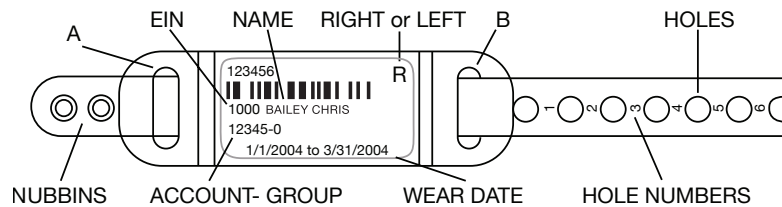


Dosimeter Finger Ring Badge Label information and wear instructions



Instructions for Fitting the Ring

1. Curl the ring into your approximate ring size and snap the ring together.
2. Try the ring on the finger you will be wearing it on.
3. Adjust the ring to the proper size, leaving a little room for expansion as you work.
4. Remember the number of the ring size (between the nubbins). This will help you quickly size the new ring you receive for the next wear period.
5. Remove the ring from your finger and tuck the strap end under the pouch through hole "A".
6. The ring is now ready to wear and can be easily adjusted during the month, if this is required. When the wear period is complete, simply return the whole ring (strap included) to RDC.
7. To quickly size the ring for the next wear period, simply adjust the strap to the correct ring size and tuck the strap under before putting it on.

An accredited dose monitoring program is usually required to demonstrate compliance with state and/or federal regulations. The thermoluminescent dosimeter (TLD) finger ring service is designed to provide valuable dose information for a facility's personnel monitoring program. To obtain accurate and meaningful data from our service, it is important to observe the following:

1. The ring should be worn at all times when there is a potential for occupational exposure to ionizing radiation.
2. Rings not in use should be stored in an area without radiation from your activities to prevent non-work related exposure.
3. Remember to return the control dosimeter with the rest of the dosimeters for the same wear period. The control is used to monitor background radiation and any dose received by the rings in transit. Any dose recorded is subtracted from the other rings, so it should not be used for any other purpose.
4. The TLD detector chip is located under the name label, so the ring should be worn with the label facing the source of radiation. If gloves are used, then the ring should be worn under the gloves. When the correct size adjustment is determined, the excess snap can be inserted under the TLD pouch.
5. The TLD pouch is heat sealed and water resistant. The plastic ring is able to be sterilized but is not resistant to all solvents.
6. Care should be taken to use only the ring marked with the proper individual's identification. Right and left hand rings are designated by "RU" or "LU" in the upper right corner of the label. If it is necessary to use an unassigned ring, or to reassign a ring, please provide the necessary information on a "Badge Reassignment Form" and enclose it when returning the rings. Do NOT cover or write on the pouch, as doing so may damage the chipstrate.
7. Replacement rings for the next monitoring period will be sent before the beginning of the next monitoring period. Begin wearing the rings on the date printed on the dosimeter. Return worn rings from the previous monitoring period, along with their control, in the self-addressed mailing envelope or box provided.
8. To avoid an unreturned dosimeter charge, return badges promptly after each wear period.

Find out more at thermofisher.com/dosimetrieservices