

Low Volume Handling Suggestions for Iodine- 125 Labeled Compounds

Attention

Due to the high concentration of this compound, this vial may contain <10 μ L of volume.

When pipetting such small volumes we recommend the following steps to insure the correct amount of activity is attained:

For assays using the total product volume (one use of per vial)

- 1) Before the first use and the lid still secured, shake the vial 3 times in a downward motion (as if for a thermometer)
- 2) If possible, use the NENSure vial as your assay vessel
- 3) If your total assay volume is larger than the NENSure vial can hold, use your assay buffer to dilute the compound, then transfer contents to your assay vessel

For assays using only partial product volume (multiple uses from the same vial)

- 1) Before the first use and the lid still secured, shake the vial 3 times in a downward motion (as if for a thermometer)
- 2) Select the appropriate pipet, recommended for volumes <10 μ l
- 3) If possible, dilute the product in a buffer that will maintain the radiochemical purity of the compound. Dispense single use aliquots into suitable containers and store at an appropriate temperature. Refer to the product's technical data sheet available at perkinelmer.com/COA

For help with any aspect of your assay, contact PerkinElmer [Technical Support](#).