

## Clinical centrifuges

# Using Thermo Scientific compact centrifuges for the effective and efficient processing of clinical samples

## Authors

Richard Sicard, Thermo Fisher Scientific, Asheville, North Carolina, USA

Romana Hinz, Senior Global Product Manager for Compact and Micro Centrifuges at Thermo Fisher Scientific

## Keywords

Compact centrifuge, Platelet-poor plasma (PPP), Urinalysis, Blood chemistry analysis

## Introduction

Clinical laboratory testing plays an important role in the detection, diagnosis, and treatment of disease. The use of centrifugal force for separation of blood and urine is a very crucial step within the process. It is very important not only to perform this step correctly, but also efficiently so as to derive the best results for the patient sample.

Blood is typically separated to yield platelet-poor plasma (PPP), which in turn is used for most coagulation tests. Ideally, PPP should have a platelet count of less than  $10 \times 10^3$  per  $\mu\text{L}$ <sup>1</sup>.

Similarly, the analysis of urine is also very important in the clinical laboratory. Information derived from this testing can predict many important medical concerns, one of which is the function of kidneys or associated infections. This common test measures various compounds such as white blood cells, red blood cells, epithelial cells, bacteria, calcium oxalate crystals, mucous threads, yeast, and hyaline casts that pass through the urine<sup>2,3</sup>.

Finally, blood chemistry analysis is also a very common test in the clinical laboratory. These results are used to evaluate a variety of components such as potassium, aspartate aminotransferase (AST), phosphorus, and lactate dehydrogenase (LDH)<sup>4</sup>.

The objective of this note is to illustrate the efficiency and effectiveness of Thermo Scientific™ small benchtop centrifuges in critical centrifugation steps of the following clinical testing: (A) platelet-poor plasma (PPP) preparation, (B) urinalysis, and (C) blood chemistry analysis.

Thermo Scientific™ C Series Centrifuges (see Figure 1) can be used with the Thermo Scientific™ TX-150 swinging bucket rotor for high-speed horizontal processing. The TX-150 rotor accommodates up to 24 x 5/7 mL blood collection tubes or 8 x 50 mL conical tubes at 4,500 rpm (3,260 x g) but when configured with the Thermo Scientific™ HIGHConic™ rotor, it can achieve up to 8,700 rpm (10,155 x g) with 6 tubes. While a lower capacity than the standard swinging bucket rotor, this rotor's higher speed does make it a very important rotor in efficiently and effectively processing PPP.

These small benchtop centrifuges offer sample protection with the Thermo Scientific™ ClickSeal™ biocontainment lid system certified by Public Health England, Porton Down, UK. They also feature Thermo Scientific™ Auto-Lock™ Rotor Exchange which allows rotors to be installed and removed from the chamber of the centrifuge with the push of a button—no tools required.



Figure 1. Thermo Scientific™ C1T, C1TR and C1FR Centrifuges

## Methods

Platelet-poor plasma (PPP) preparation was done by performing centrifugation of whole blood using 2.7 mL BD Vacutainer™ Plus plastic citrate tubes. The initial platelet concentrations from whole blood as well as for each PPP preparation were determined with the aid of a hematology analyzer.

Urinalysis is composed of two main steps. In the first step urine was centrifuged using 12 mL KOVA™ tubes in the TX-150 swinging bucket rotor to separate sediment from urine which was then examined in the second step for crystals, casts, cells, yeast, and bacteria under a microscope.

Blood chemistry analysis was performed by collecting blood into 7 mL BD Vacutainer SST Plus tubes, followed by centrifugation (TX-150 and HIGHConic rotors). The serum was tested for potassium, aspartate aminotransferase (AST), phosphorus, and lactate dehydrogenase (LDH) using a chemistry analyzer.

## Results

Clinically acceptable results were derived for all centrifugation steps performed using a Thermo Scientific small benchtop centrifuge with the TX-150 swinging bucket rotor or HIGHConic fixed angle rotor. Acceptable results for PPP were accomplished at 4,500 rpm for 10 minutes using the TX-150 rotor and 1 minute at 8,700 rpm using the HIGHConic rotor. Similarly, it was at 1,700 rpm for 4 minutes using the TX-150 rotor, whereby acceptable results were achieved for the Urinalysis testing. And finally, a 2-minute centrifugation time at 4,500 rpm using the TX-150 rotor for blood samples were the centrifugation parameters suitable for stat clinical blood chemistry testing. Subsequently, centrifugation of peripheral blood was also successfully performed at 7,200 rpm for 1 minute using the HIGHConic rotor.

