Thermo Fisher S C I E N T I F I C

2023 ISBER, PJ-01

Investigating the labeling workflow and incorporation of automation for increased collaboration

Amy Armstrong¹, Wilfredo Marin², Elana S. Novotny³, Thermo Fisher Scientific

1. 168 Third Ave, Waltham, MA 02451 2. 275 Aiken Road, Asheville, NC 28804 3. 75 Panorama Creek Dr, Rochester, NY 14625

Introduction

Biobanks have long since served as foundational facilities, storing varying amounts of biomaterial and data for translational research efforts. In order to improve biobanking operations, innovation is required in how sample information is captured to facilitate sample identification, viability, and sharing with collaborators. It may seem easiest to continue handwritten labeling methods, but the possibility of mistakes that render a sample useless is too high of a risk. Here we test a protocol that includes a barcode scanner intended to increase speed, security, and support of persistent tagging for the traceability of biological material and associated

Results

Handwriting

Figure 2. Statistics of the handwriting study. This handwriting test was imperfect because the same person could get the same tube more than once. However, since the scientists were told speed was important, the information was likely not memorized between rounds.



Automation

Figure 5. The interface of VisionMate HSX software with the different parameters that can be optimized for convenient export of data. The VisionMate HSX High Speed Barcode Reader comes with software for exporting data from tubes into a workable format, such as with Excel software. While straightforward, a 15 min training session was arranged to make sure users understood the minutiae of how to export plate data into Excel software.



Conclusions

Scaling up

Figure 8. Comparing handwriting versus automation for data entry. Writing information by hand on tubes requires more time, on average, per tube than using automation (data entry of information into Excel software in this case). Actual reading of the tubes using automation is fast (~1 sec), though there is some initial time needed to install, learn the software, and set up how exports will occur. The potential for time savings and fewer mistakes is apparent when looking at scaling up to a realistic number of tubes, such as 30,000 per day.

data in a biobank [1].

Once implemented as a standard practice, biobankers can maintain confidence through the chain of custody that accompanies many samples. Incorporating sample information into a laboratory information management system (LIMS) helps increase the value of the well-organized biomaterial without burdening the vial with potentially identifiable information. With an initial investment and potential plans to move towards automation, most facilities can recover these costs in a short period from the ability to handle higher volumes of samples, increasing capacity at scale [2].

Materials

Tubes: Thermo Scientific[™] Matrix[™] ScrewTop Tri-coded Tube (Cat. No. 3741-WP1D-BR)



Automation reader: Thermo Scientific[™] VisionMate[™] HSX High Speed Barcode Reader

- 7.9 in. x 5.9 in. x 6.3 in. (L x W x H); 6 lb
- Reading area: 5 in. x 3.39 in.



Figure 3. Comparing the writing on the tubes. The information in red is where errors were introduced during the rounds. In general, no one specific type of data was misrepresented more than others.

0218973	0975283	7182004	6900128	0048917
PL2	URN	CRD	SAL	TIS
0001151	0001151	0002837	0007735	0001151
2/24/2001	6 /30/2002	1/17 /2019	7/11/ 1995	6/25/1984
12/14/2022	11/28/2022	11/29/2022	12/18/2022	12/19/2022
3980117	1153277	3399557	5170228	5098270
AMN	RBC	TER	U24	CLN
0002837	0001151	0007735	0003278	0002837
4/19/1992	11/23/1970	10/23/1960	8/30/1972	5/13/1980
11/11/2022	11/21/2022	10/31/2022	10/27/2022	11/30/2022

Figure 4. Example tubes showing mistakes, and comments made by scientists during the study.



"Holding onto the cap was more comfortable and made it easier to write on the tube, but being left-handed meant the orientation of the label is different."

-Scientist 1



Figure 6. A successful (top) and unsuccessful (bottom) read of the 96-tube rack of the Matrix tubes. These images illustrate some of the capabilities of the automation software, such as alerting the user if the number of samples is not as expected.











Automation can save biobanks time and lead to less



Procedure

Scientists were provided with 5 simulated samples of information to mimic the last stage of labeling a tube for storage in a biobank (Figure 1). Their job was to copy all information onto the tube while being as fast and as accurate as possible. All 15 tubes were then mixed, and each scientist took 5. They had to recopy the information onto new tubes, to simulate transferring a sample. This process was repeated once more. The tubes were then analyzed for accuracy.

To compare to automation, these data were aligned with 15 tubes using Microsoft[™] Excel[™] software and the VisionMate HSX High Speed Barcode Reader. To better simulate processes at a biobank, the data were typed into Excel software.

Figure 1. The data provided for marking the tubes in this exercise. Each tube was marked with (1) sample ID, (2) aliquot sample type [3], (3) clinic ID, (4) collection date, and (5) date of birth. The data were all made-up but generated based on a realistic system



Observation: Tubes that were written by a left-handed person (holding the cap with the right hand) were more convenient for a right-handed person to pick out of the rack and read. However, the two left-handed people in this study did not write in the same orientation. These differences should be considered at each specific biobank.

"The tubes roll so it is hard to hold them and transfer information from one to another."

