

<b>Quanterix</b>	<b>Form</b>	
	<b>Document No:</b> FRM-0081	<b>Revision No:</b> 03
	<b>Name:</b> Homebrew Bead Filling Record	<b>Effective:</b> 07 May 2018 <b>Page 1 of 5</b>

**1. BATCH INFORMATION**

Component name	Homebrew Helper Bead (779) 1.5g
Item number	103208
Lot Number (Vialled Reagent)	825608
Lot Number (Source Material/ Concentrate lot)	825605
Expiration date (DDMMYYYY)	13 Mar 2019
Bead Concentration (beads/ml)	$1.219 \times 10^9$
Quantity (# of units to be filled)	16
Vial / bottle size	4mL
Fill Volume Target	1.8g
Filling date (DDMMYYYY)	24 Sep 2018
Filling operator	BJ <sup>2</sup>
Checked by	Rajina Manandhar

**2. PURPOSE**

The form is used to record the required information for Bead filling operations. This document is used alongside Formulation and Label documentation to record filling for better traceability.

<b>Quanterix</b>	<b>Form</b>	
	<b>Document No:</b> FRM-0081	<b>Revision No:</b> 03
	<b>Name:</b> Homebrew Bead Filling Record	<b>Effective:</b> 07 May 2018 <b>Page 2 of 5</b>

**3. REFERENCES**

Document Number	Document title
FRM-0113	Line Clearance Checklist
SOP-0028	Quanterix Glossary
SOP-0033	Balance Operation
SOP-0034	Gowning
SOP-0039	WandMixer
SOP-0040	Cleaning Labware
SOP-0042	Zebra Label Printer Operation (ZM400)
SOP-0043	Fill Volume Information for Simoa Components
SOP-0049	Manual Filling Operation
PROC-0004	Quanterix Chemical Hygiene Plan
PROC-0003	Quanterix Biosafety Manual

**4. SAFETY AND ENVIRONMENTAL INFORMATION****4.1. Waste Handling and Disposal**

- 4.1.1 Solutions containing chemical and biological substances must be decontaminated according to the Chemical Hygiene Plan and Biosafety Manual prior to disposal.

**5. EQUIPMENT**

Document Title	Equipment #
Pipetting equipment	PS41736
Weighing equipment	28804486

**6. MATERIALS**

Description	Item #
Vial, 2 mL, Clear, PP, Sterile	100442
Closure, 2 mL Clear HDPE, Sterile	100443
Bottle, 4 mL, Amber, PP	100444
Closure, 4mL, White, PP	103086
Bottle and Cap, 8 mL, Clear	101394

<b>Quanterix</b>	<b>Form</b>	
	<b>Document No:</b> FRM-0081	<b>Revision No:</b> 03
	<b>Name:</b> Homebrew Bead Filling Record	<b>Effective:</b> 07 May 2018
		<b>Page</b> 3 of 5

## 7. PERFORMING THE FILL

- Refer to SOP-0049 for the manual filling process.

### 7.1. Fill Bulk Solution into Reagent Bottles

7.1.1 Print labels then adhere to the required number of reagent bottles.

7.1.2 Fill the bulk into the labeled reagent bottles then cap. Verify the fill weight throughout the filling event as specified below.

- If the total number of bottles is less than 300, identify 3 bottles for weight verification check. If the total number of bottles is greater than 300, use 5 bottles for weight verification. The bottles should be from the beginning, middle, and end of the filling episode.
- Record the tare weight of each bottle with a cap and record in the table below.
- Record serial number of the bottle used for weight check if applicable, if not then identify the bottle with a number.
- Once a weight verification bottle is filled, stop and weigh the sample to ensure it is in the specified range.
- If bottle passes specified range continue the filling process. If bottle fails the specified range, stop the filling process and contact immediate supervisor. Document course of action in the comment section below.

Released

<h1>Quanterix</h1>	<b>Form</b>			
	Document No: FRM-0081		Revision No: 03	
	Name: Homebrew Bead Filling Record		Effective: 07 May 2018 Page 4 of 5	

<input type="checkbox"/> N/A	4-mL Amber Vial				
Item: Dye Encoded Homebrew Target: 1.8g - 2.0g	Beg. Bottle <u>1</u>	Middle Bottle 1 <u>8</u>	Middle Bottle 2 <u>    </u>	Middle Bottle 3 <u>    </u>	End Bottle <u>16</u>
Gross weight	4.00	3.99			4.02
Tare weight	2.19	2.19			2.19
Net weight	1.81	1.80			1.85 <small>BF 24 Sep</small> 1.83
Acceptance	<input checked="" type="checkbox"/> Pass <input type="checkbox"/> Fail	<input checked="" type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input checked="" type="checkbox"/> N/A	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input type="checkbox"/> N/A	<input checked="" type="checkbox"/> Pass <input type="checkbox"/> Fail

