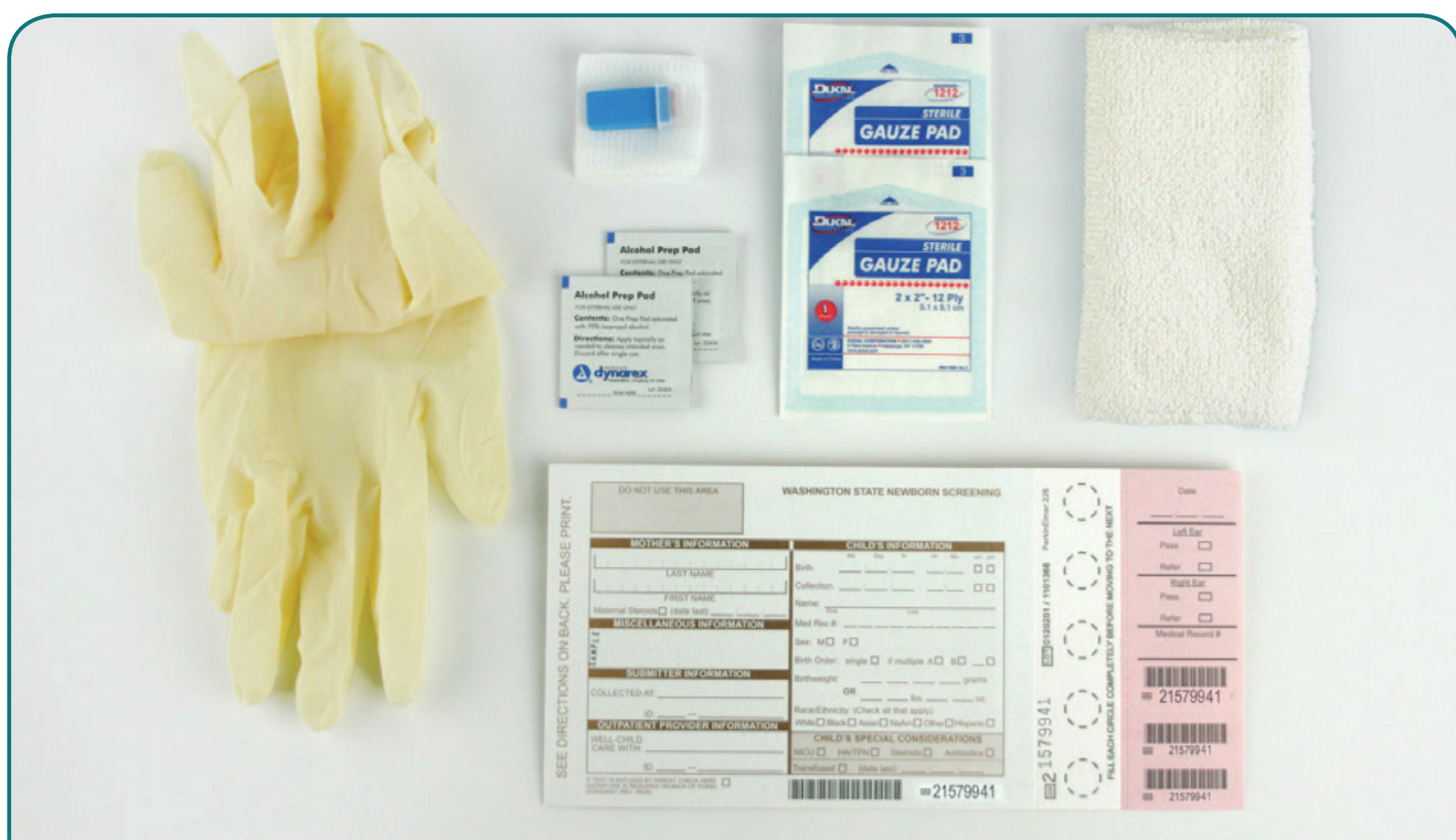
