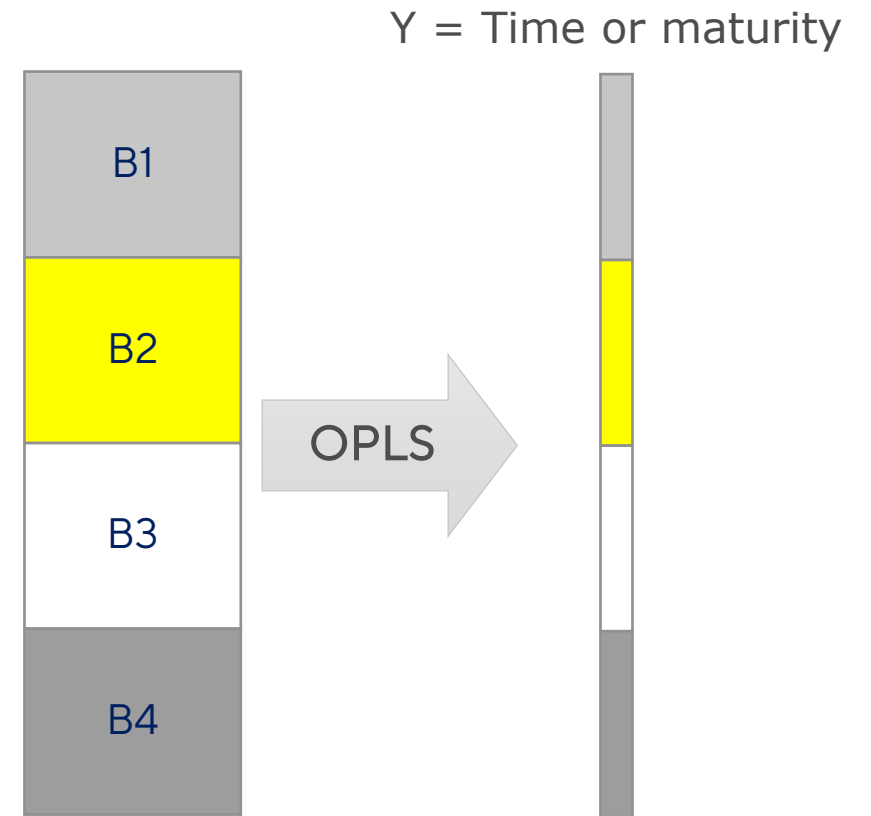
- Data are usually divided into 2 parts
 - Process data (from process DB) and Batch condition data (from LIMS/ lab DB)
- Process/ Evolution data table
 - Each row is one time point
 - One column containing batch ID
 - One column containing time
 - One column for each measured variable
- Batch condition data table
 - One row for each batch
 - Includes both initial (start) conditions and final batch results

BatchID	Innoc	QP1	QP2	Amount	Yield
Ba					
Ca					
Ia	1126,44	82	74	6368	0,493462
Ma	1050,56	89	82	6658	0,479405

BatchID	Batch Age (h)	Ethano	Temp	Molasses	NH3	Air	Level	pH
	TAGID Age001	TAGID E001	TAGID T001	TAGID FM001	TAGID FN001	TAGID FA001	TAGID L001	TAGID P001
Ba	0	0,00748	35,745	476,196	23,7236	1642,39	41,739	4,74916
Ba	0,1	0,00589	35,6153	1088,33	79,4786	214,35	41,8553	4,72612
Ba	0,2	0,022471	31,4477	110,36	61,1111	2342,21	41,9245	5,03004
Ba	0,3	0,38293	30,3056	1103,91	81,0109	2516,66	42,0182	5,04351
Ba	0,4	0,52197	30,1153	1101,82	80,1929	2704,02	42,0332	5,11628
Ba	0,5	0,53908	30,3526	1125,21	80,3073	2878,89	42,194	5,10217
Ba	0,6	0,62089	30,3014	1125,25	82,2556	3059,62	42,2775	5,05596
Ba	0,7	0,62848	30,432	1223,36	83,2116	3250,84	42,3894	5,11921
Ba	0,8	0,6288	30,3547	1200,65	84,4421	3361,14	42,4392	5,06234
Ba	0,9	0,64116	30,3445	1221,71	85,291	3469,93	42,5393	5,0456
Ba	1	0,63781	30,3887	1335,22	102,406	3648,03	42,7231	5,06487
Ba	1,1	0,6239	30,4464	1454,89	117,635	3894,62	42,8332	5,11161
Ba	1,2	0,63635	30,4195	1505,33	119,847	4081,53	42,9684	5,07076
Ba	1,3	0,63682	30,4062	1561,77	124,496	4288,1	43,1022	5,0702
Ba	1,4	0,66693	30,3954	1591,76	121,286	4396,04	43,2432	5,136
Ba	1,5	0,61144	30,4209	1656,4	122,312	4561	43,386	5,11634
Ba	1,6	0,70631	30,4195	1691,32	134,849	4734,59	43,5261	5,13043
Ba	1,7	0,74649	30,3981	1700,39	136,08	4905,19	43,6769	5,08862
Ba	1,8	0,7561	30,3927	1752,02	133,969	5041,69	43,8232	5,14039
Ba	1,9	0,78523	30,3834	1788,94	143,72	5235,29	43,9784	5,09226
Ca	0	0,06621	23,8236	667,187	47,4239	2306,88	41,2399	3,91324
Ca	0,1	0,44434	23,8293	3974,65	72,453	2495,62	41,5049	4,62487
Ca	0,2	0,39177	23,6177	1165,16	86,2716	2672,74	41,9365	4,46631
Ca	0,3	0,73487	23,559	1087,61	80,4525	2857,77	41,4882	4,80647
Ca	0,4	1,04285	23,4681	988,638	73,6315	3033,6	41,5867	5,00197
Ca	0,5	1,2538	23,5914	68,3734	68,3734	3204,5	41,6309	5,07865
Ca	0,6	1,38865	23,5349	892,362	66,2413	3386,03	41,7156	4,97938
Ca	0,7	1,50922	23,4454	871,689	65,8538	3533,8	41,7866	5,07898
Ca	0,8	1,59636	23,524	893,85	65,6356	3700,41	41,9517	4,91969
Ca	0,9	1,56373	23,559	805,478	65,8137	3870,66	41,93	5,02699
Ca	1	1,57155	23,4253	942,716	72,3848	4029,19	42,0022	5,07226
Ca	1,1	1,58124	23,4309	874,833	74,3663	4181,43	42,018	5,00927
Ca	1,2	1,59309	23,5737	1006,21	78,5567	4362,48	42,114	5,05587
Ca	1,3	1,58734	23,5122	1051,41	82,8364	4527,13	42,2471	5,09192
Ca	1,4	1,59889	23,5082	1086,24	84,301	4690,71	42,3415	5,06636
Ca	1,5	1,56685	23,452	1141	86,9122	4857,94	42,4225	5,08
Ca	1,6	1,55185	23,452	1196,39	93,0529	5026,86	42,5321	5,09196
Ca	1,7	1,52985	23,5095	1235,22	99,2319	5185,84	42,6327	5,06988
Ca	1,8	1,51146	23,555	1242,52	103,239	5339,45	42,7396	5,10555
Ca	1,9	1,48927	23,5903	1334,05	109,377	5532,02	42,8474	5,11779
Ca	2	1,46971	23,6289	1458	115,043	5693,43	42,9747	5,09786
Ia	0	0,00755	32,7412	95,489	67,4109	1454,73	41,7391	5,25271
Ia	0,1	0,00927	32,244	646,314	141,272	2060,92	41,7195	5,28144
Ia	0,2	0,02917	31,2095	700,168	141,332	2246,25	41,8504	4,40458
Ia	0,3	0,0696	30,2919	750,195	147,08	2433,64	41,9145	5,09611
Ia	0,4	0,02148	23,8236	774,388	147,196	2610,63	41,984	5,86605
Ia	0,5	0,04886	23,3532	812,629	147,411	2801,66	42,0623	5,17497
Ia	0,6	0,08038	30,0381	863,725	146,967	3009,14	42,199	5,20184
Ia	0,7	0,0883	30,0468	932,47	166,019	3263,19	42,2805	5,21192
Ia	0,8	0,12259	30,0214	987,423	16,9872	3436,25	42,3598	4,40948
Ia	0,9	0,16033	23,3825	1046,43	54,3661	3602,73	42,4461	4,18415
Ia	1	0,18011	30,004	1092,97	79,9821	3772,63	42,5336	4,23087
Ia	1,1	0,19878	30,008	1143,9	79,7954	3938,46	42,6518	4,33524
Ia	1,2	0,19174	23,3939	1240,44	79,6046	4107,3	42,7597	4,3309
Ia	1,3	0,1039	30,0173	1374,9	79,4123	4276,67	42,8709	4,39864
Ia	1,4	0,1909	30,0254	1436,27	79,1488	4443,29	42,9795	4,43225
Ia	1,5	0,15617	18,9385	1517,36	78,7337	4616,03	43,1051	4,38884
Ia	1,6	0,17027	30,0635	1747,34	78,319	4786,82	43,2636	4,36235
Ia	1,7	0,20691	30,147	1856,78	87,7843	4958,15	43,3963	4,34843
Ia	1,8	0,20695	30,147	1941,75	91,6167	5141,44	43,5263	4,24185
Ia	1,9	0,22396	30,1742	1935,29	91,2748	5434,43	43,7355	4,34213
Ia	2	0,25643	30,1175	2040,33	120,347	5368,98	43,8976	4,37778
Ma	0	0,17042	31,1189	69,2015	0,11429	1463,82	41,477	4,60889
Ma	0,1	0,27795	28,4127	284,471	61,5214	1916,99	41,9499	5,04947
Ma	0,2	0,37221	31,6595	806,234	65,3679	2250,15	41,5207	3,75869
Ma	0,3	0,66193	31,1009	903,244	70,3214	2450,6	41,634	3,91439
Ma	0,4	0,83476	30,7486	900,763	64,9643	2645,95	41,6714	4,05918
Ma	0,5	1,02756	30,6104	845,771	68,2179	2840,12	41,7223	4,19384
Ma	0,6	1,11933	30,6614	881,461	69,0072	3044,12	41,7999	4,17463
Ma	0,7	1,17445	30,7898	901,657	72,3359	3140,25	41,909	4,22025
Ma	0,8	1,21193	30,8706	940,624	75,4286	3350,11	41,9143	4,30142
Ma	0,9	1,22247	30,8706	952,2	76,7679	3518,6	42,0396	4,27682
Ma	1	1,20914	30,859	938,195	80,9929	3662,91	42,0955	4,43667
Ma	1,1	1,18651	30,7768	1177,85	94,8357	3853	42,1953	4,5057
Ma	1,2	1,17126	30,7459	1090,76	89,15	4015,53	42,2967	4,57611
Ma	1,3	1,11846	30,7239	1003,28	102,393	4215,11	42,3136	4,62339
Ma	1,4	1,07127	30,7486	1235,47	111,119	4366,53	42,525	4,68668
Ma	1,5	1,03262	30,7553	1382,33	116,629	4482,46	42,6695	4,78641
Ma	1,6	0,93695	30,754	1425,5	122,825	4637,2	42,6911	4,87114
Ma	1,7	0,84588	30,7767	1553,31	134,543	4836,48	42,8604	4,97266
Ma	1,8	0,90702	30,7982	1586,11	136,654	5022,28	42,8647	5,06116
Ma	1,9	0,89667	30,8278	1674,8	143,8	5192,63	43,119	5,07939
Ma	2	0,89611	30,8498	1755,47	149,468	5356,97	43,2785	5,11682

