Automated Wort and Beer Quality Control Analysis

Kane, H.,¹ O'Shaughnessy, C.,² and Suoniemi-Kähärä, A.³ ¹Thermo Fisher Scientific, UK; ²Campden BRI, UK; ³Thermo Fisher Scientific, Finland





Introduction

We describe a system capable of producing fast and reliable results from several analytes by combining automated colorimetric detection and solid phase extraction techniques. The analyzer, Thermo ScientificTM GalleryTM Plus Beermaster, can determine bitterness and simultaneously do other colorimetric determinations (e.g. SO₂, FAN, pH, color, polyphenol and beta-glucan) from beer or wort samples. In addition many water quality parameters can be measured using the same analyzer.

In this study we present a comparison of bitterness measurement from beer and wort samples between a new automated method and the iso-octane extraction method. We also present method comparison studies from beer and wort samples for pH, color, FAN and SO₂.

The bitterness method uses a solid-phase extraction column integrated into an automated photometric analyzer. In the automated bitterness method, samples are first acidified then passed through the extraction column which binds bittering substances. The sample matrix is washed out and bittering substances are eluted and measured at 275 nm. Bitterness units are automatically calculated from absorbance results.

Materials and Methods

Instruments

Thermo Scientific Gallery Plus Beermaster, an automated discrete photometric analyzer, Thermo Fisher Scientific Oy, Vantaa, Finland

Methods

Methods are fully automated using bar-coded traceable system reagents. Bitterness calibration was performed using samples with known bitterness values. Colorimetric methods were calibrated with either a water based standard solution or samples with known values. Reference methods were European Brewery Convention (EBC) 9.8 for bitterness, EBC 8.10 and 9.10 for FAN, EBC 9.6 for Color, EBC 9.25.1 for SO₂ and EBC 9.35 and 8.17 for pH.

Samples

To establish robustness over a range of alcohol and color values, ten small pack beer samples and ten worts were analyzed in duplicate using the Beermaster and reference methods for, pH, Color, Bitterness, Free Amino Nitrogen (FAN), Sulphur Dioxide (SO_2) (Note: SO_2 was measured in beer sample only.)

The worts (Table 1) were frozen prior to analysis to ensure product stability. They were gently defrosted overnight at room temperature and then centrifuged at 3,600 rpm for 5 minutes to remove any cold trub that had formed upon freezing. Ten small pack beers with alcohol content ranging from 0.5 to 9% (Table 2), and covering a range of color values were analyzed in duplicate. All the beers were degassed by leaving them overnight on the bench in a conical flask. In the case of SO₂ analysis a fresh can of sample was opened immediately before analysis.

Table 1.	Worts	used	in	this	study	V
----------	-------	------	----	------	-------	---

Code	Туре
WA	1038 Ale
WB	1040 Ale
WC	1040 Ale
WD	Adjunct
WE	1050 Best Bitter
WF	16 Plato Lager
WG	11 Plato Lager
WH	11 Plato Lager
WI	1050 Best Bitter
WJ	16 Plato Lager

Table 2. Beers used in this study

Code	Туре	Declared ABV (%)
BA	Lager	0.00
BB	Lager	2.30
BC	Stout	2.80
BD	Ale	3.50
BE	Lager	3.80
BF	Stout	4.20
BG	Lager	4.80
BH	Ale	5.20
BI	Ale	6.60
BJ	Lager	9.00

Results and Discussion

Tables 3-11 summarize the mean and precision data for analyses of wort and beer samples using the Beermaster as compared to the reference methods.

