

# Comparability Study of Manual and Automated Particle Characterization with MFI

*Dan Bach Kristensen, Senior Pharmaceutical Scientist and Trine Meiborg-Sloth, Pharmaceutical Technician  
Takeda Pharma A/S, Langebjerg, Denmark*

*Scott Mack, Senior Development Scientist, Angelica Olcott, MFI Product Manager and Erik Gentalen, Director of Applications  
ProteinSimple, Santa Clara, CA, USA*

## Introduction

Micro-Flow Imaging (MFI) has become a standard application for particle analysis of protein formulations<sup>1</sup> because of its ability to easily detect particle size and morphology for a diverse range of particle contaminants, including translucent protein fragments and silicone micro-droplets. Recently, the US Pharmacopeial Convention proposed changes to the USP <787> standard, recommending the use of particle morphology detection through technologies such as Micro-Flow Imaging for sub-visible particle characterization<sup>2,3</sup>.

Higher throughput and standardization are needed to meet the demand for biopharmaceutical testing of sub-visible particle characterization. The MFI 5000 series instruments can be automated with the addition of the Bot1 autosampler which reduces the hands-on time for sample evaluation. Because the automated mode requires a different sample presentation, the method parameters are often modified in order to provide the same results in automated mode as it does in manual mode. Here we outline the process for transferring a manual protocol to an automated format with the Bot1 autosampler and the MFI 5000 series, with regard to sample volume and mixing parameters.

To demonstrate the comparability of running the MFI system in manual and automated modes, we ran a model protein system comprised of 1% BSA at multiple sites. Each sample was run at least eight times in both the manual and automated modes at two different laboratories. No statistically relevant differences were observed between manual and automated runs at either facility.



## Materials and Methods

### SAMPLE PREPARATION AND INSTRUMENT OPERATION

- 1% BSA sample was prepared, heated at 60 °C overnight (at least 12 hours) with shaking at 1200 rpm at the ProteinSimple facility. The sample was then aliquotted and stored at –80 °C. Frozen aliquots were shipped to the Takeda facility on dry ice and stored at –80 °C until use.
- Flow cell was washed with 1 mL of 10% Triton-X 100 in distilled, deionized (DDI) water (0.2 micron filtered) and rinsed 4X with 1 mL water before each run.
- A fourth stirring step was added to the protocol to ensure particles did not settle during the sampling process on the Bot1 (see **Table 1**).
- Manual pipetting steps were performed by turning off the Bot1 autosampler and manually pipetting sample into the MFI 5200 system via the sample inlet port. The same analysis method was used for both the automated and manual protocols (**Figure 1**).

### METHOD TRANSFER

Translating a manual method to an automated method requires modification of the protocol to ensure comparable results due to differences in sampling handling between the two methods.

Method transfer considerations include:

- **Changes in sample introduction** — manual pipetting or stirrer with syringe barrel mix differently than an automated pipettor. The user can control the duration and speed of mixing with the automated pipettor.
- **Suitability of sample type** — manual or automated sample requirements are similar; highly viscous or extremely concentrated samples may not be suited for automation.
- **Sample volumes** — sample and purge volumes may change due to new labware and fluid path.
- **Ambient temperature** — samples should be stable at room temperature during an automated run.

OPERATION	LIQUID	VOLUME (ML)
Flush	Filtered water or filtered buffer	0.90
Flush	Filtered water or filtered buffer	0.90
Flush	Filtered water or filtered buffer	0.90
Dry system		
Flush	Filtered water or filtered buffer	0.90
Optimize Illumination	Filtered water or filtered buffer	0.22
<b>Baseline</b>	Filtered water or filtered buffer	0.70
Flush	Filtered water or filtered buffer	0.90
Dry system		
Stir 4 cycles, speed 5	Sample (well 1)	0.90
Flush	Sample (well 1)	0.50
Stir 4 cycles, speed 5	Sample (well 1)	0.90
Optimize Illumination	Sample (well 1)	0.22
Stir 4 cycles, speed 5	Sample (well 1)	0.70
<b>Analysis Run</b>	Sample (well 1)	0.70
Flush	Filtered water or filtered buffer	0.90
Flush	Filtered water or filtered buffer	0.90
Flush	Filtered water or filtered buffer	0.90

