

### CERTIFICATE OF RADIOACTIVITY

#### Internal Standard Kit for Liquid Scintillation Counting

<sup>3</sup>H FOR AQUEOUS SOLVENTS (<sup>3</sup>H-W)

Product No. 1210-121A

This kit includes 40 internal standard capsules containing a tritium-labeled aqueous compound in solid form. The labeled compound, D-[5-<sup>3</sup>H(N)]-Glucose is produced by Revvity.

The absolute activity of the capsule is calibrated by comparison with the reference standards of tritiated toluene supplied by the National Institute of Standards and Technology (NIST), USA. The Standard Reference Material (SRM) No. 4947C is certified to have an estimated accuracy of ± 1.2%.

The mean value of the absolute activity having a ±0.5% error in relation to NIST SRM No. 4947C and the % C.V. of the distribution of activity among the capsules are based on results of a 5% random samples of the capsule batch.

<b>Activity</b>	199000	<b>DPM</b>	0.0896	<b>mCi</b>	3.3167	<b>kBq</b>
<b>% CV</b>	1.30%					
<b>Reference Date / Lot No.</b>	December 5, 2023					

**The total activity of the kit is:** 3.5856 mCi 132.6667 kBq

#### Precautions on Storage and Use

The chemical stability of the tracer compound and the press-through package provide stability better than 1% of the absolute activity for at least one year from the date of calibration. **Expiry Date:** December 5, 2025

The contents of these standards are exempt from NRC or Agreement State licensing requirements. "Radioactive Material - Not for Human Use - Introduction into Foods, Beverages, Cosmetics, Drugs or Medicinals or into Products Manufactured for Commercial Distribution is prohibited. Exempt Quantities Should Not Be Combined."

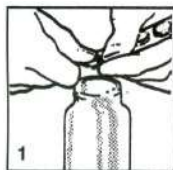
Approved by: Frank Passamonte  
Chemist

Date: 12/5/2023

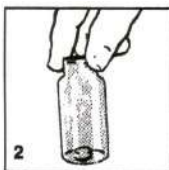
#### Instructions for use

This kit includes 40 Internal standard capsules with labeled solid state organic compound. The capsules are in a press-thru strip package and ready-for-use with organic scintillation systems.

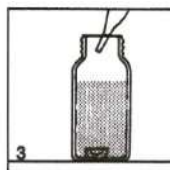
Ref. M, Reunanen and E. Soini: "New Method for obtaining the Quench Correction Curve in Liquid Scintillation Counting", Liquid Scintillation Counting, Vol. 3. ed. M. A. Crook, Heyden 1974.



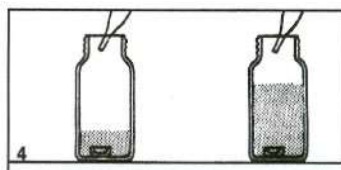
1 Press the desired, radioactive standard capsule into an empty sample vial. The capsules are delivered in press-thru strips. All standard 20ml LSC vials are acceptable.



2 Take care that the capsule is facing upwards in the vial as shown in fig. 2. The capsule can be turned by shaking the vial.



3 Samples soluble in organic solvents  
Pipette the whole scintillation cocktail into the sample vial. Shake the vial with a tight cap on. Let the radioactive standard dissolve ( few seconds ). NOTE Do not remove the glass capsule.



4 Samples soluble in water  
First pipette the total amount of water needed in the actual experiment into the vial. Let the radioactive standard dissolve for two minutes. Shake the vial with a tight cap on. Then add the scintillation liquid and shake again to mix the cocktail properly.

**CERTIFICAT DE RADIOACTIVITE**

Kit standard interne pour scintillation liquide de comptage

$^3\text{H}$  pour les solvants aqueux ( $^3\text{H-W}$ )

**Numéro de produit. 1210-121A**

Ce kit comprend 40 capsules internes standard contenant un composé marqué au tritium aqueux sous forme solide. Le composé marqué, D-[5- $^3\text{H}$  (N)]-glucose est produit par Revvity.

L'activité absolue de la capsule est étalonné par comparaison avec les normes de référence de toluène tritiée fournies par le National Institute of Standards and Technology (NIST), États-Unis. Le 4947C Matériaux de Référence n° (SRM) est certifié pour avoir une précision évaluée à  $\pm 1,2\%$ .

La valeur moyenne de l'activité absolue à une erreur de  $\pm 0,5\%$  par rapport à 4947C NIST n° SRM et le CV% de la répartition des activités entre les capsules sont basées sur les résultats d'un échantillon aléatoire de 5% du lot capsule.

<b>Activité</b>	199000	<b>DPM</b>	0.0896	<b>mCi</b>	3.3167	<b>kBq</b>
<b>% C.V</b>	1.30%					
<b>Date / de référence Numéro de lot</b>	5 décembre 2023					

<b>L'activité totale du kit est</b>	3.5856	<b>mCi</b>	132.6667	<b>kBq</b>
-------------------------------------	--------	------------	----------	------------

**Précautions sur le stockage et l'utilisation**

La stabilité chimique du composé marqueur et la presse-par paquet d'assurer la stabilité meilleure que 1% de l'activité absolue pour au moins un an à compter de la date de l'étalonnage. : **Date d'expiration: 5 décembre 2025**

Le contenu de ces normes sont exemptés du CNRC ou exigences de l'Accord de licence d'État. "Matières radioactives - Non à usage humain - Introduction dans les aliments, boissons, cosmétiques, médicaments ou Médicinal ou en produits manufacturés pour la distribution commerciale est interdite. Quantités exemptées devraient pas être combinés. "

**Approuvé par:** Frank Passamonte **Date** 5 décembre 2023  
**Chimiste**