N/A Comment:

---



---



---

<input checked="" type="checkbox"/> N/A	8-mL Clear Vial				
Item: 103207 Target: 5.0 - 5.2g	Beg. Bottle <u>    </u>	Middle Bottle 1 <u>    </u>	Middle Bottle 2 <u>    </u>	Middle Bottle 3 <u>    </u>	End Bottle <u>    </u>
Gross weight					
Tare weight					
Net weight					
Acceptance	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input type="checkbox"/> N/A	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input type="checkbox"/> N/A	<input type="checkbox"/> Pass <input type="checkbox"/> Fail

N/A Comment:

---



---



---

Quanterix	Form	
	Document No: FRM-0081	Revision No: 03
	Name: Homebrew Bead Filling Record	Effective: 07 May 2018 Page 5 of 5

<input checked="" type="checkbox"/> N/A	2-mL Clear Vial				
Item: 103206	Beg. Bottle	Middle Bottle 1	Middle Bottle 2	Middle Bottle 3	End Bottle
Target: 1.5g - 1.7g	_____	_____	_____	_____	_____
Gross weight					
Tare weight					
Net weight					
Acceptance	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input type="checkbox"/> N/A	<input type="checkbox"/> Pass <input type="checkbox"/> Fail <input type="checkbox"/> N/A	<input type="checkbox"/> Pass <input type="checkbox"/> Fail

N/A Comment:

---



---



---

7.1.3 Document number of vials filled, QC samples, and amount for inventory.

Vials filled	QC Samples	Inventory (Filled-QC)
16	0	16

7.1.4 Store filled reagent bottles at 2–8°C.

**8. DOCUMENTATION/RECORDS MANAGEMENT**

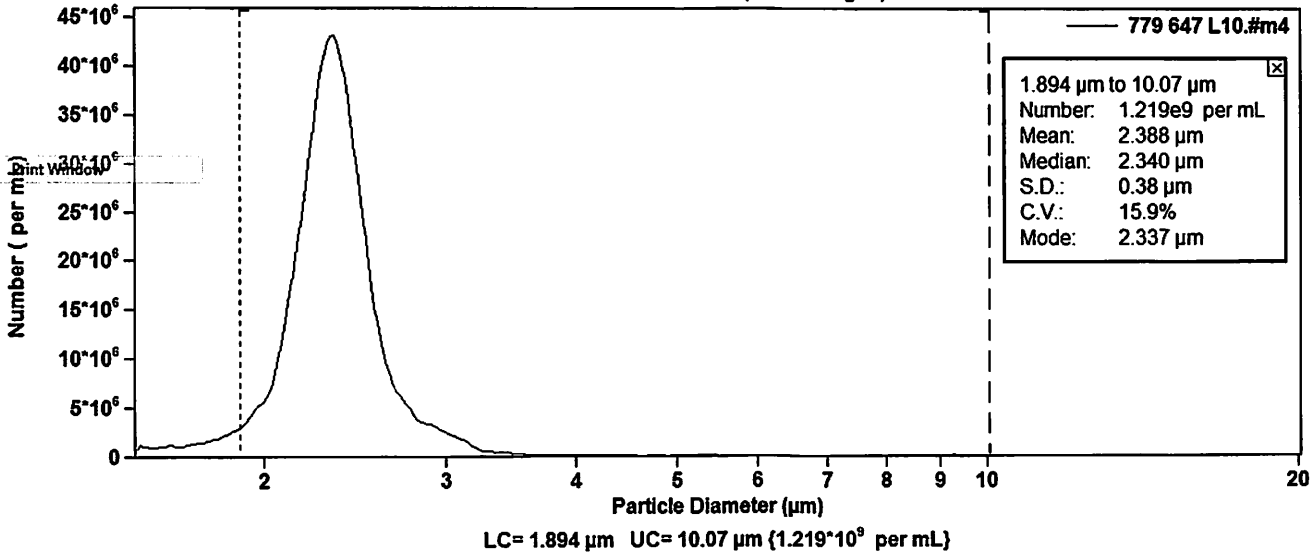
8.1. Completed records must be submitted to Quality Control for review and archival.