Table 3. Summary of pH analysis results for ten different wort samples

		Measured ¹		Thermo S	Scientific	Method			Re	eference	Method			
Analysis	Sample	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
			5.42	5.00	F 44	0.025	0.025	4.00	4.00	4.00	0.000	0.000	0.005	
рн	WA	5.22	5.13	5.08	5.11	0.035	0.025	4.86	4.86	4.86	0.000	0.000	0.065	
	WB	5.22	5.22	5.21	5.22	0.007	0.005	5.12	5.12	5.12	0.000	0.000	0.033	
	WC	5.15	5.21	5.21	5.21	0.000	0.000	5.14	5.12	5.13	0.014	0.010	0.079	
	WD	5.07	5.08	5.06	5.07	0.014	0.010	4.96	4.96	4.96	0.000	0.000	0.058	
	WE	4.87	5.11	5.12	5.12	0.007	0.005	4.95	4.95	4.95	0.000	0.000	0.019	
	WF	5.26	5.25	5.25	5.25	0.000	0.000	5.22	5.23	5.23	0.007	0.005	0.126	
	WG	5.30	5.28	5.28	5.28	0.000	0.000	5.23	5.24	5.24	0.007	0.005	0.070	
	WH	5.31	5.26	5.25	5.26	0.007	0.005	5.22	5.22	5.22	0.000	0.000	0.090	
	WI		5.09	5.09	5.09	0.000	0.000	4.90	4.88	4.89	0.014	0.010	0.032	
	WJ	5.22	5.23	5.23	5.23	0.000	0.000	5.21	5.18	5.20	0.021	0.015	0.258	

¹These are the reference method mean readings for the fresh pre-frozen worts.

Tuble 4. Outlinning of privanaryois results for terrainerent beer sumple
--

		Measured ²		Thermo S	cientific	Method			Re	eference	Method		
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
	CB44	4.08	4 16	4 16	4 16	0.000	0.000	4 16	4 03	4 10	0.092	0.065	0 500
нα	BA	4.00	4.12	4.12	4.12	0.000	0.000	3.98	3.92	3.95	0.032	0.030	0.111
•	BB		4.37	4.34	4.36	0.021	0.015	4.28	4.21	4.25	0.049	0.035	0.212
	BC		5.03	5.02	5.03	0.007	0.005	3.93	3.96	3.95	0.021	0.015	0.009
	BD		4.21	4.32	4.27	0.078	0.055	4.00	4.12	4.06	0.085	0.060	0.241
	BE		4.39	4.30	4.35	0.064	0.045	4.26	4.29	4.28	0.021	0.015	0.379
	BF		4.62	4.63	4.63	0.007	0.005	3.98	4.03	4.01	0.035	0.025	0.026
	BG		4.12	4.11	4.12	0.007	0.005	4.00	4.05	4.03	0.035	0.025	0.176
	BH		4.20	4.20	4.20	0.000	0.000	3.81	3.80	3.81	0.007	0.005	0.008
	BI		4.27	4.33	4.30	0.042	0.030	4.25	4.28	4.27	0.021	0.015	0.486
	BJ		4.10	4.09	4.10	0.007	0.005	4.06	4.09	4.08	0.021	0.002	0.426

²This is the QC sample from a commercially available lager. The value is the mean of readings developed as a reference method between February 2013 and July 2013.

The standard deviations of the duplicate samples showed that the precision of the Beermaster relative to the reference pH method was similar for both beers and worts. Overall the precision of the Beermaster in measuring pH proved better than that of the reference method and was within acceptable levels for such an instrument. Statistical analysis using the two-sample *t*-test and the one-way ANOVA test suggested that in the majority of cases and based on current data there is no statistically significant difference (p-value >0.05) in pH measurements for beer and wort when using the Beermaster versus the reference pH method.