**TABLE 1.** Description of automated Protocol D used for 1% BSA solution. Original version of Protocol D uses 0.90 mL for Flush step. Steps from 'Flush' after Baseline through 'Analysis Run' are repeated for each sample in the run. Stir of four cycles was used prior to each sample analysis to ensure particles remained in suspension.

Sample volumes should be chosen based on particle concentration in the sample, as described in the next section. Sample volumes should also be adjusted further to accommodate the sample delivery format of the instrument configuration. Thus, moving from a manual to automated protocol may require changes to sample volume and flush volume initially to achieve accurate results. This is due to the fixed sample delivery format which includes 1–2 mL deep well plates, the 1 mL pipette tips and the interface to the inlet port (see **Table 2**).

### Sample Analysis Report

<b>Method Identification</b> Name: For Bot1-0.7_0.0 Type: Sample Analysis and Report  Root Folder  Hardware Configuration: Set Point: 3      Flow Cell Model: 100µm(SP3) LED Slider Switch: NA Stir Speed(rpm) - NA Magnification Dial: NA Direction - NA Diaphragm Control: NA Stir before Run(min) - 0.0		<b>Image Frames</b> Storage Trigger: Capture all images Max Frames: 20000      Storage Space: 0.4334 GB Discard Empty Frames: <input type="checkbox"/> Seconds between Frames:																																																																																			
<b>Run Setup</b> Termination Type: Sample Dispensed(ml)      Edge Particle Rejection: <input type="checkbox"/> Total Available Volume(ml): 0.90      Fill Particles: <input checked="" type="checkbox"/> Purge Volume (ml): 0.00      Consecutive Runs: 1 Analyzed Volume(ml, approximate): 0.77		<b>Report</b> Chart Type: Histogram <table border="1"> <thead> <tr> <th>Axis</th> <th>Parameter 1</th> <th>Parameter 2</th> <th>Min</th> <th>Max</th> <th>Step</th> </tr> </thead> <tbody> <tr> <td>X</td> <td>ECD(µm)</td> <td>-</td> <td>1.00</td> <td>100.00</td> <td>1.00</td> </tr> <tr> <td>Y Left</td> <td>Count(#)</td> <td>-</td> <td>0.00</td> <td>MAX</td> <td>-</td> </tr> <tr> <td>Y Right</td> <td>-</td> <td>-</td> <td>-</td> <td>-</td> <td>-</td> </tr> </tbody> </table>		Axis	Parameter 1	Parameter 2	Min	Max	Step	X	ECD(µm)	-	1.00	100.00	1.00	Y Left	Count(#)	-	0.00	MAX	-	Y Right	-	-	-	-	-																																																										
Axis	Parameter 1	Parameter 2	Min	Max	Step																																																																																
X	ECD(µm)	-	1.00	100.00	1.00																																																																																
Y Left	Count(#)	-	0.00	MAX	-																																																																																
Y Right	-	-	-	-	-																																																																																
<b>Filters</b> <table border="1"> <thead> <tr> <th>Parameter</th> <th>Condition</th> <th>Value</th> <th>AND/OR</th> <th>Condition</th> <th>Value</th> </tr> </thead> <tbody> <tr> <td>ECD(µm)</td> <td>≥</td> <td>1</td> <td>AND</td> <td>≤</td> <td>100</td> </tr> </tbody> </table>		Parameter	Condition	Value	AND/OR	Condition	Value	ECD(µm)	≥	1	AND	≤	100	<b>Report Options</b> # of Chart Data pages Included: 10 Include Method Summary: <input checked="" type="checkbox"/> Include Particle Statistics and Images in report: <input checked="" type="checkbox"/>  <b>ECD (µm)</b> <table border="1"> <thead> <tr> <th></th> <th>1.00</th> <th>2.00</th> <th>5.00</th> <th>10.00</th> <th>15.00</th> <th>25.00</th> <th>40.00</th> <th>50.00</th> <th>70.00</th> </tr> </thead> <tbody> <tr> <td>Min Value</td> <td>1.00</td> <td>2.00</td> <td>5.00</td> <td>10.00</td> <td>15.00</td> <td>25.00</td> <td>40.00</td> <td>50.00</td> <td>70.00</td> </tr> <tr> <td>Max Value</td> <td>2.00</td> <td>5.00</td> <td>10.00</td> <td>15.00</td> <td>25.00</td> <td>40.00</td> <td>50.00</td> <td>70.00</td> <td>100.00</td> </tr> <tr> <td># Images</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> </tr> <tr> <td>Min Value</td> <td>1.00</td> <td>2.00</td> <td>5.00</td> <td>10.00</td> <td>15.00</td> <td>25.00</td> <td>40.00</td> <td>50.00</td> <td>3.00</td> </tr> <tr> <td>Max Value</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> <td>100.00</td> </tr> <tr> <td># Images</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> <td>10</td> </tr> </tbody> </table>			1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00	Min Value	1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00	Max Value	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00	100.00	# Images	10	10	10	10	10	10	10	10	10	Min Value	1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	3.00	Max Value	100.00	100.00	100.00	100.00	100.00	100.00	100.00	100.00	100.00	# Images	10	10	10	10	10	10	10	10	10
Parameter	Condition	Value	AND/OR	Condition	Value																																																																																
ECD(µm)	≥	1	AND	≤	100																																																																																
	1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00																																																																												
Min Value	1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00																																																																												
Max Value	2.00	5.00	10.00	15.00	25.00	40.00	50.00	70.00	100.00																																																																												
# Images	10	10	10	10	10	10	10	10	10																																																																												
Min Value	1.00	2.00	5.00	10.00	15.00	25.00	40.00	50.00	3.00																																																																												
Max Value	100.00	100.00	100.00	100.00	100.00	100.00	100.00	100.00	100.00																																																																												
# Images	10	10	10	10	10	10	10	10	10																																																																												

