



ISOFLUX™ System

SAMPLE SUBMISSION INSTRUCTIONS

Checklist for completion of sample submission:

Pre-collection	
<input type="checkbox"/>	Signed agreement in place to process samples (Purchase Order, Research Agreement, etc.)
<input type="checkbox"/>	As much advanced notice as possible is provided to Fluxion before collection (minimum of 1 day)
Sample Collection	
<input type="checkbox"/>	Samples are collected, packed, and sealed per guidelines
<input type="checkbox"/>	Shipment is sent with appropriate courier to exact address below (MORNING DELIVERY IS REQUIRED)
<input type="checkbox"/>	Sample Submission Form (paper copy) is included with shipment
<input type="checkbox"/>	Sample Submission Form (electronic copy) is emailed to support@fluxionbio.com

Fluxion Biosciences, Inc.
Attn: Sample Processing (Job #XXXX)
385 Oyster Point Blvd. #3
South San Francisco, CA 94080
USA

+1 (650) 241-4777
www.fluxionbio.com

IsoFlux Specimen Preparation Information for Fresh Whole Blood Samples

1. PREPARATION

- Prepare to ship samples the same day they are collected.
- Samples are accepted for delivery Monday – Friday.
- Please confirm with your shipper that the sample will arrive during these days.
- Materials required (these are suggested part numbers through VWR, you can use comparable alternatives, please comply with your institution's and federal shipping guidelines):
 - 10cc EDTA Blood Collection Tube: BD #366643 or similar (available from VWR as P/N BD-366643)
 - Styrofoam shipping container:
 - For 2 tubes (Foam Mailer for Two 5/7/10 mL Tubes, VWR #73320-066)
 - For 8 tubes (Foam Mailer for Eight 5/7/10 mL Tubes, VWR #73320-086)
 - Plastic biospecimen outer bags (For multiple blood tubes, VWR # 89400-274)
 - Outer shipping box (Assembled Foam Unit in Corrugated Carton, Interior Dimensions 9 1/2 x 7 1/2 x 5 5/8", VWR # 89413-472)

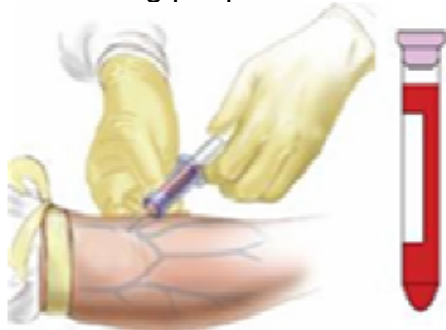
2. SAMPLE COLLECTION

a. Before collecting the sample:

- Label tubes with patient ID number(s).
- Complete the Sample Shipment Form.

b. Collecting the sample:

- Draw blood into 10cc EDTA Collection Tube using standard venipuncture procedures for collecting peripheral venous blood. Try to fill as much of the tube possible (>7mL).



- Invert the tube gently 8-10 times.

c. Packing the sample:

- The samples should be packed/shipped at Room Temperature. Do NOT refrigerate or freeze samples at any time.
- Place the collection tube(s) in the small Styrofoam shipping box.



- Place the small box into the plastic specimen bag and seal it using the seal-sealing strip



- Place the sealed bag into the insulated shipping box.



- Add a layer of bubble wrap on top to fill empty space.

- Enclose the Sample Shipment Form inside the cardboard box and seal up the box for shipment.

d. Precautions:

- **Do not expose to temperatures less than 4°C or greater than 37°C.**
- Do not freeze.
- Ship sample immediately (the day of collection).
- Handle sample with care – avoid shaking or hard inversion of the tubes.
- Hemolysis of the sample will affect cell isolation procedures. Overtly hemolyzed samples may be rejected at Fluxion's discretion.

3. TRANSFER TO FLUXION BIOSCIENCES

Prepare to ship the sample same day as collection for same day (local) or next day delivery. Shipping companies such as FEDEX, UPS, and local medical couriers all have same day or overnight delivery capabilities. Please contact Fluxion if you have any questions regarding shipments (650) 241-4777.

IMPORTANT: When shipping overnight, please select a morning delivery option (often called 'Priority Overnight' or equivalent). Samples that arrive late in the day may not be able to be processed.

Please use the exact shipping address as printed below

**Fluxion Biosciences
Attn: Sample Processing (Job # XXXXX)
385 Oyster Point Blvd.
Suite #3
South San Francisco, CA 94080
USA
+1 (650) 240-4777**

Please also include the following e-mail address in your shipping courier's electronic shipment notification update: support@fluxionbio.com

Adhere to D.O.T. and international shipping guidelines with respect to transport of biological materials. Please contact your courier or Fluxion for information if you are uncertain about the regulations for shipping various materials. Pack the samples in the appropriate materials as suggested by the guidelines included in this document.

IMPORTANT: Please inform your courier of the type of biological sample you are shipping and any associated details such as dry ice, cold packs, or potentially hazardous materials. Ensure any required forms or paperwork are filled out ahead of time to avoid delay.

IMPORTANT: Please ensure that your shipment will arrive to Fluxion's address on a weekday (Monday – Friday). Fluxion cannot be liable for shipments held up for delivery over the weekend.