Data Structure In the BEM

- OPLS between process time points and time gives model representing process evolution in data set
- When based on “good” or “normal” batches the model represents normal batch evolution

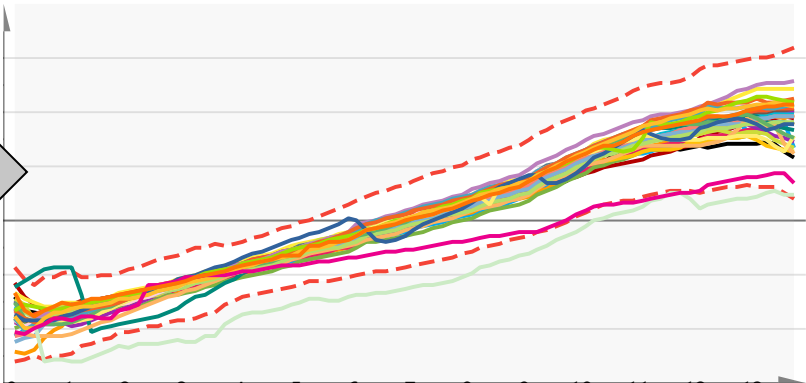


Observations, From Variables to Scores

- Observations are now described by "scores"
- Scores are linear combinations of original variables
 - Why are observations close to each other?
 - What is the reason between dissimilarities?