FIGURE 1. Analysis method for manual and automated samples at both ProteinSimple and Takeda sites. This analysis method was used with the batch protocol for modified Bot1 Protocol D. All samples were analyzed under the same conditions, with an analysis volume of 0.77 µL, and Optimize Illumination volume of 0.22 µL. Particle baseline was established with 0.2 µm-filtered DDI water.

	MANUAL OPERATION	BOT1 PROTOCOL D	MODIFIED BOT1 PROTOCOL D
Bot1 flush with sample	N/A	0.90 mL*	0.50 mL
Sample volume dispensed (defined in Method)	0.90 mL	0.70 mL	0.77 mL
Sample purge volume (defined in Method)	0.20 mL	0.00 mL	0.00 mL
Optimized Illumination volume and liquid type	0.22 mL of buffer (MFI 5200)	0.22 mL of sample (MFI 5200)	0.22 mL of sample (MFI 5200)
Dead volume	0.10 mL	0.03 mL	0.03 mL
<b>Total Sample Volume</b>	<b>1.20 mL</b>	<b>1.85 mL</b>	<b>1.52 mL</b>

\*Volume may be reduced by optimizing the batch protocol for a specific sample type. Optimize Illumination step is equivalent to a purge step, so additional purge step is not required.

TABLE 2. Method comparison of manual to modified Bot1 Protocol D. The Bot1 Protocol D was optimized further to reduce flush volume, resulting in a total sample volume of 1.52 mL for Modified Bot1 Protocol D, which was the automated protocol used in this study.