–Scientist 2



"One of the things I was worried about when writing down the information was smearing the ink." -Scientist 3



Figure 7. Examples of data entry using Excel software. Depending on how the database is set up, much more information can be incorporated with each tube using automation and Excel software. Although one mistake occurred copying the data into Excel software, it was caught and corrected before moving on, because of having more time.

Observation: Specific biobanks will dictate which LIMS is best for their

purposes and what information they need recorded with respect to each

sample. The capabilities of the VisionMate reader enable it to be flexible

and meet many different needs.

	0		6						
A B Tube 3540907564				B 3540907564			(BEMT): Online financial pla		
		2	Sample ID		7182004			sustainability <i>Biopreserva</i>	
Sample ID 7182004		3	Aliquot Sample Type		CRD			odotamaomity	Dioprocorvatio
Aliquot Sample Ty	4	Clinic ID		0002837			3 Batsou E at al (2018) Stand		
Clinic ID	0002837	5	Collection Date		1/17/2019				
Collection Date 1/17/2019		6	Date of Birth		11/29,	11/29/2022		3.0. Biopreservation and Bio	
Date of Birth	11/29/2022	7		unte Tune Definition	Candhla				
		8	Allquot Sar	hple Type Definition					
		10			AAAAAA	AAAA			edaments
	11	Additional Samples Tube Numbers		BBBBBBBBBB			/		
		12							a thank Carina F
		13	-					vve would like t	o inank Sanna E
		14	Addition	al Sample Information:	YYYYYY	Y V		participation du	ring the handwri
		15			111111	1			C
	A		В	С			D	E	F
9	Tube	35409	07564	3540909779	9	354	0907557	3540907556	3540909771
Sample ID		718	2004	0975283		2381904		1039273	0218973
Aliquot Sample Type		CRD		URN	BON		BON	BMA	PL2
	Clinic ID	000	2837	0001151		00	07735	0009889	0001151
Collection Date		1/17	/2019 6/30/2002		e.	10/15/2005		1/10/1944	2/24/2001
Date of Birth		11/29	/2022	2022 11/28/2022		12/15/2022		12/15/2022	12/14/2022
Aliquot Sample Type Definition		Cord blood		Urine, random ("spot")		Bone		Bone marrow aspirate	Plasma, double spun
Additional Samples Tube		XXXXXXXXXX		22222222		GGGG	GGGGGGG	НННННННН	VVVVVVVV
		ΑΑΑΑΑΑΑΑΑ		DDDDDDDDD				PPPPPPPPP	ККККККККК
		BBBBBBBBBB		ZZZZZZZZZZ					тттттттт
				LLLLLLLL					
3									
Additional Sample Information:		YYYYYYY YYYYYYYY YYYYYYYY		YYYYYYYY YYYYYYYY		YYYYYYY		ΥΥΥΥΥΥΥ	YYYYYYY
									ΥΥΥΥΥΥΥ
5									

- erroneous data. Automation can be more scalable.
- Incorporation of automation can help position biobanks for collaborations.

References

- 1. International Organization for Standardization (2018) Biotechnology—Biobanking—General requirements for biobanking (ISO 20387:2018). https://www.iso.org/standard/67888.html.
- 2. Odeh H et al. (2015) The Biobank Economic Modeling Tool Online financial planning to facilitate biobank ability. Biopreservation and biobanking 13(6):421–429.
- F et al. (2018) Standard PREanalytical Code Version preservation and Biobanking 16(1):9–12.

wledgments

like to thank Sarina Bellows and Colin Wright for their on during the handwriting trial.

realistic system.			different viewing angles
Scientist 1	Scientist 2	Scientist 3	
0218973	6900128	0000197	
PL2	SAL	BLD	
0001151	0007735	0002837	
2/24/2001	7/11/1995	12/8/1956	
12/14/2022	12/18/2022	11/15/2022	
1039273	0048917	1153277	
BMA	TIS	RBC	
0009889	0001151	0001151	
1/10/1994	6/25/1984	11/23/1970	
12/15/2022	12/19/2022	11/21/2022	
2381904	5098270	3399557	"Loo
BON	CLN	TER	
0007735	0002837	0007735	
10/15/2005	5/13/1980	10/23/1960	
12/15/2022	11/30/2022	10/31/2022	
0975283	3980117	2270019	
URN	AMN	CEL	
0001151	0002837	0003278	
6/30/2002	4/19/1992	9/11/1945	
11/28/2022	11/11/2022	10/27/2022	
7182004 CRD 0002837 1/17/2019 11/29/2022	4078291 ZZZ 0003278 3/14/1999 11/15/2022	5170228 U24 0003278 8/30/1972 10/27/2022	Observ during t easier a would lil

'Looking through to copy the data from one tube to another was hard.' –Scientist 3

> "My neck feels strained even after only copying down 5 tubes." -Scientist 1

servation: The data were provided with dates using a slash (/), but ng the experiment the scientists found that using a dash (-) was ier and was no more likely to be misinterpreted as a slash. This Id likely be a convenience set by the specific biobank.



Science at a scan Scan the QR code on the right with your mobile device to download this and many more scientific posters.



*

Learn more at thermofisher.com/biobanking

This product is intended for General Laboratory Use. It is the customer's responsibility to ensure that the performance of the product is suitable for customer's specific use or application. © 2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Microsoft and Excel are trademarks of Microsoft Corporation. COL27646 0423