

Standard Process for Whole Blood to Plasma or Serum

Parameters:

- **Storage Temperature:** -80°C
- **Final Product Volume:** 0.5mL
- **Final Product Vial:** 1.0 Matrix cryovial (Thermo Fisher catalog number 3740)

Procedure:

1. Spin whole blood tubes at 2000xg for 15 minutes at 4°C.
 - a. Set acceleration = 20% of centrifuge maximum
 - a. Deceleration = 20% of centrifuge maximum
2. Aseptically pipette 500 µL of plasma or serum into a labeled 1.0mL Matrix cryovial.
3. Store aliquots at -80°C until shipment.