CRITERIA	DESCRIPTION
Volume	Minimum 900 $\mu\text{L}$ *
Throughput	>10 samples/day (or unattended operation)
Labware	1 or 2 mL deepwell plates
Mixing format	Aspirate and dispense steps using automated pipettor
Ambient temperature	Samples can be tested at room temperature (ambient)
Viscosity	Some viscosity, not highly viscous ( $\leq 10$ cP units)
Concentration	>50–150 mg/mL for protein formulations
MFI 5000 Series	Bot1 is used with MFI 5000 series

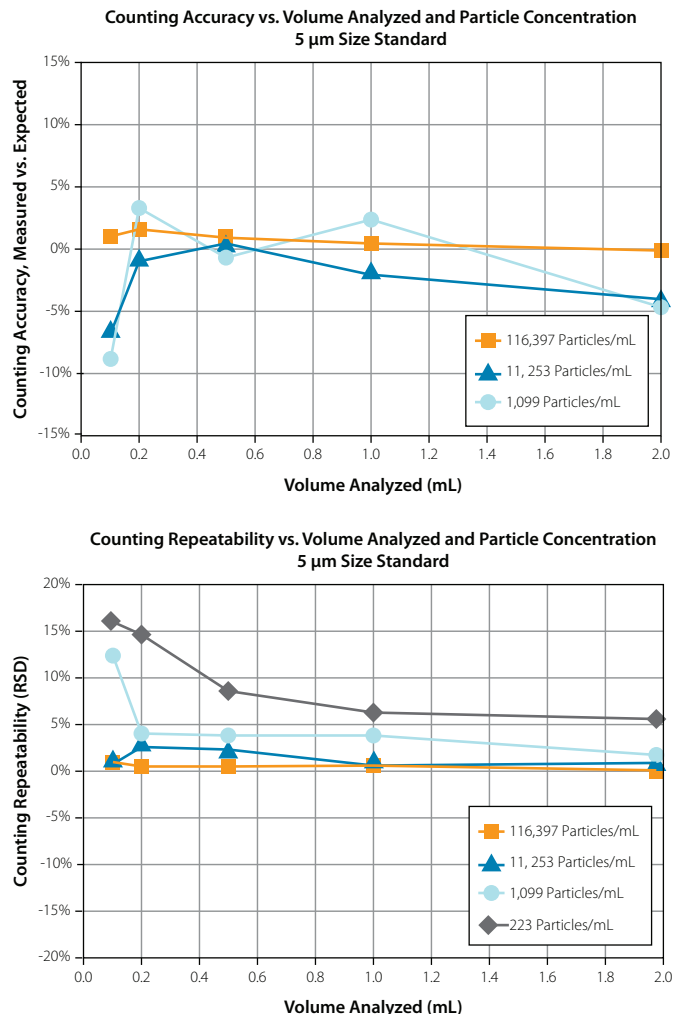
\*Required sample volume may be reduced following method optimization.

**TABLE 3.** Recommended criteria for MFI 5000 series with Bot1 autosampler.

Optimization of the automated protocol for the 1% BSA solution was performed as follows. We used a larger volume initially in Protocol D and then modified pipetting parameters as described in the modified Protocol D to reduce the volume requirements (Table 2). Additional stirring is required to counteract the effects of settling time during the automated process, so that samples are well mixed at the time of analysis. Sample types used with the MFI system in manual or automated mode should also be evaluated for suitability, as outlined in Table 3.

### GUIDELINES FOR SAMPLE VOLUME

Particle concentration in the sample can aid in determining required volume for accurate sizing and concentration measurements. The plots in Figure 2 show the impact of particle concentration on the appropriate volume. Optimal measurements are achieved for values at the 0.00% line of each graph. For example, in samples with particle concentrations of 10,000 P/mL or greater, volumes of at least 500-900  $\mu\text{L}$  should be used. Note that in samples with particle concentrations of 1,000 P/mL, volumes of 1 mL or greater should be used.



**FIGURE 2.** Impact of volume on counting accuracy for differing particle concentrations (top). Impact of volume on counting reproducibility for differing particle concentrations (bottom).

