

Finger stick sample requirements for DBS NAD assay

Recommendations to aid in collection

- 1) Run hands under warm water for a few minutes prior to collection.
- 2) Hydrate several hours before collection.

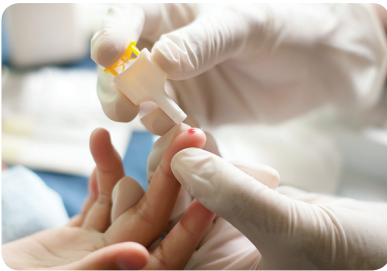


Disinfect and prick finger

1. Wash hands with soap and warm water. Dry with clean paper towel.
2. Select a finger and wipe end of finger

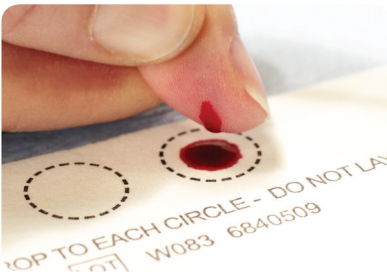
Puncture using lancet device

3. Prick finger using a sterile single-use lancet according to manufacturer protocol:
 - Twist cap in either direction (as indicated by arrows) to remove.
 - Press end into skin until lancet is released.
4. Dry wipe first spot of blood from finger with gauze before spotting.

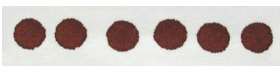


Wipe away first drop of blood. Let second drop fall onto circle. Do not fill circle.

5. Spot blood onto filter paper card in the center of the circle outlines.
 - Only ONE spot of blood per circle (blood volume should not exceed 125 μ L per spot according to manufacturer's instructions). DO NOT ATTEMPT TO FILL CIRCLE.
 - Do not spot on top of an existing blood spot.
 - Do not touch skin to the filter paper. If contact with skin occurs, wash hands immediately after collection.
 - Blood spot must soak through to back of card.
 - Back of filter paper should not be in contact with any surface while spotting or drying.
6. Allow blood spots to dry completely before storage (approximately 3 hours).
7. Place dry DBS card in biohazard bag with desiccant, seal, and store or ship.



Acceptable specimen



Unacceptable specimens



No blood

Insufficient blood

Supersaturated blood

Storage & shipping requirements

- DBS samples collected for NAD assay on Whatman® FTA mini filter paper cards may be stored at room temperature for up to 7 days or at $\geq 20^{\circ}\text{C}$ for up to 4 months.
- Samples should be shipped overnight on dry ice once dried or immediately after removal from storage.

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

Prepare paperwork and obtain sample

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

Drying collected blood spots

5. Allow blood specimen to air dry horizontally at ambient temperature of 18° to 25°C (64° to 77°F), on nonabsorbent 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.
 - Store in cool place until shipping (avoid heat and direct sunlight).
 - Do not stack.

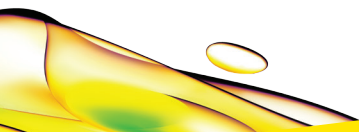
Prepare for shipping

6. Ship DBS card, test requisition form, and consent form as soon as test strip is fully dried.
7. 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 mail 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, 8:00 AM - 8:00 PM ET) or on our website at www.revvity.com/category/clinical-services

Disclaimer: This testing service has not been cleared or approved by the U.S. Food and Drug Administration. Testing services may not be licensed in accordance with the laws in all countries. The availability of specific test offerings is dependent upon laboratory location. The content of this pamphlet is provided for informational purposes only, not as medical advice. It is not intended to substitute the consultation, diagnosis, and/or treatment provided by a qualified licensed physician or other medical professional.



revvity