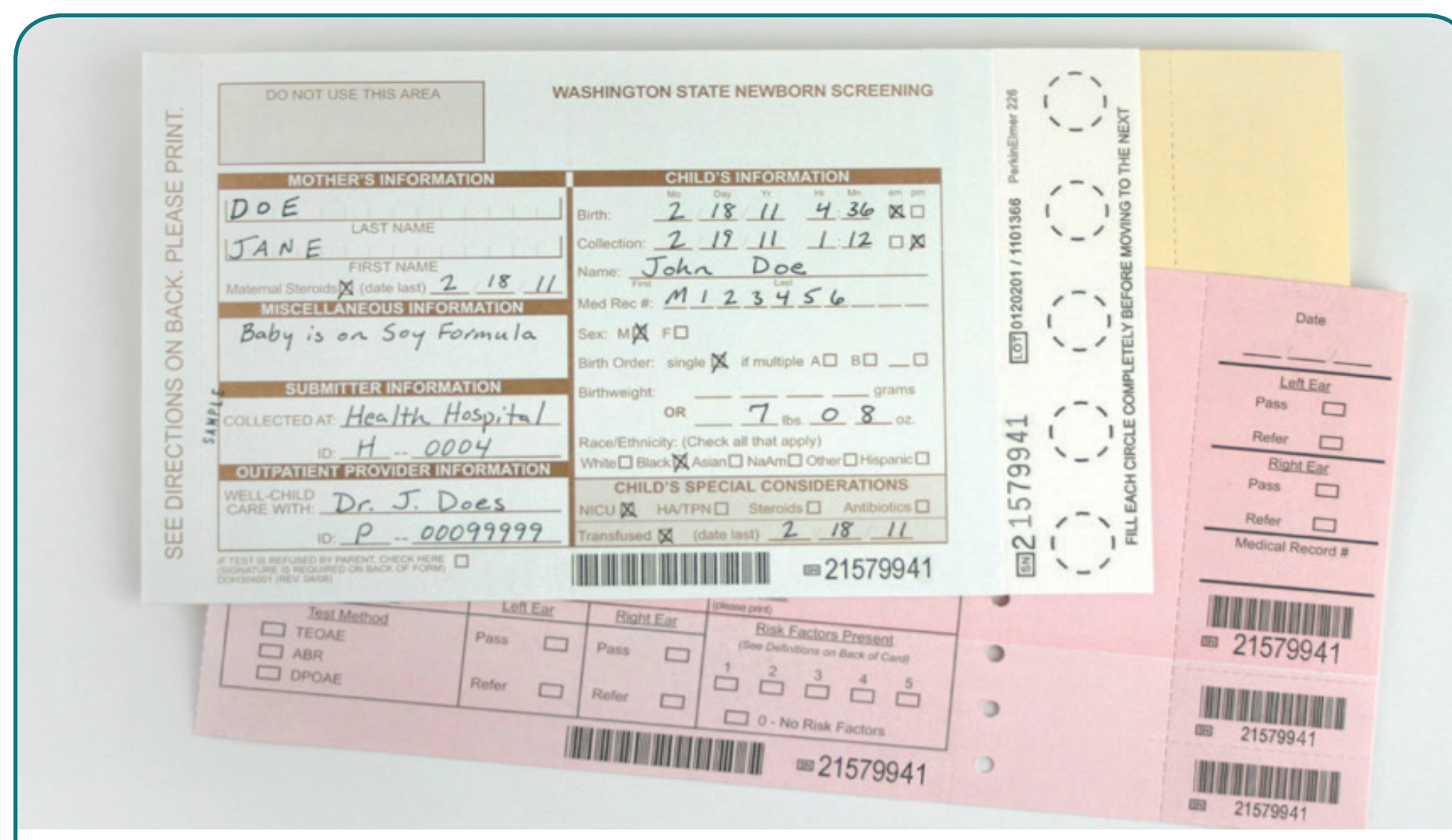