### ACCOUNTING FOR DIFFERENCES IN SAMPLE HANDLING AND VOLUME REQUIREMENTS

The automated protocol uses 1.52 mL of sample as compared to the manual protocol’s 1.20 mL. Volume and mixing steps for the automated protocol were optimized to ensure consistent counts and concentrations between replicates. As shown in Table 2, the initial total volume requirements of 1.85 mL for Protocol D were reduced to 1.52 mL in the modified version by adding automated mixing steps and reducing the sample flush volume. To reduce the total volume required, the stir speed was increased slightly from a setting of 3 to 5 (users can control the

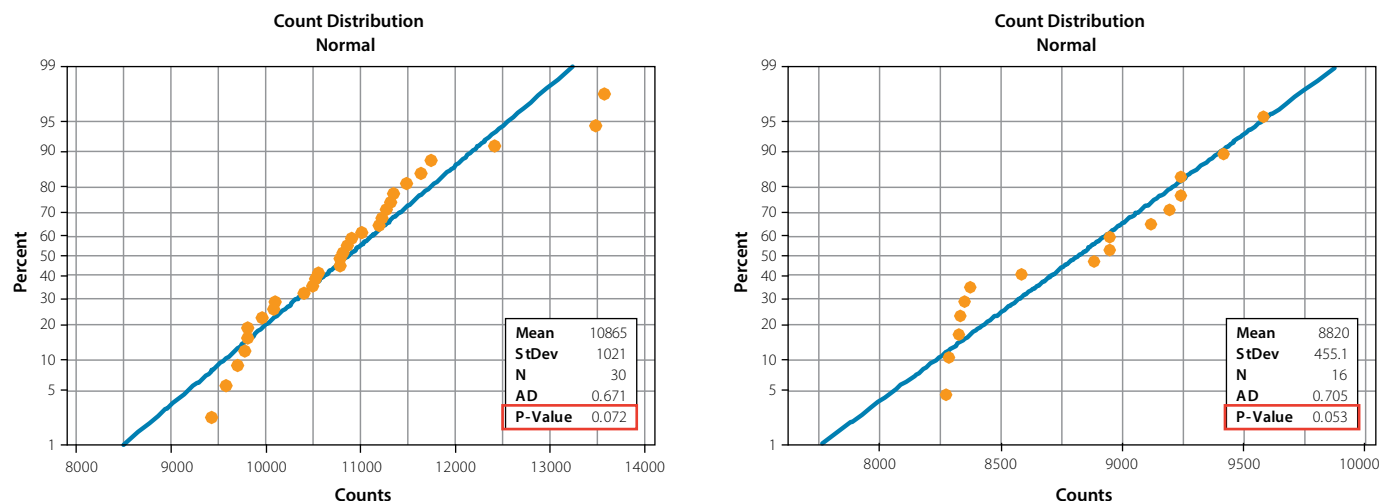


FIGURE 3. Data distribution (counts) at ProteinSimple (left) and Takeda (right). Anderson-Darling normality test was applied to the data distribution of counts. Both data sets were found to have a normal distribution, with a P-value > 0.05.

speed of this setting from 117 to 700  $\mu\text{L}/\text{sec}$ ). With this change, the actual sample flush volume required was reduced to 500  $\mu\text{L}$  from 900  $\mu\text{L}$  with no additional variation due to precipitation or aggregation. In each case, two runs of four replicates each were averaged and analyzed for count and concentration. Results from each method were analyzed using standard statistical tests as outlined in the next section.

## Results

At least eight manual and automated runs were performed at each site, using the same manual and automated protocols for pipetting and analysis (see Figure 1).

Data were analyzed to assess normality of data set, and were not found to significantly deviate from normal distribution. P-values were 0.072 and 0.053 for ProteinSimple and Takeda, respectively. Variances between automated and manual methods were not statistically different by f-test (Figure 3).

Employing a standard analysis for statistical variance (ANOVA) indicated that only between 0 and 4% of variance in particle counts can be attributed to method type (automated versus manual, Figure 4), which is not significant. Regression analysis showed no relationship between other input variables and variance in the data (Figure 5).