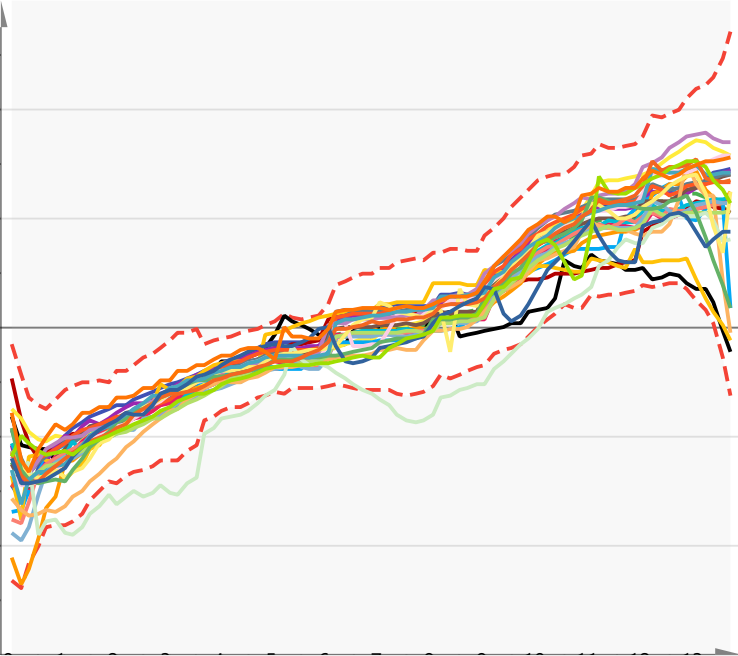
Batch	Observation	1	2	3	4	5	6	7	8	9	10	11	12	13	14
1	1	0.16667	0.53995	30.8995	832.241	4.81927	1849.75	42.4759	9.4295	832.541	4.81927	1849.75			
2	2	0.33333	0.27134	30.7574	688.828	8.35432	2171.85	42.7382	4.4249	1221.37	12.7726	4115.6	0.624008		
3	3	0.5	0.36416	30.6462	754.033	0.26431	2365.03	42.7909	4.67838	1975.4	18.0389	4445.43	0.110101		
4	4	0.66667	0.43617	30.6109	778.134	0.08213	2581.89	42.8518	4.87028	2792.34	18.1196	9026.39	0.178701		
5	5	0.83333	0.49184	30.2048	814.789	0.09287	2794	42.9077	4.88195	3870.33	13.2119	11748.6	0.231899		
6	6	1	0.52271	30.2043	849.971	34.2959	2907.04	42.9773	4.17424	4420.3	47.8079	14675.6	0.301490		
7	7	1.16667	0.46687	30.2761	1049.20	81.2448	3003.03	43.2078	4.49489	5005.25	346.149	20040.7	0.4312		
8	8	1.33333	0.51481	30.2862	946.48	72.3172	3257.42	43.1246	4.35114	6059.37	185.203	21024.4	0.448799		
9	9	1.5	0.49872	30.2807	999.92	76.6829	3420.47	43.1998	4.43836	7285.49	264.885	24446.9	0.523701		
10	10	1.66667	0.46687	30.2761	1049.20	81.2448	3003.03	43.2078	4.49489	5005.25	346.149	20040.7	0.4312		
11	11	1.83333	0.44191	30.1494	1097.74	85.5041	3764.42	43.2015	4.55808	9405.99	631.683	31050.3	0.7087		
12	12	2	0.42349	30.1099	1162.24	80.0974	3930.96	43.4708	4.43812	10540.1	921.79	35736.3	0.803001		
13	13	2.16667	0.41017	30.1729	1251.6	97.8488	4088.18	43.5778	4.70239	11420.7	433.48	38824.1	0.906		
14	14	2.33333	0.40891	30.3445	1371.2	108.91	4241.15	43.6956	4.82419	13192	727.43	44085.4	1.0198		
15	15	2.5	0.40241	30.3979	1494.94	118.643	4427.14	43.8124	4.93381	14689	846.273	48132.7	1.1376		
16	16	2.66667	0.44336	30.2949	1491.48	120.53	4492.06	43.9461	5.12019	14840.8	976.303	51024.3	1.2723		
17	17	2.83333	0.50911	30.1897	1797.51	139.184	4779.32	44.0337	5.09892	18046	1114.49	57882.1	1.4079		
18	18	3	0.5999	30.1884	1843.33	149.613	4917.81	44.2389	5.14818	19909.3	1244.1	62799.9	1.5431		
19	19	3.16667	0.71672	30.2078	1941.63	156.428	5096.81	44.4049	5.18841	21850.9	1420.72	67896.7	1.7291		
20	20	3.33333	0.68078	30.2788	1989.89	160.847	5248.08	44.4778	5.14844	23843.4	1383.87	73146.4	1.9022		
21	21	3.5	0.90086	30.3058	2047.35	165.365	5422.43	44.7663	5.15047	25880					
22	22	3.66667	1.02034	30.2894	2097.58	169.444	5566.73	44.9223	5.1312	27959					
23	23	3.83333	1.23016	30.189	2146.44	173.71	5732.48	45.1149	5.14971	30236					
24	24	4	1.37332	30.1703	2199.38	177.927	5831.44	45.298	5.14449	32334					
25	25	4.16667	1.52179	30.2943	2265.47	182.293	6140.46	45.4747	5.14941	34424					
26	26	4.33333	1.46049	30.2824	2416.4	186.494	6300.23	45.7021	5.14842	37404.3	2441.87	107952	3.0273		
27	27	4.5	1.83416	30.2974	2534.91	200.904	6214.95	45.9083	5.14468	39578.2	2854.77	114467	3.2325		
28	28	4.66667	2.03994	30.2894	2487.87	216.284	6303.27	46.1305	5.13784	42236.1	3075.05	120470	3.4547		
29	29	4.83333	2.30149	30.2441	2744.42	224.57	6390.24	46.3478	5.14688	44020.5	3280.57	124841	3.6982		
30	30	5	2.55879	30.2292	2906.43	234.661	6374.27	46.5971	5.18376	47927.1	3534.13	133235	3.9213		
31	31	5.16667	2.87899	30.2145	2960.12	234.209	6484.27	46.8781	5.19489	50982.2	3740.34	139713	4.2023		
32	32	5.33333	3.05003	30.1863	3260.15	241.374	6510.44	47.0344	5.19087	52443.4	3939.72	146209	4.3388		
33	33	5.5	2.77264	30.1520	3348.73	242.669	6510.04	47.1351	5.03062	53487.1	3932.41	152705	4.4599		
34	34	5.66667	2.42789	30.2020	3385.1	246.072	6510.4	47.2593	4.99441	54850.4	4020.48	159260	4.5833		
35	35	5.83333	3.05003	30.2108	3432.23	249.241	6509.48	47.2833	4.91897	56237.7	4127.78	165791	4.7073		
36	36	6	4.67491	30.2180	3478.93	252.419	6519.99	47.4549	4.80719	57994.6	4242.49	172318	4.8291		
37	37	6.16667	1.36933	30.2735	3529.87	256.582	6478.5	47.6649	4.84497	60224.4	4399.07	178794	4.9521		
38	38	6.33333	1.1484	30.2494	3520.42	259.21	6510.81	47.8721	4.8885	62754.9	4570.5	185807	5.1463		
39	39	6.5	1.05865	30.2468	3644.24	263.799	6524.46	48.0553	5.023	65413.2	4768.1	193184	5.4958		
40	40	6.66667	0.94013	30.2044	3508.14	268.409	6504.8	48.2491	5.08711	66527.3	4886.5	198339	5.5933		
41	41	6.83333	0.96235	30.1122	3424.33	269.032	6497.94	48.4297	4.97749	69961.7	5014.94	204384	5.7639		
42	42	7	0.98921	30.2364	3616.22	247.214	6514.81	48.6429	4.87849	71854.8	5460.78	211841	6.1467		
43	43	7.16667	1.23849	30.2414	3620.52	251.501	6514.75	48.8266	5.18894	75536.3	5395.28	217843	6.3508		
44	44	7.33333	1.01002	30.2189	3543.15	245.459	6509.81	49.2015	5.10982	77770.5	5540.79	224373	6.5237		
45	45	7.5	0.72084	30.2189	3574.1	249.301	6510.81	49.4076	5.13146	80306.5	5703.04	230891	6.7142		
46	46	7.66667	0.63094	30.2743	3778.38	252.939	6505.86	49.6476	5.14407	83331.9	5891.94	237397	6.9718		
47	47	7.83333	0.54516	30.2819	3973.48	259.891	6522.44	49.9308	5.1802	86005.3	6091.93	243951	7.2121		
48	48	8													
49	49	8.16667													
50	50	8.33333													

BakersYeast.M / Scores [Comp. 1]



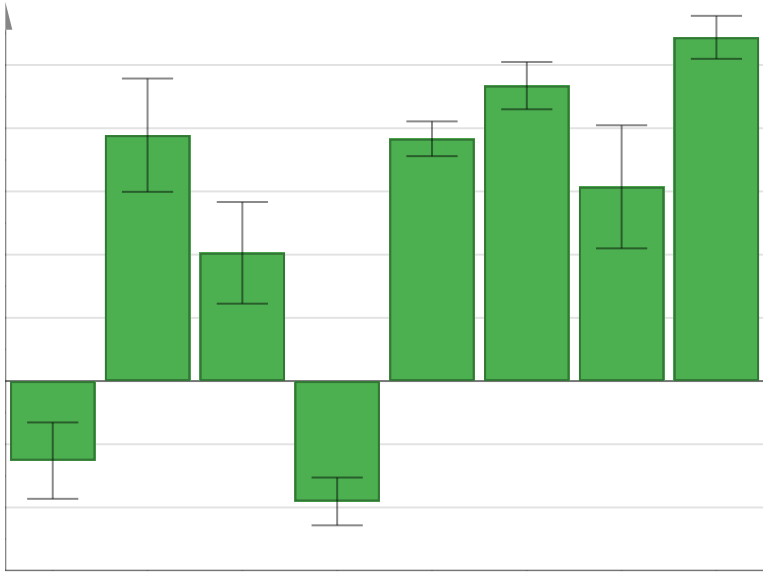
Example – Baker's Yeast

BakersYeast.M1
Scores [Comp. 1]



Observations - HOW

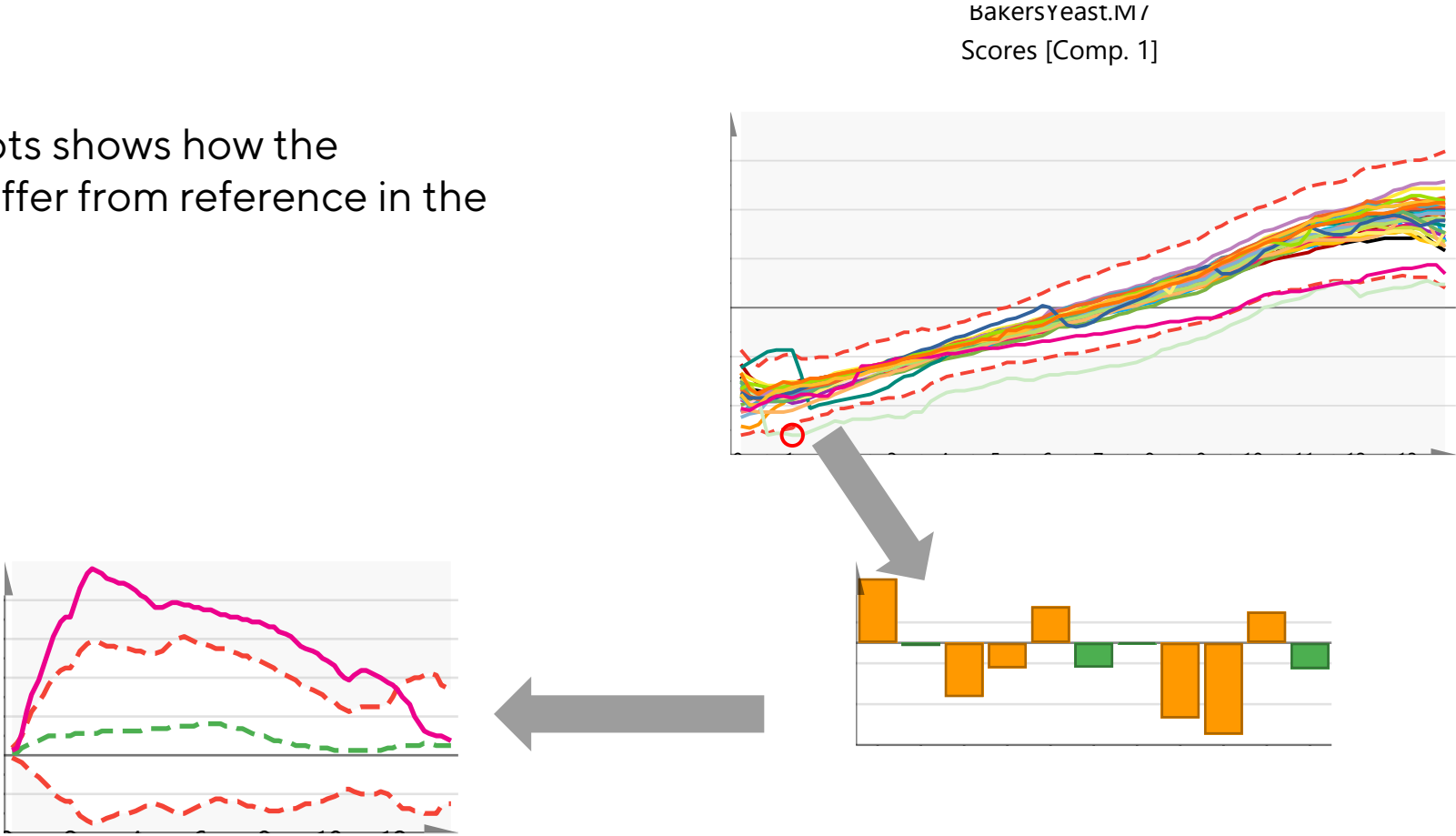
BakersYeast.M1 (OPLS), First model excl low QP2
Normalized to unit length