		Measured ¹		Thermo S	cientific	Method			Re	eference	Method		
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
Colour	WA		25.78	26.05	25.91	0.190	0.140	27.48	27.50	27.49	0.014	0.010	0.055
(EBC)	WB	38.00	31.09	31.12	31.10	0.021	0.015	33.50	33.55	33.53	0.035	0.025	0.008
	WC	35.00	31.15	30.95	31.05	0.141	0.100	33.63	33.65	33.64	0.014	0.010	0.025
	WD	12.00	11.90	11.88	11.89	0.014	0.010	12.85	12.88	12.87	0.021	0.015	0.012
	WE	37.00	36.77	36.74	36.76	0.021	0.015	38.65	38.70	38.68	0.035	0.025	0.010
	WF	20.00	19.40	19.44	19.42	0.028	0.020	20.15	20.18	20.17	0.021	0.015	0.021
	WG	11.00	11.48	11.48	11.48	0.000	0.000	11.55	11.60	11.58	0.035	0.025	0.164
	WH	13.00	13.10	13.12	13.11	0.014	0.010	13.33	13.33	13.33	0.000	0.000	0.029
	WI		39.59	39.55	39.57	0.028	0.020	40.98	41.05	41.02	0.049	0.035	0.018
	LM	21.00	19.28	19.32	19.30	0.028	0.020	20.48	20.50	20.49	0.014	0.010	0.012

	Table 5. Summary	v of color analy	vsis results for te	n different wort sam	ples
--	------------------	------------------	---------------------	----------------------	------

¹These are the reference method mean readings for the fresh pre-frozen worts.

|--|

		Measured ²		Thermo S	cientific	Method			Re	eference	Method		
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
	CB44	8.88	8.50	8.52	8.51	0.014	0.010	8.93	8.93	8.93	0.000	0.000	0.015
Colour	BA		7.76	7.77	7.77	0.001	0.005	8.13	8.10	8.12	0.021	0.015	0.029
(EBC)	BB		5.08	5.09	5.09	0.007	0.005	5.38	5.38	5.38	0.000	0.000	0.011
	BC		165.77	165.66	165.71	0.080	0.055	181.30	184.50	182.90	2.263	1.600	0.059
	BD		14.46	14.46	14.46	0.001	0.001	15.15	15.20	15.18	0.035	0.025	0.022
	BE		8.12	8.11	8.12	0.007	0.005	8.68	8.70	8.69	0.014	0.010	0.012
	BF		91.77	91.99	91.88	0.156	0.110	102.30	103.00	102.65	0.495	0.350	0.022
	BG		7.36	7.32	7.34	0.028	0.020	7.58	7.58	7.58	0.000	0.000	0.053
	BH		52.69	52.94	52.82	0.177	0.130	58.80	59.30	59.05	0.354	0.250	0.029
	BI		13.39	13.40	13.39	0.007	0.005	14.08	14.15	14.12	0.049	0.035	0.031
	BJ		13.12	13.11	13.11	0.007	0.005	13.68	13.73	13.71	0.035	0.025	0.027

²This is the QC sample from a commercially available lager. The value is the mean of readings developed as a reference method between February 2013 and July 2013.

The standard deviation of the duplicate samples showed that the precision of the Beermaster relative to the reference color method was similar for both beers and worts. Interestingly in all cases the color readings from the Beermaster were slightly lower than that of the reference method. As a result, statistical analysis using the two-sample *t*-test and the one-way ANOVA test suggested that based on the current data there was a statistically significant difference (p-value >0.05) in color measurements for beer and wort when using the Beermaster versus the reference color method.

To assess whether inclusion of a factor would compensate for these differences, the statistical tests were re-run using the original Beermaster values multiplied by a factor of 1.05. By applying a factor of 1.05 to the Beermaster data, statistical analysis using the two-sample *t*-test and the one-way ANOVA test suggested that in the majority of cases and based on current data there is no statistically significant difference (p-value >0.05) in color measurements for beer and wort when using the Beermaster versus the reference color method.