Comparison of data showed results at Takeda were approximately 2000 counts/mL lower than those at

One-way ANOVA: Counts versus Method

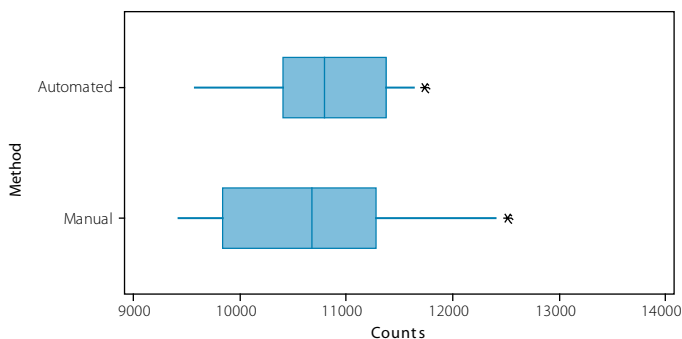
Source	DF	SS	MS	F	P
Method	1	203214	203214	0.19	0.667
Error	28	30009739	1071776		
Total	29	30212953			

S = 1035 R-Sq = 0.67% R-Sq(adj) = 0.00%

Level	N	Mean	StDev
Automated	14	10953	952
Manual	16	10788	1102

Pooled StDev = 1035

Boxplot of Counts vs. Method



One-way ANOVA: Counts versus Method

Source	DF	SS	MS	F	P
Method	1	315900	315900	1.58	0.229
Error	14	2790735	199338		
Total	15	3106636			

S = 446.5 R-Sq = 10.17% R-Sq(adj) = 3.75%

Level	N	Mean	StDev
Automated	8	8960.2	492.1
Manual	8	8679.1	395.6

Pooled StDev = 446.5

Boxplot of Counts vs. Method

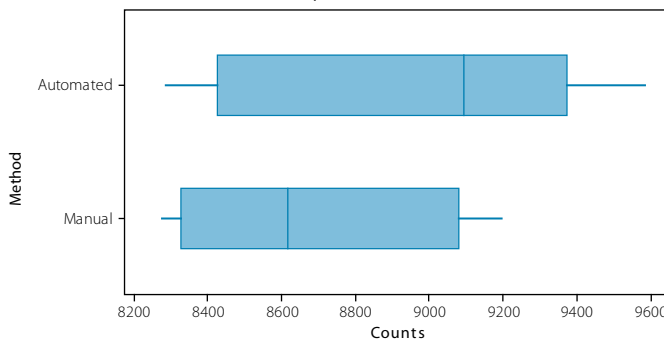


FIGURE 4. Impact of method type on counts at ProteinSimple (top) and Takeda (bottom). Analysis using one-way ANOVA did not show a significant effect for method type on particle counts, with an R-Sq(adj) of 0.0% for ProteinSimple and an R-Sq(adj) of 3.75% for Takeda.

Predictor	Coef	SE Coef	T	P
Constant	10534.6	519.9	20.26	0.000
Automated	172.3	410.2	0.42	0.678
Set 2	279.3	404.0	0.69	0.496
Day 2	224.6	410.2	0.55	0.589
Run 2	107.0	553.1	0.19	0.848
Run 3	-266.9	553.1	-0.48	0.634
Run 4	164.7	605.9	0.27	0.788

S = 1106.28 R-Sq = 6.8% R-Sq(adj) = 0.0%

PRESS = 48120295 R-Sq(pred) = 0.00%

Analysis of Variance

Source	DF	SS	MS	F	P
Regression	6	2064274	344046	0.28	0.940
Residual Error	23	28148679	1223856		
Total	29	30212953			

Source	DF	Seq SS
Automated	1	203214
Set 2	1	585203
Day 2	1	463322
Run 2	1	153389
Run 3	1	568725
Run 4	1	90420