**9. REVISION HISTORY**

Revision	Detailed Description	Date	Originator
1.0	Initial Release	07Jan2016	B. Flaherty
02	DCR-16-1150: The fill range for part # 100451 was changed from "1.8 – 2.0 g" to "1.5-1.7g" to accurately reflect the fill range of this product.	18Jan2017	R. Fregeolle
03	DCR-18-0474: Update to new HB part numbers.	07May2018	B. Pink

*End of Document*

Differential Number (Smoothing=7)



BTM sp5

**1. Component or Kit Information**

Component Name:	Honeybrow Helper Bead 779 1.5mL
Part Number:	105208
Lot Number:	825608
Expiration Date:	13 Mar 2019
Storage Temperature:	2-8°C
Number of labels requested:	18
Completed By:	BS
Date:	24 Sep 2018
Verified by:	RM
Date:	24 Sep 2018

# of Labels Printed	Printed by	Date
18	RM	24 Sep 2018

**2. Label Inspection, completed by Verifier**

Print Quality	Item #	Lot #	Expiration Date	Storage Temp.	Barcode Check & Scan
<input checked="" type="checkbox"/> Pass	<input checked="" type="checkbox"/> Pass	<input checked="" type="checkbox"/> Pass	<input checked="" type="checkbox"/> Pass	<input checked="" type="checkbox"/> Pass	<input type="checkbox"/> Pass
<input type="checkbox"/> Fail	<input type="checkbox"/> Fail	<input type="checkbox"/> Fail	<input type="checkbox"/> Fail	<input type="checkbox"/> Fail	<input type="checkbox"/> Fail
<input type="checkbox"/> N/A	<input type="checkbox"/> N/A	<input type="checkbox"/> N/A	<input type="checkbox"/> N/A	<input type="checkbox"/> N/A	<input checked="" type="checkbox"/> N/A

Labels Verified by: BS

Date: 24 Sep 2018

Homebrew Helper Bead Vial (779) 1.5ml

REF 103208 2\*18° Conc: 1.22 x10<sup>8</sup> beads/ml  
 LOT 825608  
 13-Mar-2019 Quanterix Corp  
 For Research Use Only LBT-181-220 Rev 1.0

*BF 24 Sep 18*

Verifier attach label above, initial and date label

3. Labeling Line Clearance (✓ each box)

Area clear of unrelated material	Material lot #, paperwork & labels match	Work area clean and clear of debris
✓	✓	✓

Performed by	Date
<i>BF</i>	<i>24 Sep 18</i>

4. Label Accountability

- A. # of Bottles/Kits Labeled: 16
  - B. # of Labels on Form: 1
  - C. Total # of Labels (A+B): 17
  - D. # of Labels Requested: 18
  - E. Difference (D-C): 18
  - F. # of Labels Destroyed: 1
  - G. Reconciliation (F-E): XC
- BF 24 Sep 18*





<h1>Quanterix</h1>	<b>Form</b>	
	<b>Document No:</b> FRM-0108	Revision No: 05 Effective: 20 Mar 2018 Page 4 of 5
	<b>Name:</b> Component Labeling and Filling Record	

7.  **N/A Filling Bulk Solution into Reagent Bottles**

Fill the bulk into the labeled reagent bottles then cap. Verify the fill weight throughout the filling event as specified below.

- If the total number of bottles is less than 300, identify 3 bottles for weight verification check. If the total number of bottles is greater than 300, use 5 bottles for weight verification. The bottles should be from the beginning, middle, and end of the filling episode.
- Record the tare weight of each bottle with a cap and record in the table below.
- Record serial number of the bottle used for weight check if applicable, if not then identify the bottle with a number.
- Once a weight verification bottle is filled, stop and weigh the sample to ensure it is in the specified range.
- If bottle passes specified range continue the filling process. If bottle fails the specified range, stop the filling process and contact immediate supervisor. Document course of action in the comment section below.
- To convert target fill and target range into grams multiply by 1.02.

Vial/Bottle Size	Target Fill	Target Range	Filling Equipment

	Begin Bottle	Middle Bottle 1	<input type="checkbox"/> N/A Middle Bottle 2	<input type="checkbox"/> N/A Middle Bottle 3	End Bottle
Gross weight					
Tare Weight					
Net weight					
Acceptance	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail	<input type="checkbox"/> Pass <input type="checkbox"/> Fail
Performed by:	Date:		Verified by:		Date:

Verification Performed by (Mfg): *Ben Sy*  
*ljk* 25 Sep 2018  
 Packet Reviewed by (QC): *Imah Cam*

Date: *24 Sep 2018*  
 Date: *25 Sep 2018*

**REVISION HISTORY**

Revision	Detailed Description	Date	Originator
1.0	Initial Release. Created new doc number FRM-100-0035 that replaces FRM.009. (Doc number follows convention FRM-NNN-NNNN.)	30Oct2015	K. Lerma
1.1	Add specific gravity calculation to convert from mL to grams. Formatting.	18Dec2015	K. Lerma
1.2	DCR-16-0598: Transfer documents from QMS 2.0 to R&D vault. Archive the document in QMS 2.0 once released in R&D.	28Sep2016	S. Chin
03	DCR-16-1219: Update header and footer format from M-Files to MasterControl	16Nov2016	S. Moriarty
04	DCR-18-0218: Add Line Clearance to improve the accuracy of the kitting process.	08Feb2018	D. Ahuja
05	DCR-18-0363: Delete step header, unnecessary header causing issues with use of n/a box.	16Mar2018	B. Flaherty

*End of Document*