## Methods and materials

### A. Platelet-poor plasma (PPP)

Blood was obtained from 10 unknown donors by venipuncture collection at the same day of the testing. 2.7 mL of whole blood was collected into BD Vacutainer Plus plastic citrate tubes and after 30 minutes each sample was tested to get the initial platelet count before processing. The tubes were processed using a small benchtop centrifuge. The rotors and centrifugation conditions are summarized in Table 3. Platelet counts were performed on the whole blood and PPP samples using a hematology analyzer.

### B. Urinalysis

Urine specimens were pooled from 10 unknown patients and collected at the same day of the testing into 12 mL KOVA tubes. The tubes were processed using a small benchtop centrifuge until a moderately cohesive button was formed at the bottom of

the tubes (see Table 3). The supernatant was decanted and a volume of 1 mL was left inside the tube. After resuspending the sediment a drop was poured onto a glass slide, covered with a coverslip and examined under a microscope at 10x (low power field) and 40x magnifications (high power field).

### C. Blood chemistry analysis

Blood was collected into BD Vacutainer SST Plus tubes from 6 random donors (2 tubes per donor). After waiting 30 minutes for samples to clot, the tubes were centrifuged using a small benchtop centrifuge. The rotors and spin conditions are listed in Table 3. The serum was tested for potassium, aspartate aminotransferase (AST), phosphorus, and lactate dehydrogenase (LDH) using a chemistry analyzer.

## Data analysis

### A. Platelet-poor plasma (PPP)

The values were reported as the mean and the standard deviation (SD) of the Platelet concentrations from whole blood and PPP preparation.

### B. Urinalysis

The white blood cells (WBCs), red blood cells (RBCs), and epithelial cells counts were reported as number of cells/HPF (red blood cells per high power field); the counts of casts (cylindrical particles sometimes found in urine that are formed from coagulated protein secreted by kidney cells) were reported as number of casts/10 LPF (low power field). Bacteria, calcium oxalate crystals, uric acid crystals, mucous threads and yeast were all examined at LPF and estimated as “rare”, “few”, “moderate”, or “many”.

### C. Blood chemistry analysis

The mean and standard deviation were calculated.

**Table 2. Thermo Scientific Rotor Buckets & Adapter**

Purpose	Rotor	Buckets	Adapter
PPP; Blood chemistry analysis	TX-150 rotor (PN 75005701)	Round buckets (PN 75005702)	5/7 mL blood collection tube adapter (PN 75005739) Rubber pads (PN 76003266)
	HIGHConic rotor (PN 75005709)	–	5/7 mL adapter (PN 75005804) Rubber pads (PN 005700F)
Urinalysis	TX-150 rotor (PN 75005701)	Round buckets (PN 75005702)	14 mL urine tube adapter (PN 75005738)

**Table 3. Centrifugation conditions<sup>a</sup>**

Clinical testing	Rotor	RPM	RCF [x g]	Time <sup>b</sup> [min]
PPP	TX-150	4,500	3,260	10
	HIGHConic	8,700	10,155	1
Urinalysis	TX-150	1,700	465	4
Blood chemistry analysis	TX-150	4,500	3,260	2
	HIGHConic	7,200	6,967	1

a Room temperature

b Acceleration time/deceleration time: Standard

## Results

### A. Platelet-poor plasma (PPP)

See Table 4.

**Table 4. Platelet count in whole blood and PPP preparation from 10 unknown donors**

	WB platelet count <sup>a,b</sup> [K/ $\mu$ L] (Before spin)	Platelet count <sup>a,b</sup> [K/ $\mu$ L] (After spin)
TX-150	222.9 $\pm$ 94.6	2.5 $\pm$ 2.2
HIGHConic	178.7 $\pm$ 105.8	4.3 $\pm$ 1.3

a Centrifugation conditions shown in Table 3

b Mean  $\pm$  SD

### B. Urinalysis

The number of WBCs, RBCs, epithelial cells, casts, crystals and yeast are shown in Table 5 and 6.

### C. Blood chemistry analysis

The mean and SD values of potassium, AST, phosphorus, and LDH in serum are summarized in Table 7.

**Table 5. Urinalysis from 10 unknown patients**

Patient	WBC <sup>a,b</sup>	RBC <sup>a,b</sup>	Epithelial Cells <sup>a,b</sup>	Bacteria <sup>a,c</sup>	Calcium oxalate crystals <sup>a,c</sup>	Uric acid crystals <sup>a,c</sup>
1	6 - 10	> 50	1 - 5	rare	none	none
2	1 - 5	> 50	6 - 25	moderate	none	none
3	> 50	6 - 10	1 - 5	moderate	none	none
4	0 - 1	0 - 1	1 - 5	none	moderate	none
5	1 - 5	0 - 1	6 - 25	moderate	none	rare
6	1 - 5	0 - 1	1 - 5	none	none	–
7	10 - 25	6 - 10	6 - 25	many	none	–
8	10 - 25	1 - 5	1 - 5	rare	moderate	–
9	10 - 25	1 - 5	1 - 5	moderate	none	–
10	10 - 25	1 - 5	6 - 25	moderate	none	–