Predictor	Coef	SE Coef	T	P
Constant	8885.2	249.4	35.62	0.000
Automated	398.8	245.6	1.62	0.136
Set 2	41.4	229.8	0.18	0.861
Run 2	-162.4	313.1	-0.52	0.615
Run 3	-590.6	319.1	-1.85	0.094
Run 4	-189.8	319.1	-0.59	0.565

S = 442.818 R-Sq = 36.9% R-Sq(adj) = 5.3%

PRESS = 5405327 R-Sq(pred) = 0.00%

Analysis of Variance

Source	DF	SS	MS	F	P
Regression	5	1145756	229151	1.17	0.388
Residual Error	10	1960880	196088		
Total	15	3106636			

Source	DF	Seq SS
Automated	1	401933
Set 2	1	15025
Run 2	1	21197
Run 3	1	638197
Run 4	1	69403

FIGURE 5. Linear regression analysis was used with dummy coding to evaluate relevant variables (manual method, automated method, day) as a group. Analysis shows no significant impact of system inputs (automation, day to day variance) on particle counts. The various system inputs could account for 0.0% (ProteinSimple, left) and 5.3% (Takeda, right) of the variation in the data. An R-Sq (adj) of >30% would indicate a significant impact.

MANUAL			AUTOMATED		
RUN	COUNTS	COUNTS/ML	RUN	COUNTS	COUNTS/ML
1	8783	11184	1	8463	10780
2	9745	12409	2	8444	10773
3	8530	10862	3	8483	10812
4	8238	10490	4	9003	11479
<b>Average</b>	8824.00	11236.25	<b>Average</b>	8598.25	10961.00
<b>Std Dev</b>	565.63	720.25	<b>Std Dev</b>	234.09	299.43
<b>%CV</b>	6.41%	6.41%	<b>%CV</b>	2.72%	2.73%
RUN	COUNTS	COUNTS/ML	RUN	COUNTS	COUNTS/ML
5	7918	10083	5	8840	11274
6	7703	9809	6	9124	11633
7	7815	9951	7	8905	11343
8	7403	9427	8	10572	13471
<b>Average</b>	7709.75	9817.50	<b>Average</b>	9360.25	11930.25
<b>Std Dev</b>	192.73	245.40	<b>Std Dev</b>	707.47	899.69
<b>%CV</b>	2.50%	2.50%	<b>%CV</b>	7.56%	7.54%
AVERAGE OF ALL 8 MEASUREMENTS			AVERAGE OF ALL 8 MEASUREMENTS		
<b>Average</b>	8227.83	10477.17	<b>Average</b>	9058.00	11545.67
<b>Std Dev</b>	875.47	1114.73	<b>Std Dev</b>	787.21	1001.27
<b>%CV</b>	10.64%	10.64%	<b>%CV</b>	8.69%	8.67%

TABLE 4. Count and concentration for automated versus manual format for 1% BSA, ProteinSimple.

ProteinSimple. This difference is likely related to sample quality in terms of sample stress or degradation during shipment. Neither site showed any significant difference in counts or concentration between the manual and automated method, confirming that similar results can be obtained using either format.

#### COMPARISON OF THE MANUAL VERSUS AUTOMATED METHOD, PROTEINSIMPLE

Data from tests performed at ProteinSimple are listed in **Table 4**. The automated format of 1% BSA protein solution produced consistent counts and concentration data, compared to the manual format. The modification to Protocol D reduces the sample flush volume without impacting the consistency of

results between replicates in the automated method. The standard deviation for both methods was 5% or less for all samples tested.

#### COMPARISON OF THE MANUAL VERSUS AUTOMATED METHOD, TAKEDA

Data from tests performed at Takeda are listed in **Table 5**. At the Takeda site, the automated format of 1% BSA protein solution produced consistent counts and concentration data which were also comparable to the manual format. The standard deviation for both methods was 6% or less for all samples tested. Takeda results were similar to ProteinSimple's in terms of successful transfer of the 1% BSA assay to an automated format.