Variables - WHY

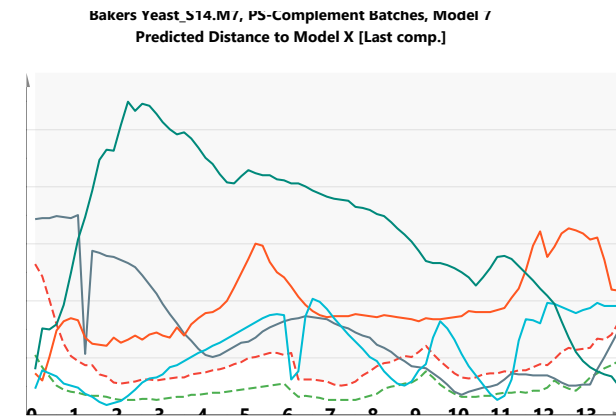
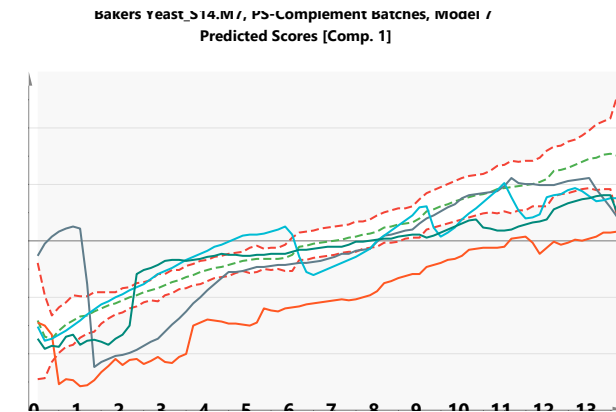
Contribution Plots - Scores

- Score contribution plots shows how the observation / group differ from reference in the model space



Validation of Batch Evolution Model (BEM)

- The poor performing and misbehaving batches are used to validate the model
- Will the model recognize batches with suboptimal behavior?
- Batches outside the ± 3 StDev limits are considered as definite deviators
- Batches outside the ± 2 StDev limits are considered as risky





Example: Cell Culture Evolution (BEM/BLM)

Modelling Cell Culture Metabolite Trajectories for Process Monitoring

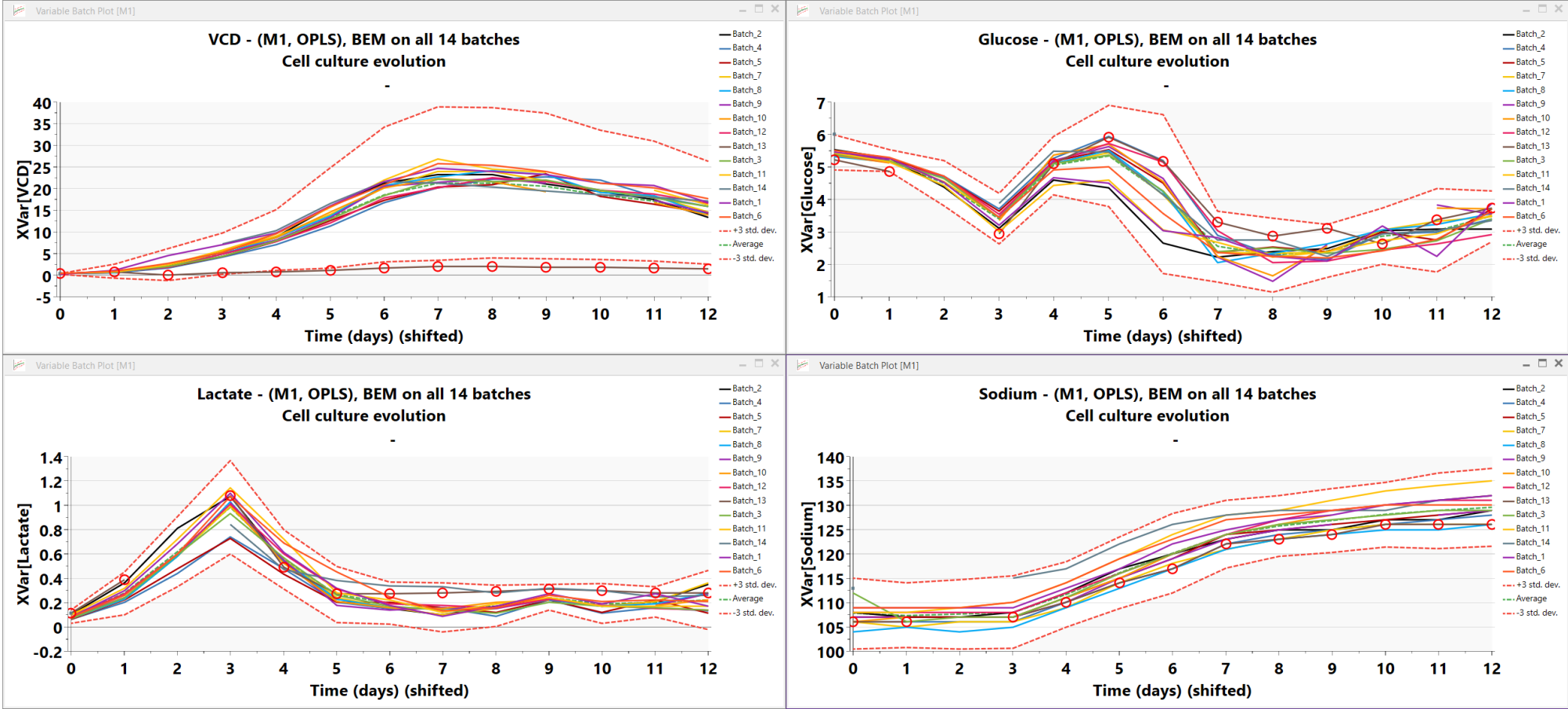
- Several fed-batch cell culture processes were performed for process development objectives.
- The experimental campaign involved processes performed under stable conditions and under unstable conditions.
- As typical parameters of the process performance metabolite measurements, cell characteristics as well as ion and gas concentration measurements were registered once per day.
- There are 14 batches. Process measurements, collected once per day, as well as initial and final batch conditions are available.

