



The safest, easiest-to-use packaging for radiochemicals.

NENSure™ Packaging System

- Features a resealable screw cap for safe and convenient access and storage of product.
- Allows convenient access to the product by pipette or by syringe, without removing the cap.
- Has a square body to prevent rolling and an octagonally shaped cap that can be easily removed using the built-in wrench on top of the outer container.
- Thaws to room temperature in less than 20 minutes.
- Contains a rubber septum coated with Teflon® to ensure compatibility with biological systems - no toxic contaminants can leach through the rubber.

Light sensitive compounds are packaged in a red NENSure vial, which contains a UV inhibitor, instead of traditional foil-wrapped vials.



Figure 1: Detailed view of the NENSure packaging system

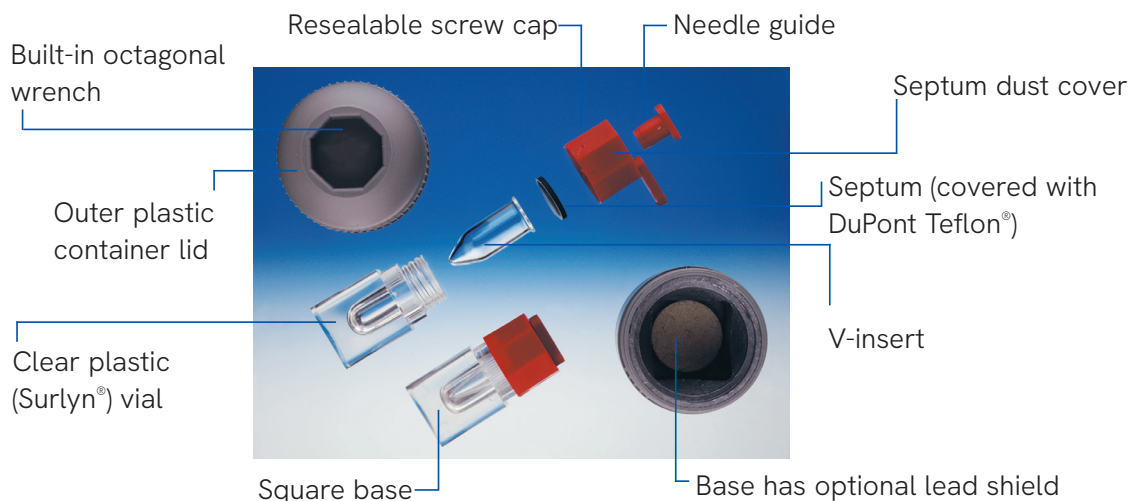


Figure 1: Detailed view of the NENSure packaging system.

First in safety and convenience NEN[®] radiochemicals from Revvity

Since 1985, the unique NENSure system has set the standard for radiochemical packaging. Based on proven performance over millions of shipments, the system provides:

- Ease of use
- Safe storage of unused product
- Reduced radiation exposure

Many NEN brand research products, including ³²P-, ³⁵S-, ³H-, ¹⁴C- and ¹²⁵I-labeled compounds, are packaged in the NENSure System, as are most radionuclides.

From radiochemicals to packaging, no one offers more to meet your research needs than Revvity.

Using a syringe to remove contents from NENSure vials

A 1.5 inch (38 mm) needle will reach the bottom of the V-insert when inserted directly through the septum of a 0.9 mL or 2.2 mL NENSure vial. A 2.25 inch (57 mm) needle is required for the 5 mL vial. The needle guide was designed to facilitate the use of microsyringes, such as the Hamilton syringe; however, it will accommodate needles of 22 gauge or finer. When using the needle guide, add 0.5 inches (12.7 mm) to the required needle length.

To expose the septum:

1. Unscrew the lid of the outer container.
2. Slide aside or remove the dust cover on the cap.
3. If required, install a charcoal trap directly through the septum.
4. Insert the syringe needle directly through the septum.

To use the needle guide:

1. Slide aside and remove the dust cover.
2. Insert and press the needle guide into the exposed opening.

3. Install a charcoal trap through the side opening of the guide until it just penetrates the septum.
4. Fit the syringe needle into the center hole of the guide and push firmly through the septum into the vial.

Using a pipette to remove contents from NENSure vials

To remove the cap:

1. Unscrew the lid of the outer container.
2. Invert the lid and position it so that the built-in octagonal wrench engages the cap.
3. Turn the wrench counterclockwise to remove the cap.

To replace the cap:

1. Position the cap on the vial.
2. Turn the wrench clockwise to screw the cap in place, ensuring a tight seal.

Installing a charcoal trap in NENSure vials

Prior to removing the vial contents or adding reagents, Revvity recommends installing a charcoal trap (NEX033T) to vent any volatiles that may have built up during shipment. This precautionary step is particularly important when opening ³⁵S-labeled compounds or ¹²⁵I nuclides.

The charcoal trap can be installed either directly through the septum or through the side opening of the needle guide (refer to figure and instructions).



| Figure 2: Charcoal trap.

revvity