MANUAL			AUTOMATED		
RUN	COUNTS	COUNTS/ML	RUN	COUNTS	COUNTS/ML
1	6570	8947	1	6641	8950
2	6833	9197	2	6477	8586
3	6263	8275	3	6093	8288
4	6344	8333	4	6965	9418
<b>Average</b>	6502.52	8688.06	<b>Average</b>	6544.21	8810.20
<b>Std Dev</b>	221.60	394.50	<b>Std Dev</b>	313.97	421.88
<b>%CV</b>	3.41%	4.54%	<b>%CV</b>	4.80%	4.79%
RUN	COUNTS	COUNTS/ML	RUN	COUNTS	COUNTS/ML
5	6748	8884	5	7043	9241
6	6198	8349	6	7043	9241
7	6670	9123	7	6261	8373
8	6235	8326	8	7373	9585
<b>Average</b>	6462.72	8670.23	<b>Average</b>	6930.14	9110.15
<b>Std Dev</b>	248.28	343.69	<b>Std Dev</b>	409.09	448.12
<b>%CV</b>	3.84%	3.96%	<b>%CV</b>	5.90%	4.92%
AVERAGE OF ALL 8 MEASUREMENTS			AVERAGE OF ALL 8 MEASUREMENTS		
<b>Average</b>	6542.31	8804.16	<b>Average</b>	6806.43	8996.05
<b>Std Dev</b>	267.33	379.17	<b>Std Dev</b>	416.26	452.95
<b>%CV</b>	4.09%	4.31%	<b>%CV</b>	6.12%	5.03%

TABLE 5. Count and concentration for automated versus manual format for 1% BSA, Takeda.

## Conclusion

The MFI 5200 produces the same high quality particle characterization in manual or automated mode. This study showed that method optimization could further reduce the sample volume required without impairing concentration accuracy. Statistical analysis confirmed that these protocols are robust and provide an example of standardization of methods across instrument configurations.

In the Takeda development laboratory, MFI was originally implemented due to its advantages over conventional techniques (HIAC-based light obscuration used according to Ph. Eur. 2.9.20/USP <788>) for measurement of sub-visible particles in the  $\mu\text{m}$  range (e.g. 2–10  $\mu\text{m}$ ).

Those advantages included:

- Lower sample consumption versus light obscuration
- 1–10  $\mu\text{m}$  particle detection
- MFI's ability to distinguish between particle types based on morphology parameters

Initially, a manual MFI 4200 system was obtained for the Takeda development work which included formulation development, primary container selection, stability studies, etc. While the advantages over HIAC-based light obscuration were obvious in terms of volume range, particle size and morphology, manual MFI use became labor-intensive and repetitive with larger development

studies containing up to 50 samples per run. Reduction of hands-on time associated with manual operation was thus a primary goal, and became the main reason for upgrading to the automated MFI 5200 system with Bot1 autosampler at an early stage. In addition, the improved reproducibility associated with the automated format further contributed to the decision to upgrade.

The option to automate provides a key advantage for particle characterization of protein formulations, offering many benefits compared to more common techniques. Automated protocols allow for much greater throughput and less hands-on time, with up to 80 samples per unattended run. Implementation of automation for higher throughput and standardization can help address demand for more rapid and consistent screening methods in particle characterization.

## References

1. Micro-flow Imaging: flow microscopy applied to sub-visible particulate analysis in protein formulations, D Sharma, D King, P Oma and C Merchant, *AAPS Journal*, 2010; 12(3).
2. In-Process Revision: <787> subvisible particulate matter in therapeutic protein injections (USP36-NF31 IS), February 2013.
3. AAPS update on USP expert committee for subvisible particle analysis, L Narhi, *Newsletter of the AAPS Aggregation and Biological Relevance Focus Group*, 2012; 3(2).



Toll-free: (888) 607-9692  
Tel: (408) 510-5500  
Fax: (408) 510-5599  
orders@proteinsimple.com  
proteinsimple.com