Process Evolution Measurements Used to Get the BEM

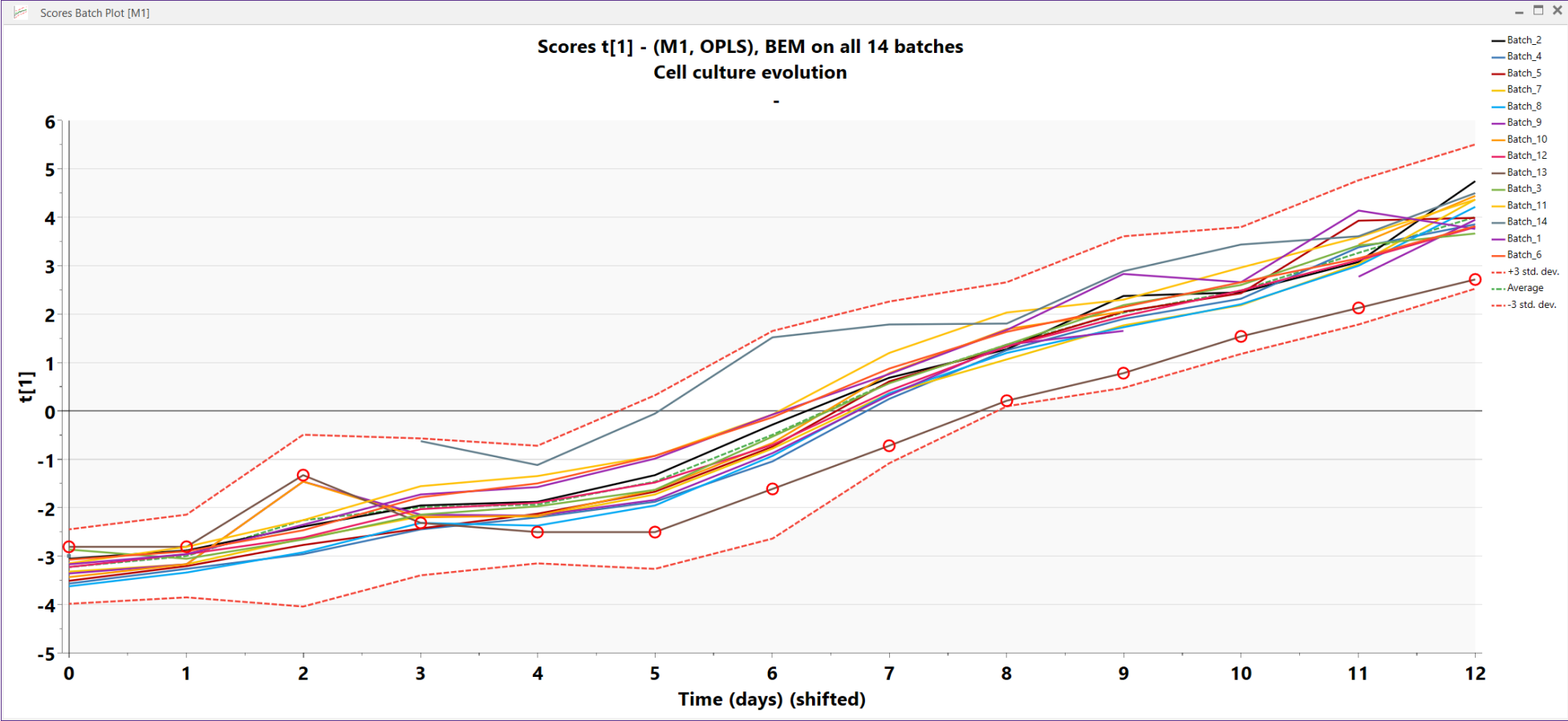
Dataset - Evolution data

1	2	3	4	5	6	7	8	9	10	11	12	13	14	
Primary ID	\$BatchID	Time (days)	VCD	Viability	Glucose	Lactate	Titer	pCO2	pH Offline	pO2	Osmolality	Potassium	Sodium	
41	40	Batch_7	0	0.324	98.2	5.4	0.07	0.0104639	49.5	7.106	91.2	306	9.4	106
42	41	Batch_7	1	0.713	98	5.19	0.23	0.0222557	48.6	7.09	92.5	306	9.3	105
43	42	Batch_7	2	2.036	99	4.56	0.58	0.0647377	36.1	7.102	90.5	304	9.1	106
44	43	Batch_7	3	5.492	99	3.45	0.99	0.165377	24.1	7.07	89.3	290	8.6	106
45	44	Batch_7	4	9.072	99	5.09	0.55	0.339043	38.9	7.094	83.8	309.5	9.4	109
46	45	Batch_7	5	14.1415	99	5.44	0.24	0.620911	58.8	7.085	69.9	319.67	9.7	114
47	46	Batch_7	6	20.1197	99	4.13	0.16	1.07201	70.3	7.09	64.9	315.5	9.4	118
48	47	Batch_7	7	23.984	98.6	2.37	0.1	1.84783	81.9	7.095	61.9	304.5	9.7	121
49	48	Batch_7	8	24.017	98.33	2.35	0.15	2.35966	81.1	7.095	55.6	308	10.9	123
50	49	Batch_7	9	21.848	96.7	2.37	0.24	2.8195	73.5	7.099	52.7	317.5	11.8	125
51	50	Batch_7	10	19.488	95	3.06	0.17	3.17237	77.9	7.103	57.7	331.5	13	126
52	51	Batch_7	11	18.207	89.13	3.32	0.21	3.43495	83.7	7.096	57.9	345	14.1	126
53	52	Batch_7	12	14.51	76.95	3.46	0.36	3.49327	78.1	7.1	77	364	15.5	126
54	53	Batch_8	0	0.31	99	5.32	0.07	0.0100164	54.4	7.064	91.4	303	9.2	104
55	54	Batch_8	1	0.63	99.1	5.15	0.23	0.0215393	46.4	7.1	89.3	300.5	9.1	105
56	55	Batch_8	2	1.829	98.9	4.66	0.58	0.0642541	35.4	7.104	90.3	298	8.9	104
57	56	Batch_8	3	4.872	99.3	3.44	1.03	0.179967	21.5	7.089	89.9	284	8.4	105
58	57	Batch_8	4	8.625	99.3	5.11	0.54	0.344157	41.3	7.07	79.8	307.5	9.2	109
59	58	Batch_8	5	13.635	99.35	5.51	0.23	0.681418	57.3	7.09	77.5	317	9.4	113
60	59	Batch_8	6	20.702	99.3	4.24	0.16	1.16472	68.8	7.09	58.3	309	9.4	117
61	60	Batch_8	7	22.644	98.06	2.05	0.15	1.76629	78.9	7.082	44.1	299.5	9.6	121
62	61	Batch_8	8	24.0935	97	2.34	0.16	2.2537	84.7	7.08	14.7	309	10.7	123
63	62	Batch_8	9	23.251	96	2.64	0.22	2.74823	77.1	7.102	47.3	317	11.8	124
64	63	Batch_8	10	19.175	93.9	3.05	0.17	3.11018	81.4	7.101	46	331.5	12.9	125
65	64	Batch_8	11	18.38	90.05	3.24	0.19	3.32718	85.7	7.095	54.8	349	14.1	125
66	65	Batch_8	12	13.958	79.2	3.56	0.28	3.49662	84	7.1	57.9	367	15.4	126
67	66	Batch_9	0	0.312	99	5.41	0.08	0.00943621	48.2	7.116	89.9	301	9.5	107