a Centrifugation conditions shown in Table 3

b hpf, high power field

c lpf, low power field

**Table 6. Urinalysis from 10 unknown patients**

Patient	Mucous threads <sup>a,b</sup>	Yeast <sup>a,b</sup>	Hyaline casts <sup>a,b</sup>	Coarse granular casts <sup>a,b</sup>	WBC casts <sup>a,b</sup>	Amorphous sediment <sup>a,b</sup>
1	none	few	none	–	–	rare
2	moderate	few	none	–	–	none
3	none	none	none	–	–	none
4	moderate	none	1 - 5	–	–	rare
5	few	few	0 - 1	–	–	rare
6	none	none	none	none	none	none
7	none	few	none	0 - 2	none	none
8	none	none	1 - 5	none	none	none
9	moderate	few	1 - 5	none	0 - 2	none
10	many	none	none	none	none	none

a Centrifugation conditions shown in Table 3

b lpf, low power field

**Table 7. Blood chemistry analysis from 6 unknown donors**

Rotor	K [mmol/L] <sup>a</sup>	AST [IU/L] <sup>a</sup>	PHOS [mg/dL] <sup>a</sup>	LDH [IU/L] <sup>a</sup>
<b>TX-150</b>	3.85 ± 0.19	19.83 ± 7.83	2.87 ± 0.27	172.17 ± 33.90
<b>HIGHConic</b>	4.35 ± 0.23	20.50 ± 4.32	3.90 ± 0.38	172.33 ± 27.98

<sup>a</sup> Mean ± SD

### Conclusion

With a spin of 10 minutes at 4,500 rpm using the TX-150 swinging bucket rotor, a compact centrifuge achieved a clinically acceptable PPP yield per laboratory standards. In addition, when this centrifuge was configured with its HIGHConic fixed angle rotor, acceptable PPP yield were achieved within 1 minute when spun at 8,700 rpm. With both rotor options the residual platelet counts in the final PPP sample were definitely < 10,000/μL commensurate of an acceptable platelet count the CLSI guidelines for PPP used for coagulation tests<sup>6</sup>. Therefore, this study has illustrated the effectiveness and efficiency by which the compact centrifuge generates not only a clinically acceptable PPP separation, but also reduces the time required for routine and prompt results.

Similarly, with urinalysis, the sample was also effectively and efficiently processed at 1,700 rpm in 4 minutes using the TX-150 swinging bucket rotor and at the end, the acceptable results were comparable with values typically obtained in the lab using other small clinical centrifuges. In addition, this same success was also seen with the results from the blood chemistry samples. With a 2 minute run at 4,500 rpm using the TX-150 swinging bucket rotor or a 1 minute run when the HIGHConic fixed angle rotor was spun at 7,200 rpm, the results were also commensurate with successful results traditionally seen by clinicians when using other clinical centrifuges.

The versatility of Thermo Scientific compact centrifuges enable them to be effective and efficient in performing the multiple centrifugation steps required in the various applications within a clinical lab.

### References:

1. Shahangian S. et al. (2005), Results of 2001 survey of hospital coagulation laboratories in the United States, Archives of pathology and laboratory medicine, 129, 47 - 60.
2. Fogazzi GB, Garigali G (2003), The clinical art and science of urine microscopy, Curr Opin Nephrol Hypertens, 12, 625 - 32.
3. Graham JC, Galloway A (2001), ACP best practice no. 167, the laboratory diagnosis of urinary tract infection. J Clin Pathol., 54, 911 - 9.
4. A. Van der Laarse (1979), Interference of the measurement of lactate dehydrogenase (LDH) activity in human serum and plasma by from blood cells, Clinica Chimica Acta, 99, 135 - 142.
5. Validation studies for Thermo Scientific small benchtop centrifuges were performed with the help of University Hospital, San Antonio, Rosemarie Martinez MT (ASCP) and Kathleen Born MT (ASCP).
6. Collection, transport and processing of blood specimens for testing plasma-based coagulation assays. Approved guidelines, 4th ed. Wayne, Pennsylvania, Clinical and Laboratory Standards Institute, 2003 (Document H21-A4).

 Learn more at [thermofisher.com/clinicalcentrifuges](https://thermofisher.com/clinicalcentrifuges)

**This is an IVD product used for clinical purposes. Availability and intended use may differ by location. It is the customer's responsibility to ensure that the performance of the product is suitable for customers' specific uses or applications.** © 2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. BD Vacutainer is a trademark of Becton Dickinson AG. KOVA is a registered trademark of Kova International Inc. **APN-12809400 1225**