Table 7. Summary of bitterness analysis results for ten different wort samples

		Measured ¹		Thermo S	cientific	Method			Re	ference	Method		
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
Bitterness	WA		41.83	42.03	41.93	0.141	0.100	33.65	33.70	33.68	0.035	0.025	0.008
(BU)	WB	34.00	24.73	22.82	23.78	1.351	0.950	29.30	28.75	29.03	0.389	0.027	0.119
	WC	36.00	25.19	24.44	24.82	0.530	0.370	29.65	29.60	29.63	0.035	0.025	0.050
	WD	26.00	32.84	31.81	32.33	0.728	0.520	23.40	23.35	23.38	0.035	0.025	0.037
	WE	51.00	56.56	63.33	59.95	4.787	3.400	55.10	56.30	55.70	0.849	0.600	0.433
	WF	34.00	39.91	38.48	39.20	1.011	0.710	32.50	33.45	32.98	0.672	0.480	0.087
	WG	35.00	41.95	42.14	42.05	0.134	0.095	30.90	31.75	31.33	0.601	0.430	0.026
	WH	32.00	40.06	41.03	40.55	0.686	0.480	27.95	30.00	28.98	1.450	1.100	0.062
	WI		51.99	53.71	52.85	1.216	0.860	51.75	53.20	52.48	1.025	0.730	0.795
	WJ	32.00	37.9	33.84	35.87	2.871	2.000	31.30	32.10	31.70	0.566	0.400	0.293

¹These are the reference method mean readings for the fresh pre-frozen worts.

Table 8. Summary of bitterness analysis results for ten different beer samples

		Measured ²		Thermo S	cientific	Method			Re	eference	Method		
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
	CB44	15.90	18.33	17.66	17.99	0.474	0.330	15.63	16.03	15.83	0.283	0.200	0.114
Bitternes	BA		22.11	21.86	21.98	0.177	0.130	17.50	17.90	17.70	0.283	0.200	0.035
(BU)	BB		15.91	15.51	15.71	0.283	0.200	13.18	13.58	13.38	0.283	0.200	0.014
	BC		20.28	18.47	19.38	1.280	0.910	23.40	24.20	23.80	0.566	0.400	0.140
	BD		25.47	24.01	24.74	1.032	0.730	20.25	20.28	20.26	0.021	0.015	0.103
	BE		15.44	14.83	15.13	0.431	0.300	12.38	12.45	12.42	0.049	0.035	0.072
	BF		26.76	24.56	25.66	1.556	1.100	24.20	24.45	24.33	0.177	0.130	0.441
	BG		20.19	22.33	21.26	1.512	1.100	19.85	19.62	19.74	0.163	0.110	0.391
	BH		23.18	23.14	23.16	0.028	0.020	26.80	26.70	26.75	0.071	0.050	0.010
	BI		17.33	16.92	17.13	0.289	0.200	17.20	17.30	17.25	0.071	0.050	0.660
	BJ		26.14	26.01	26.08	0.092	0.065	22.25	22.30	22.28	0.035	0.025	0.012

²This is the QC sample from a commercially available lager. The value is the mean of readings developed as a reference method between February 2013 and July 2013.

The standard deviation of the duplicate samples showed that the precision of the Beermaster relative to the reference bitterness method was similar for both beers and worts, and was within acceptable levels for such an instrument. Statistical analysis using the two-sample *t*-test and the one-way ANOVA test suggested that in the majority of cases and based on current data there is no statistically significant difference (p-value >0.05) in bitterness measurements for beer and wort when using the Beermaster versus the reference bitterness method.

		Measured ¹		Thermo S	cientific	Method			R	eference	Method	1	
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
FAN	WA		131.21	132.23	131.72	0.721	0.51	146.80	142.10	144.45	3.323	2.400	0.119
(mg/L)	WB	212.00	201.52	207.94	204.73	4.540	3.20	229.90	207.90	218.90	15.556	11.000	0.433
	WC	208.00	199.52	198.11	198.82	0.997	0.70	208.00	207.90	207.95	0.071	0.050	0.049
	WD	315.00	303.78	300.61	302.20	2.242	1.60	279.30	290.30	284.80	7.778	5.500	0.202
	WE	172.00	186.37	185.02	185.70	0.955	0.67	169.70	179.50	174.60	6.930	4.900	0.267
	WF	269.00	250.41	253.47	251.94	2.164	1.50	231.90	240.00	235.95	5.728	4.100	0.168
	WG	175.00	176.09	175.33	175.71	0.537	0.38	177.80	171.20	174.50	4.667	3.300	0.778
	WH	179.00	188.07	187.43	187.75	0.453	0.32	171.20	177.40	174.30	4.384	3.100	0.145
	WI		168.90	169.30	169.10	0.283	0.20	156.20	161.20	158.70	3.536	2.500	0.151
	WJ	263.00	258.55	259.64	259.10	0.771	0.54	231.80	243.40	237.60	8.202	5.800	0.168