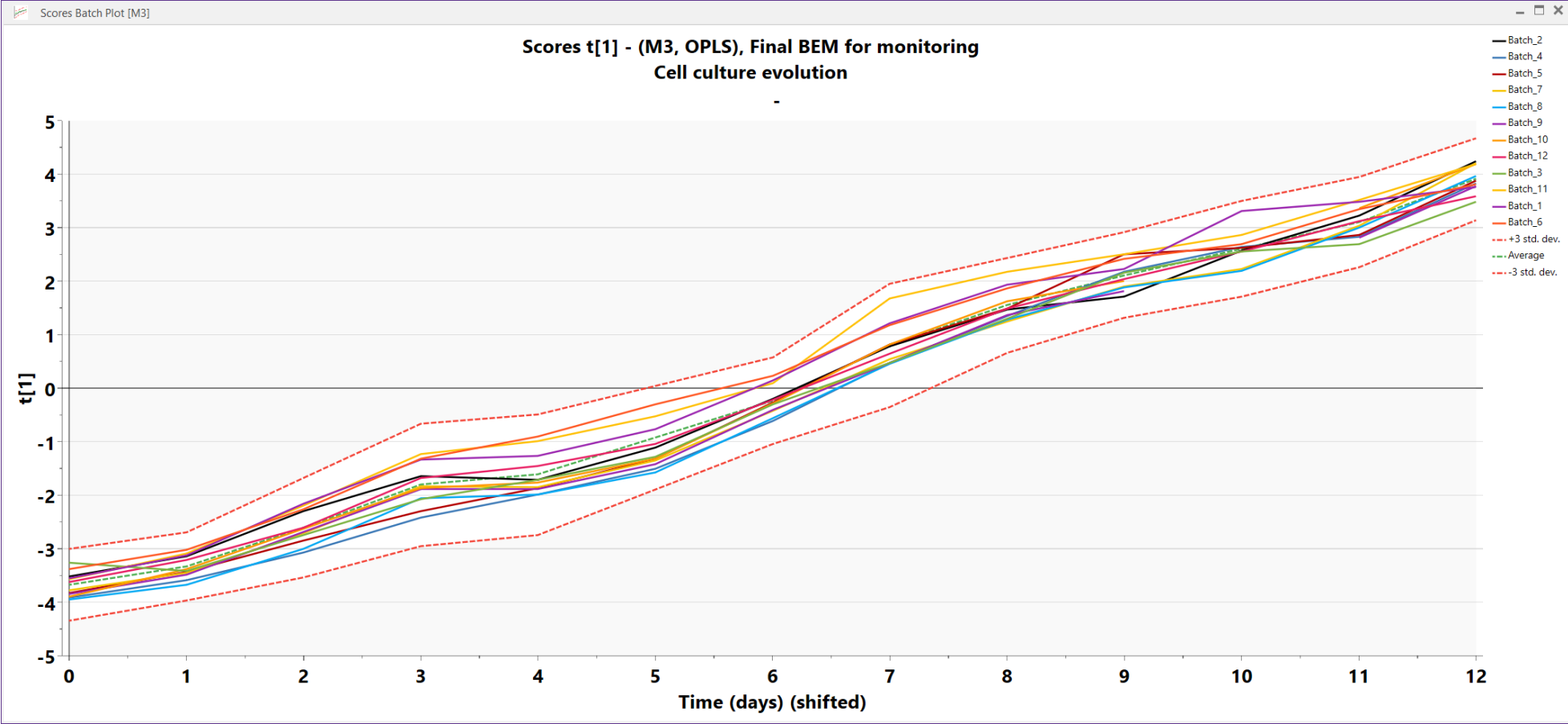
Batch Control Charts (BCC) of Individual Variables (Some Examples)



All Trajectories Summarized by the Score BCC



Result: Two Batches to Be Removed → Final BEM Representing Good Process



Batch Condition Data Used for BLMs

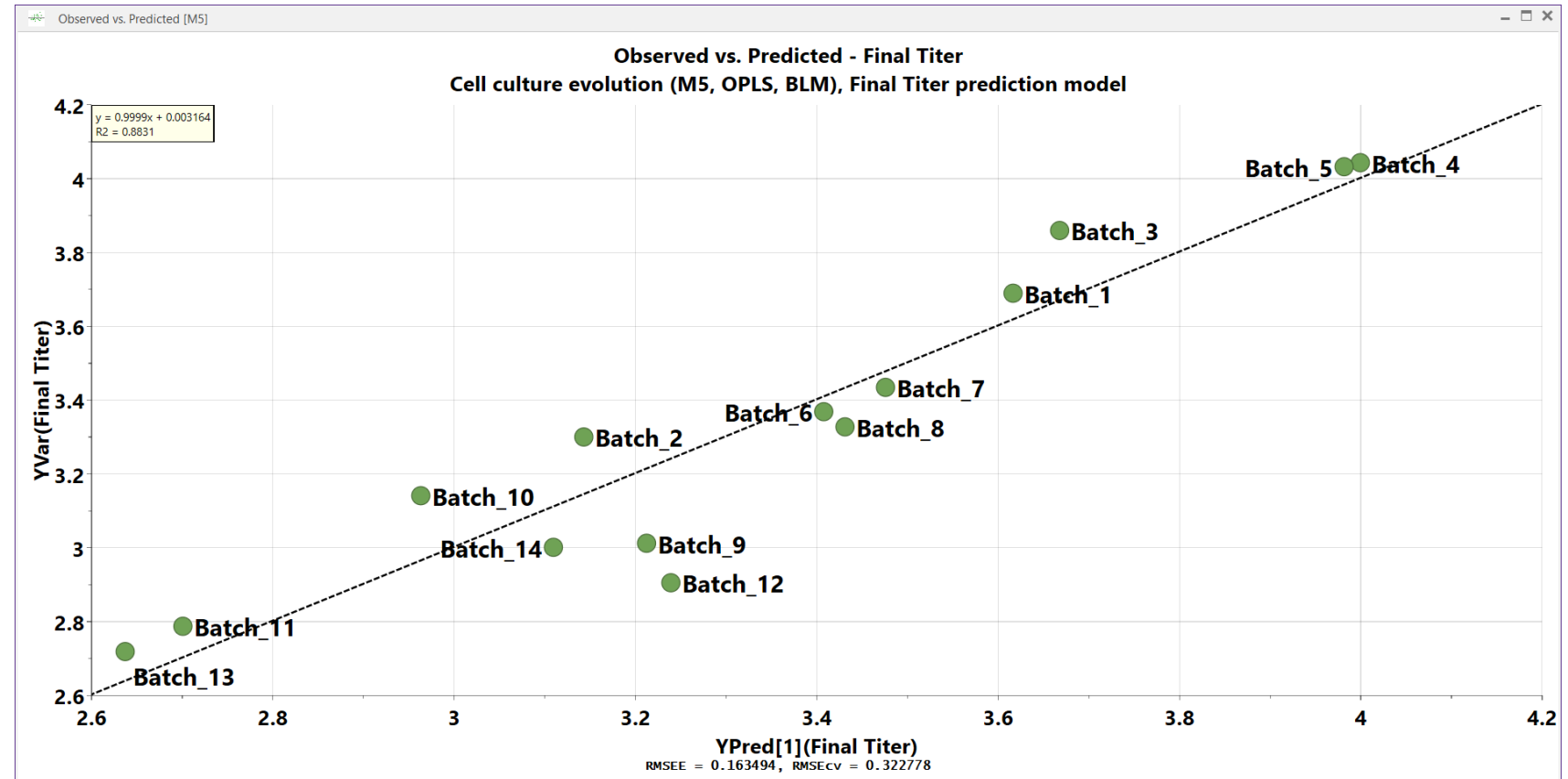
- A BLM is used to overview ALL data of completed batches
- A BLM can be used to predict final conditions
 - E.g. What process evolutions measurements correlate with final titer?
 - And when during cell culture lifetime?

	1	2	3	4	5	6	7	8	9	10	11	12
1	Primary ID	\$BatchID	Initial Glucose	Initial Lactate	Initial Osmolality	Initial Potassium	Initial Sodium	Initial VCD	Initial Viability	Initial CO2	Initial pH	Initial O2
2	\$Batch_1	Batch_1	5.45	0.08	308.5	9.7	109	0.275	99	50.3	7.1	97.4
3	\$Batch_2	Batch_2	5.53	0.1	307	9.6	108	0.3	96.5	47.7	7.109	88.7
4	\$Batch_3	Batch_3	5.38	0.07	308	10	112	0.3	97.3	52.5	7.075	83.1
5	\$Batch_4	Batch_4	5.44	0.06	307	9.5	106	0.325	99.1	40.9	7.08	88
6	\$Batch_5	Batch_5	5.42	0.06	308	9.5	107	0.325	99.1	40.4	7.08	96.5
7	\$Batch_6	Batch_6	5.5	0.08	315	9.6	109	0.3	98.2	53.4	7.084	100
8	\$Batch_7	Batch_7	5.4	0.07	306	9.4	106	0.325	98.2	49.5	7.106	91.2
9	\$Batch_8	Batch_8	5.32	0.07	303	9.2	104	0.3	99	54.4	7.064	91.4
10	\$Batch_9	Batch_9	5.41	0.08	301	9.5	107	0.3	99	48.2	7.116	89.9
11	\$Batch_10	Batch_10	5.36	0.1	300	9.4	106	0.325	99.4	49.9	7.105	88.6
12	\$Batch_11	Batch_11	5.44	0.1								108
13	\$Batch_12	Batch_12	5.38	0.08								90.5
14	\$Batch_13	Batch_13	5.22	0.11								102
15	\$Batch_14	Batch_14	6.01	0.11								94

