

## Blood sample collection instructions

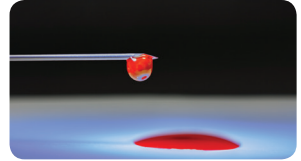
### Venipuncture



1. Disinfect patient's puncture site. Clean site with antiseptic.



2. Use standard EDTA venous collection procedures to prevent clotting.



3. Transfer one blood drop from syringe to each circle. Do not allow pipette to make contact with test strip. Be sure that each spot is completely saturated. Do not layer blood drops.

### Heel stick



1. Disinfect patient's heel. Clean heel with alcohol wipe and dry with sterile gauze.



2. Perform puncture using lancet device. Place heel upright and with lancet press down on either side of heel.



3. Wipe away first drop of blood. Let second, larger drop form. Gently touch the filter paper to the large drop and allow the blood to soak through and completely fill the circle. Apply blood to one side only. Do not allow heel to touch the sample application area, do not layer blood drops.

### Finger prick



1. Wash hands with soap and warm water. Dry with clean paper towel.

2. Disinfect patient's hand. Clean fingertip with alcohol wipe. Let hand hang down for 15-20 seconds to stimulate blood flow.



3. Perform puncture using lancet device. Place palm facing upward on hard surface and with lancet press down on finger.



4. Wipe away first drop of blood. Spot the blood drop on each circle, filling completely. Apply blood within circles, one side only. Do not allow finger to touch the sample application area, do not layer blood drops.

### Acceptable specimen



### Unacceptable specimens



No blood



Supersaturated blood

To ensure results are analyzed and returned expediently, follow these steps:

### Prepare paperwork and obtain sample

- Complete all fields on the test requisition and consent form using ball point pen.
- Complete patient name and date on the dried blood spot (DBS) card.
- Ensure the patient and/or parent(s) signature is captured on the consent form.
- Use preferred collection method to fill all spots on the DBS card (see reverse side).

### Drying collected blood spots

- Allow blood specimen to air dry horizontally at ambient temperature of 18° to 25°C (64° to 77°F), on non-absorbent open surface for approximately 3 hours.
  - Do not touch the blood spots while drying.
  - Keep protective flap open while drying.
  - Do not place test strips in front of air vents, use hair dryers, or other sources of moving air.
  - Test strips may be stored at room temperature for 7 days or at  $\geq -20^{\circ}\text{C}$  for up to 4 months.
  - Once dry, with the protective flap closed, samples can be stacked.

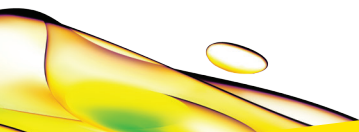
### Prepare for shipping

- Ship DBS card, test requisition form, and consent form as soon as test strip is fully dried.
- Ship by preferred method to: Revvity Omics, 250 Industry Drive, Pittsburgh, PA 15275 USA

### Frequently asked questions

- All results will be sent by portal, fax, or email to the recipient listed on the test requisition form.
- Notification of abnormal results will include a follow-up by a genetic counselor.
- For general questions on the collection and return of sample results, please contact Revvity at +1 (866) 354-2910 (Monday-Friday, 8:00AM - 8:00PM ET) or on our website at [www.revvity.com/category/clinical-services](http://www.revvity.com/category/clinical-services)

Information intended for use by healthcare professionals only.



revvity