Table 9. Summary of FAN analysis results for ten different wort samples

¹These are the reference method mean readings for the fresh pre-frozen worts.

Table 10. Summary of FAN analysis results for ten different beer samples

		Measured ²		Thermo S	cientific l	Method	thod Reference Method						
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
	CB44	76.68	69.82	69.46	69.64	0.255	0.180	81.33	79.72	80.53	1.138	0.81	0.048
FAN	BA		60.83	60.96	60.89	0.092	0.065	57.10	58.40	57.75	0.919	0.650	0.130
(mg/L)	BB		28.96	30.30	29.63	0.949	0.670	31.80	31.30	31.55	0.354	0.250	0.227
	BC		12.26	11.76	12.01	0.354	0.250	19.60	19.80	19.70	0.141	0.100	0.022
	BD		34.49	35.61	35.05	0.792	0.560	36.20	35.60	35.90	0.424	0.300	0.409
	BE		96.01	95.09	95.55	0.651	0.460	92.10	90.30	91.20	1.273	0.900	0.145
	BF		71.23	71.71	71.47	0.339	0.240	75.10	73.70	74.40	0.990	0.700	0.157
	BG		58.04	59.06	58.55	0.721	0.510	58.00	57.40	57.70	0.424	0.300	0.387
	BH		26.50	26.61	26.55	0.077	0.055	41.10	40.60	40.85	0.354	0.250	0.011
	BI		112.56	109.59	111.07	2.101	1.500	110.70	108.30	109.50	1.697	1.200	0.561
	BJ		107.97	108.92	108.45	0.674	0.480	105.10	102.50	103.80	1.838	1.300	0.184

²This is the QC sample from a commercially available lager. The value is the mean of readings developed as a reference method between February 2013 and July 2013.

The standard deviation of the duplicate samples showed that the precision of the Beermaster in measuring FAN proved better than that of the reference method (especially in the case of worts) and was within acceptable levels for such an instrument. Statistical analysis using the two-sample t-test and the one-way ANOVA test suggested that in the majority of cases and based on current data there is no statistically significant difference (p-value >0.05) in FAN measurements for beer and wort when using the Beermaster versus the reference FAN method.

		Measured ²		Thermo S	cientific	Method							
Analysis	Sample				Mean	Std Dev	SE Mean			Mean	Std Dev	SE Mean	P-Value
	230	5.60	4.81	4.94	4.88	0.092	0.065	5.60	5.12	5.36	0.339	0.240	0.302
SO2	BA		3.10	3.00	3.05	0.071	0.050	4.80	4.96	4.88	0.113	0.080	0.033
(mg/L)	BB		5.10	5.10	5.10	0.000	0.000	5.76	5.44	5.60	0.226	0.160	0.197
	BC		0.10	0.10	0.10	0.000	0.000	1.12	1.44	1.28	0.226	0.160	0.086
	BD		1.30	1.30	1.30	0.000	0.000	1.60	1.76	1.68	0.113	0.080	0.132
	BE		11.90	12.10	12.00	0.141	0.100	12.32	12.80	12.56	0.339	0.240	0.227
	BF		0.80	0.90	0.85	0.071	0.050	2.40	2.08	2.24	0.226	0.160	0.076
	BG		4.00	4.10	4.05	0.071	0.050	2.91	3.01	2.96	0.071	0.015	0.004
	BH		2.40	2.40	2.40	0.000	0.000	2.24	1.92	2.08	0.226	0.160	0.295
	BI		1.80	1.80	1.80	0.000	0.000	1.60	1.28	1.44	0.226	0.160	0.266
	BJ		15.10	19.70	17.40	3.253	2.300	16.32	16.80	16.56	0.339	0.240	0.778