	1	2	3	4	5	6	7
1	Primary ID	\$BatchID	Peak Titer	Peak VCD	Final Titer	Final VCD	Final Viability
2	\$Batch_1	Batch_1	3.6895	24.65	3.6895	20.7	90.8
3	\$Batch_2	Batch_2	3.42983	23.3	3.29939	17.5	89.3
4	\$Batch_3	Batch_3	3.99882	22.225	3.86063	17.8	92.4
5	\$Batch_4	Batch_4	4.26145	22.7	4.04247	18.2	94
6	\$Batch_5	Batch_5	4.25435	23.425	4.03154	16.4	93
7	\$Batch_6	Batch_6	3.62	25.7	3.37	20.2	92.7
8	\$Batch_7	Batch_7	3.49327	23.975	3.43495	18.2	89.13
9	\$Batch_8	Batch_8	3.49662	24.075	3.32718	18.4	90.05
10	\$Batch_9	Batch_9	3.23808	22.4	3.01178	17.325	91.45
11	\$Batch_10	Batch_10	3.40568	22.375	3.14113	16.65	89.8
12	\$Batch_11	Batch_11	2.93629	26.825	2.78789	19.6	91.8
13	\$Batch_12	Batch_12	3.08995	22.5	2.90643	18.8	94.3
14	\$Batch_13	Batch_13	2.92	1.925	2.72	1.6	93.4
15	\$Batch_14	Batch_14	3.18	21.625	3	18.1	92.4944

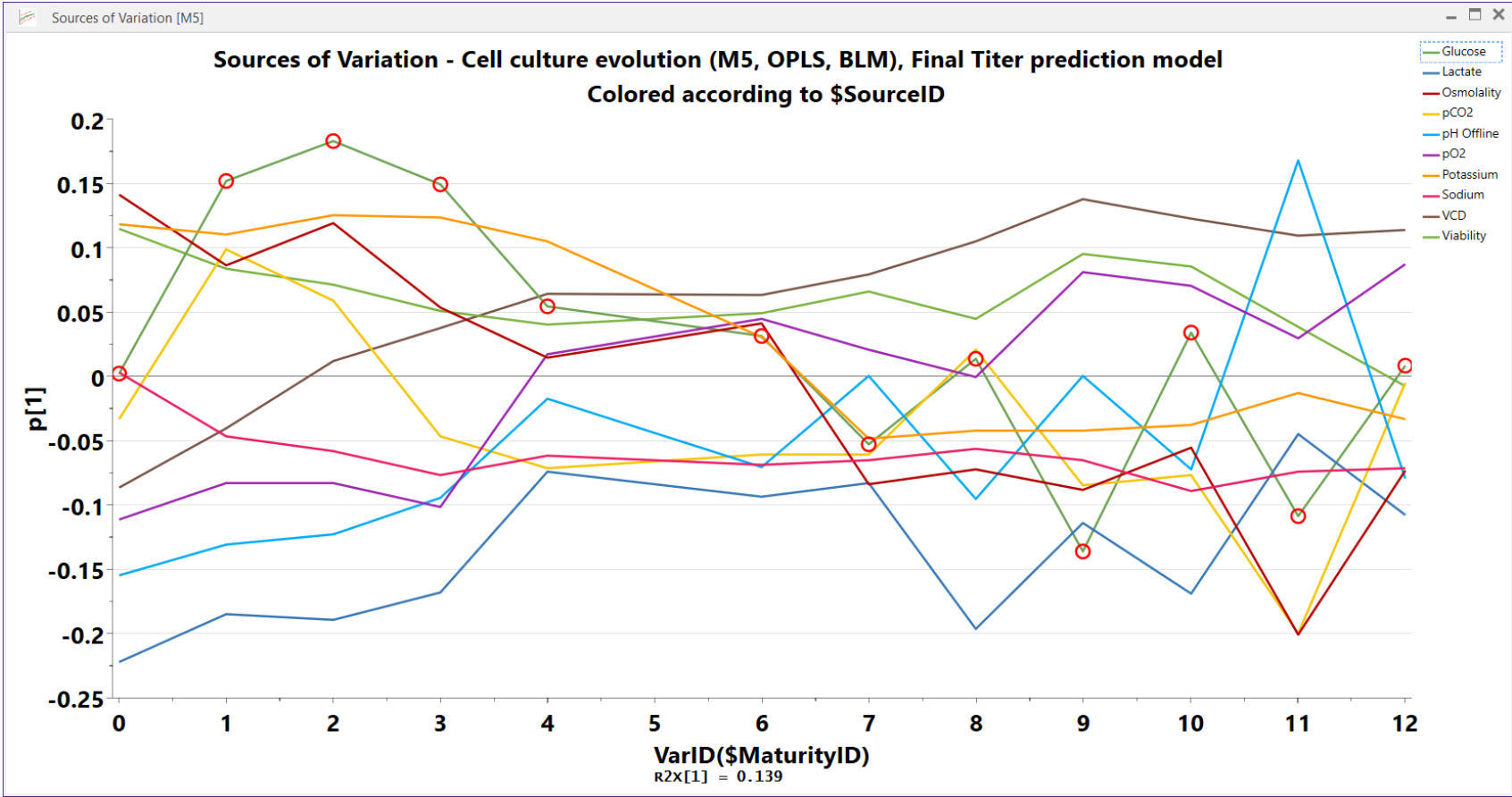
Example of BLM: Final Titer Prediction Model

- Strong agreement between measured final titer and model predictions



Sources of Variation Plot to Reveal Influence of Process Parameters

- The influence of the process parameters across lifetime of cell culture process visualized
- Glucose highlighted as an example
- Possibility to sharpen model by dropping non-influential parameters and/or time segments





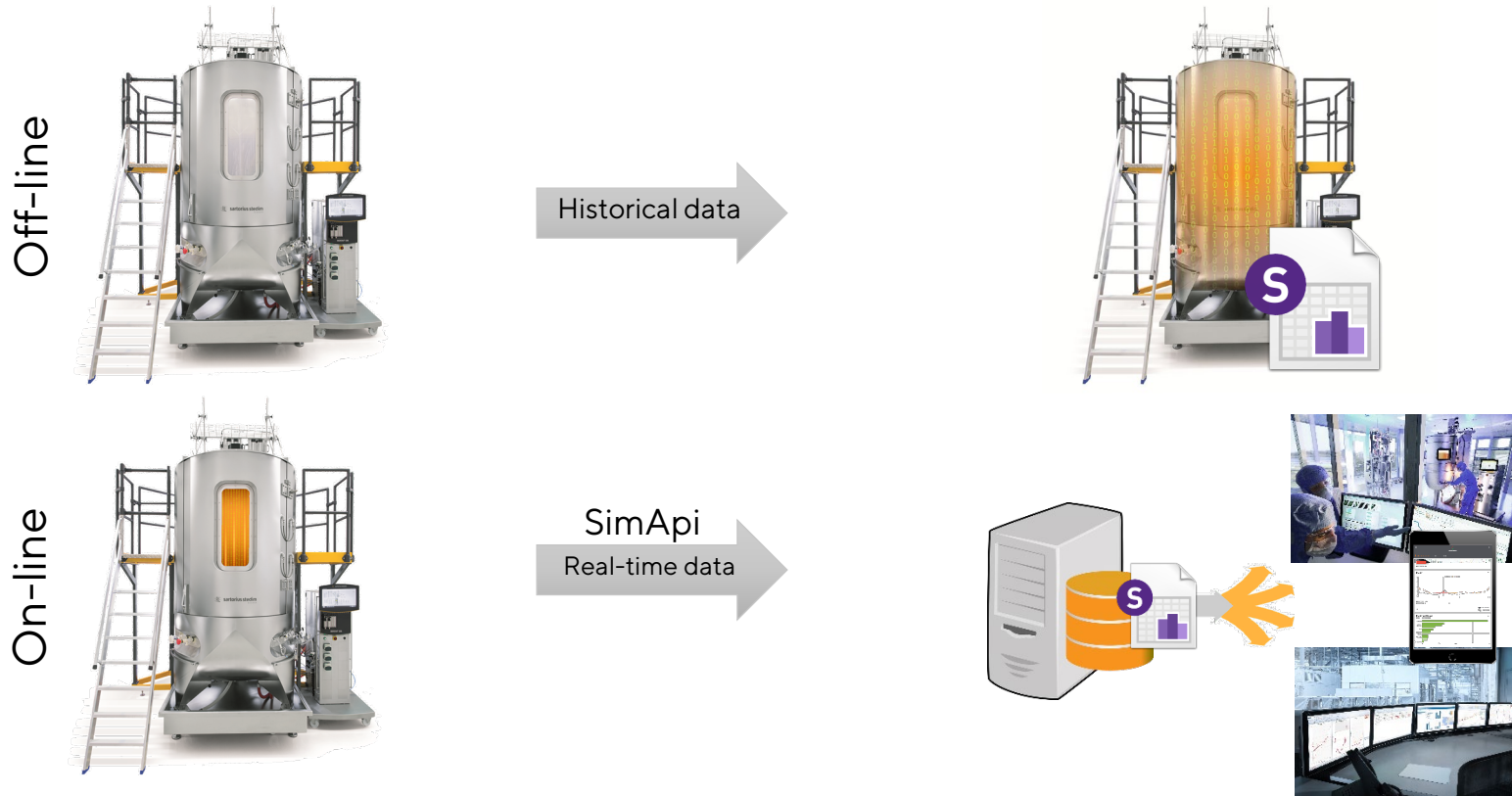
Demo (Cell Culture Dataset)

CONNECTION
ANALYSIS
DATA
SEARCHING
VERIFICATION

Conclusions From Demo Example

- This example shows a standard strategy to build multivariate models for cell culture processes with the long-term goal of enabling real-time process monitoring.
- The strategy is based on two types of models, the batch evolution model (BEM) and the batch level model (BLM).
- The objective of the BEM is to allow process monitoring of a cell culture process as it is evolving and as early as possible flag any deviation from a normal operating condition.
- The objective of the BLM is to summarize all data (initial conditions, process evolution measurements, final conditions,...) for completed batches to explore possibilities for further process optimization and process enhancement.
- The different models developed in this example suggest high quality data and a strong association between final titer and the way the cell cultures were progressing.