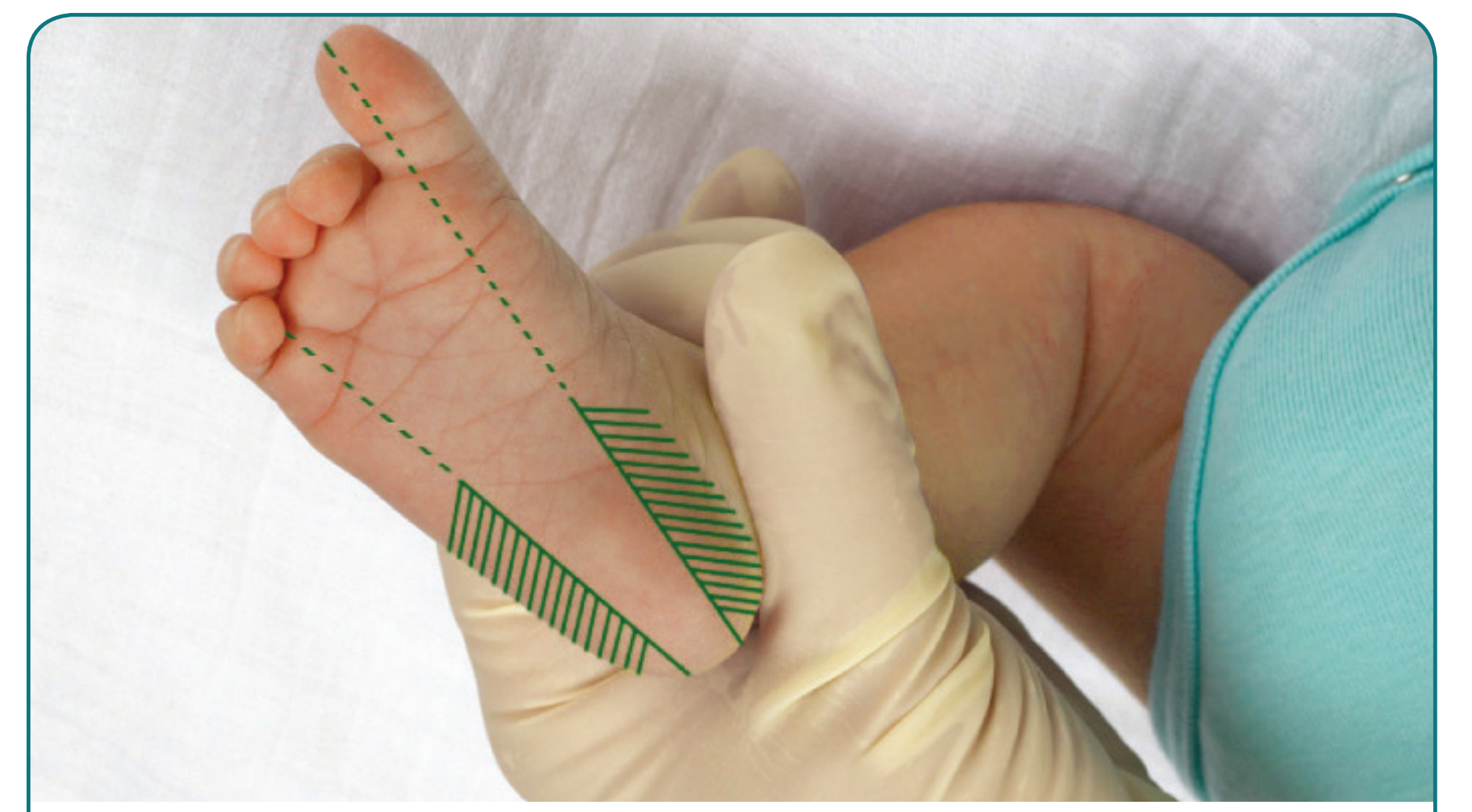
Newborn Screening: Sample Collection and Handling procedure