	Table 11. Summar	v of SO ₂ ana	lysis results for	r ten different k	beer samples
--	------------------	--------------------------	-------------------	-------------------	--------------

²This is the QC sample from a commercially available lager. The value is the mean of readings developed as a reference method between February 2013 and July 2013.

The standard deviation of the duplicate samples showed that the precision of the Beermaster relative to the reference SO₂ method was similar. Overall the precision of the Beermaster in measuring SO₂ proved better than that of the reference method and was within acceptable levels for such an instrument. Statistical analysis using the two-sample t-test and the one-way ANOVA test suggested that in the majority of cases and based on current data there is no statistically significant difference (p-value >0.05) in SO₂ measurements for beer and wort when using the Beermaster versus the reference SO₂ method.

Ten replicates of a commercially available canned lager were analyzed for pH, color, bitterness, FAN and SO₂ using the Beermaster and traditional reference methods. Table 12 summarizes the mean and precision data for these analyses.

Table 12. Summary of pH, color, bitterness, FAN and SO₂ analysis results for ten replicates of a single brand of beer

	рН		Color (EBC)		Bitterness (BU)		FAN (mg/L)		SO ₂ (mg/L)	
	Beermaster	Reference	Beermaster	Reference	Beermaster	Reference	Beermaster	Reference	Beermaster	Reference
Mean	4.14	4.03	7.32	7.58	22.16	19.74	58.16	57.75	3.97	2.96
SD	0.016	0.024	0.032	0.021	0.831	0.185	0.596	0.706	0.106	0.253
95% Confidence Interval for Mean	4.13, 4.15	4.01, 4.04	7.30, 7.35	7.56, 7.59	21.56, 22.75	19.60, 19.87	57.73, 58.59	57.25, 58.26	3.89, 4.05	2.78, 3.14

The precision of the Beermaster, as expressed in the standard deviation of the ten replicates showed with the current data and this brand of beer, precision was greatest for pH measurement and least for bitterness. The current data also suggested that the Beermaster was more precise in the measurement of pH, FAN and SO₂ compared with the reference methods used.

Conclusion

Based on the data obtained during this study, the Thermo Scientific Gallery Plus Beermaster analyzer has been shown to provide comparable performance to established methods in the measurement of pH, bitterness, FAN and SO2 In the case of color measurement, the Gallery Plus Beermaster had similar precision to the reference method but consistently gave slightly lower results suggesting that a factor of 1.05 be included to compensate for these differences in the methodology. When this factor was included, the Gallery Plus Beermaster was comparable to the established method. Analysis of ten replicates of the same brand of beer showed that the Gallery Plus Beermaster had greater precision in the measurement of pH, FAN and SO₂ compared to the reference methods used in this study. Precision values for all analyses are within acceptable levels for spectrophotometers in the brewing industry. The Gallery Plus Beermaster has been proven to provide faster results when compared to time consuming traditional methods. The low reagent and water volumes required for analysis not only reduce reagent costs but also reduce the amount of waste produced, thereby providing analysis with low environmental impact. The new bitterness measurement uses environmentally safe reagents without the requirement to use harmful iso-octane in routine analysis. The analyzer is very straightforward to use and requires minimal training or skills to run and maintain.

www.thermoscientific.com

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details



A Thermo Fisher Scientific Brand

North America Tel: 1-800-232-3342 | email: sales.industrial@thermofisher.com Tel: +358 9 329 100 | email: sales.cdx.fi@thermofisher.com Europe