General Functionality of SIMCA Family



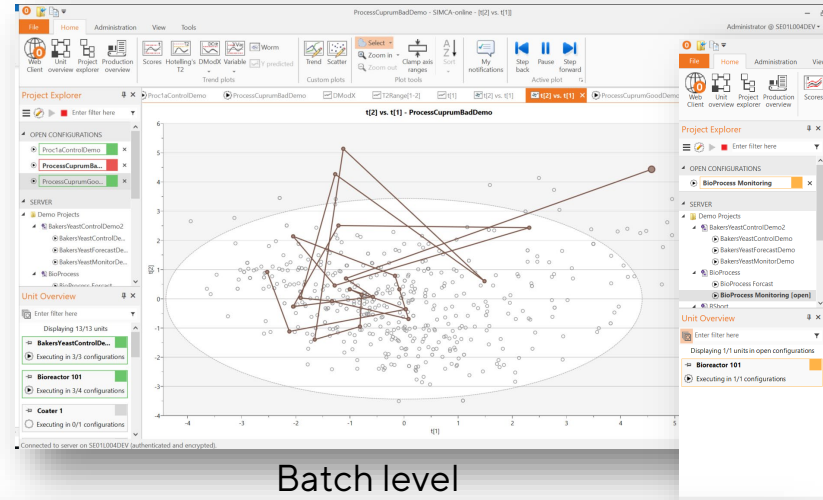
- SIMCA®
 - Creates models from historical data
 - Model objectives include
 - Finding deviating batches
 - Predicting a quality parameter
 - Resulting models are stored in a USP (Umetrics® SIMCA® Project) file

- SIMCA®-online
 - Samples process data at user specified frequency
 - Uses USP files as reference to new data
 - Visualizes the current state of the process

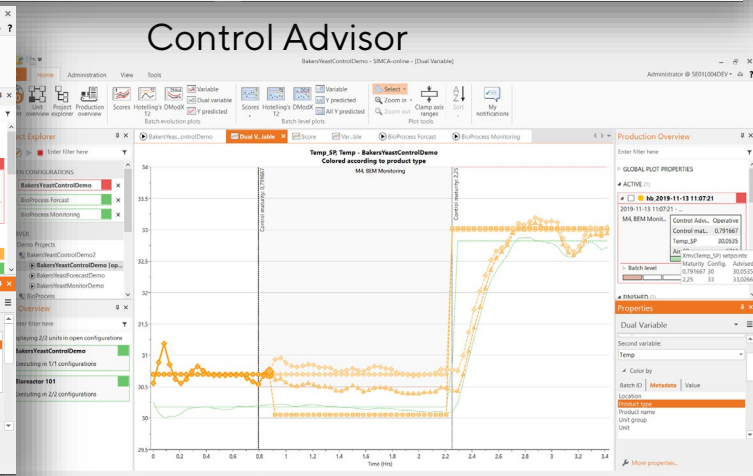
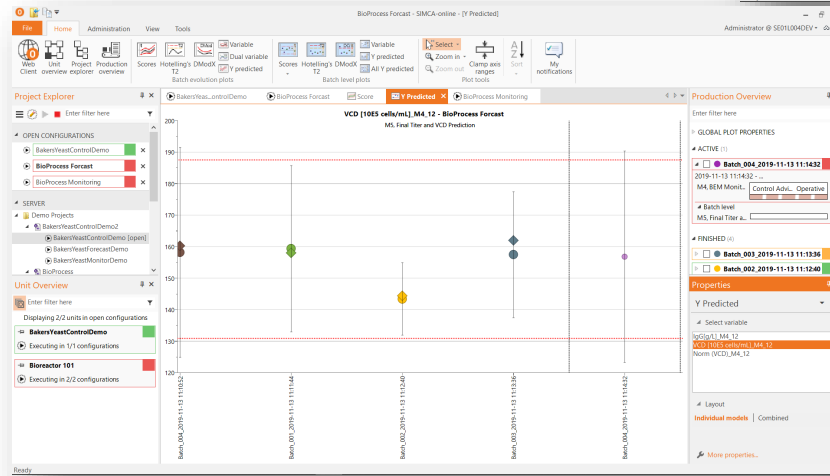
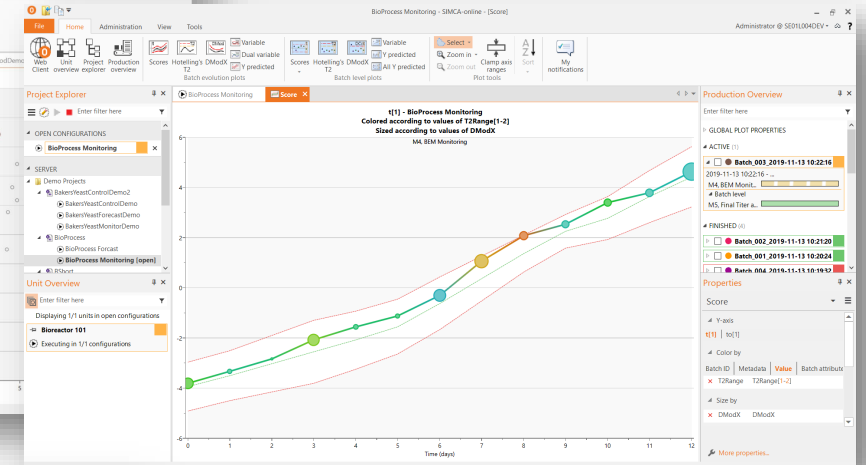
What can SIMCA®-online Do?

- **Monitor** – in real time
 - Create an ideal model of your process
 - Compare your process with the model in real time
- **Predict** – with confidence
 - Final quality prediction
 - Removes expensive measurements
 - All predictions has validity measures
- **Control** – at a glance
 - Forecast quality predictions
 - Learn how to optimize your process

Continuous



Batch evolution



Upcoming Webinars

(<https://www.sartorius.com/en/company/exhibition-conferences>)

The screenshot displays a web interface for Sartorius webinars. At the top, there are navigation tabs: 'UPCOMING EVENTS' (selected), 'PAST EVENTS', 'CALENDAR' (with a calendar icon), and 'SCHEDULE' (with a plus icon). Below the tabs is a search bar with a magnifying glass icon and the text 'Search'. The main content area is divided into two columns of event cards. Each card features a green square icon with a white checkmark, a 'STANDARD' tag, the event title, and the date and time. A vertical ellipsis menu is located to the right of each card. The events listed are:

Event Title	Date	Time
Design of Experiments (DOE) for the Beginner	TUE, JAN 26, 2021	03:00 PM - 04:00 PM CET
Multivariate Data Analysis (MVDA) for the Beginner	THU, JAN 28, 2021	03:00 PM - 04:00 PM CET
Lean-and-clean DOE using One-click analysis	TUE, FEB 16, 2021	03:00 PM - 04:00 PM CET
OPLS® in process modeling	THU, FEB 18, 2021	03:00 PM - 04:00 PM CET
Robust optimization made easy	TUE, MAR 2, 2021	03:00 PM - 04:00 PM CET
Analyzing batch process data, a step-by-step guide	THU, MAR 4, 2021	03:00 PM - 04:00 PM CET
From Design of Experiments to Design Space Estimation	TUE, MAR 23, 2021	03:00 PM - 04:00 PM CET
Multiblock Orthogonal Component Analysis (MOCA) – A Novel Tool for Data Integration	THU, MAR 25, 2021	03:00 PM - 04:00 PM CET