- Before collecting the sample, the following materials must be gathered:
- Sterile lancet with tip less than 2.4mm long
 - Sterile 70% alcohol pads
 - Sterile gauze pads
 - Warm, moist towel or compress
 - Fully completed, in-date filter paper blood collection form
 - Sterile gloves



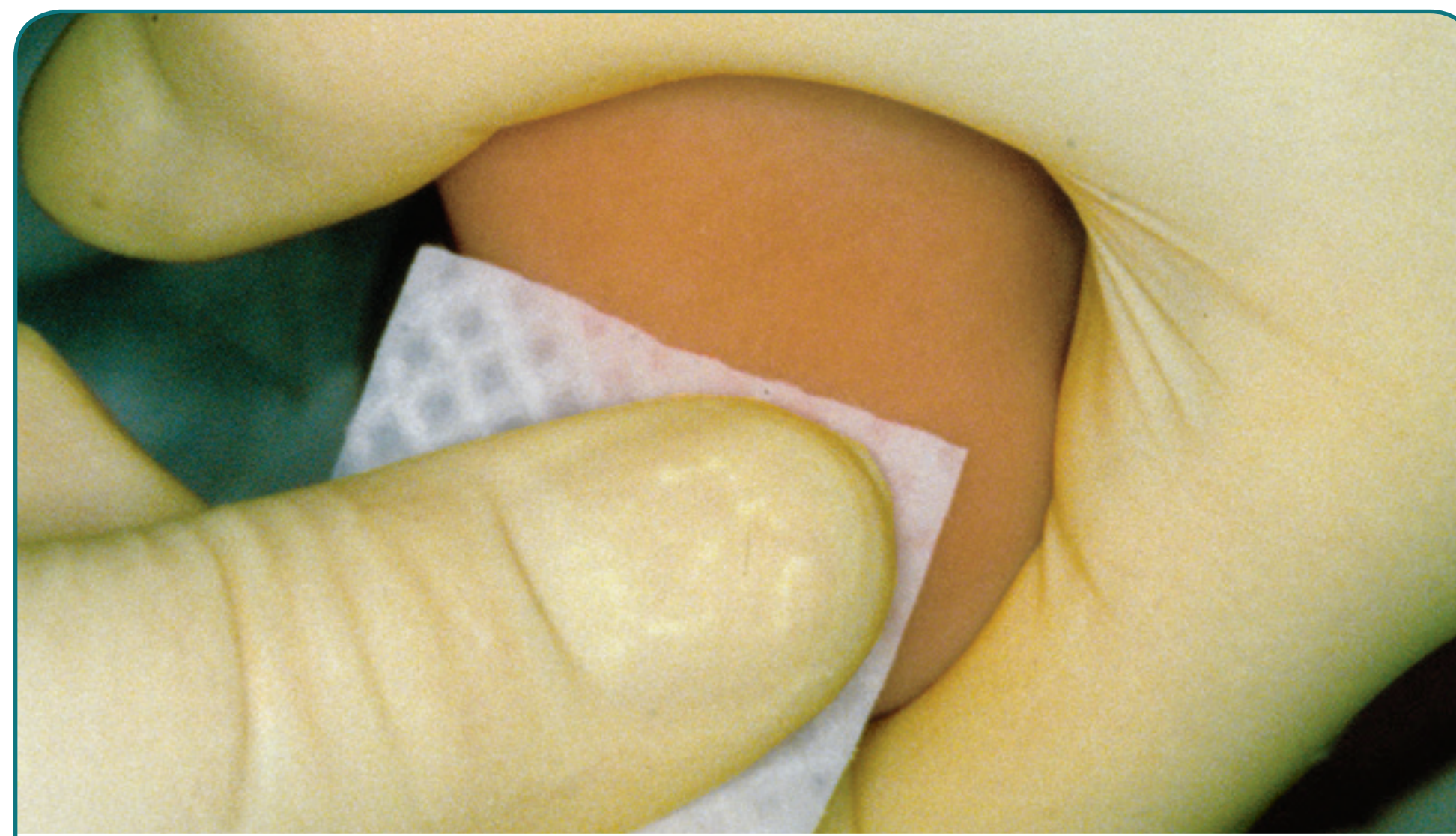
- 1** All fields on the form must be completed. Take steps to ensure that the filter paper is not touched or contaminated in any way during this process. If a copy of the form, such as a hearing copy, is required to be removed, do so at this time.



- 2** Areas on the side of the heel containing hatch marks are acceptable puncture sites.



- 3** Warm puncture site for 3-5 minutes with a warm, moist towel no hotter than 42°C.



- 4** Cleanse the puncture site with a sterile alcohol pad and allow to air dry.



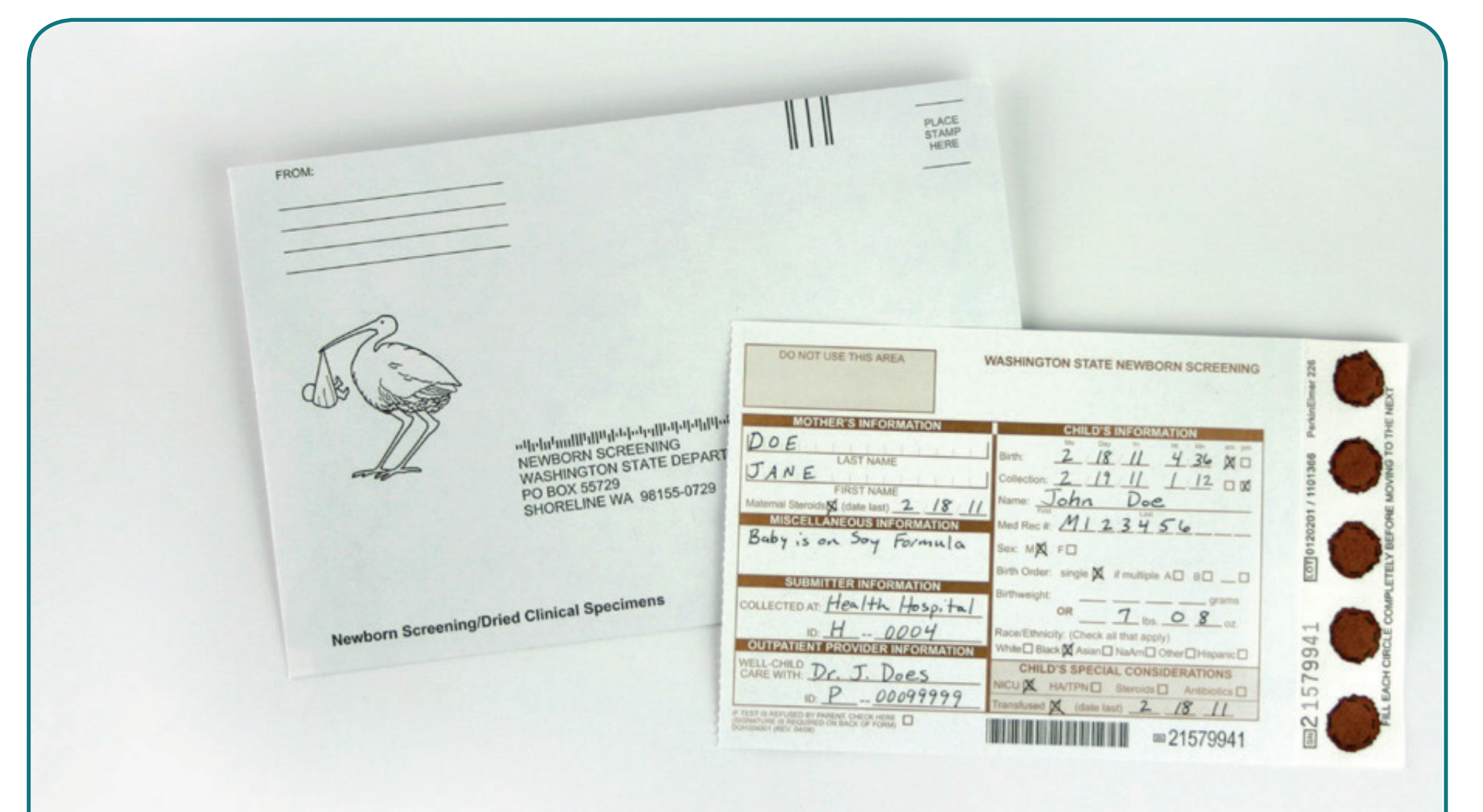
- 5** Puncture skin and wipe away the first drop of blood. Allow a second large drop to form.



- 6** Lightly touch the filter paper to the second drop of blood. Allow the blood to soak through and fill the preprinted circle. DO NOT APPLY MORE THAN ONE DROP TO A CIRCLE. Fill all remaining circles in this manner. If blood flow is diminished, repeat steps 3-6 with sterile equipment.



- 7** Allow the specimen to dry in a horizontal position for at least three hours, away from direct contact with surfaces, heat, and sunlight.



- 8** Send the specimen to the appropriate location no more than 24 hours after collection.

