

Warfarin, [<sup>14</sup>C]-

Product Number: NEC839

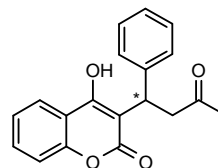
**LOT SPECIFIC INFORMATION**

Lot Number: 3112943

Specific Activity: 58.9 mCi/mmol

2179.3 MBq/mmol

Production Date: 21 Jun 2023



M.W. 308.33

C<sub>19</sub>H<sub>16</sub>O<sub>4</sub>

CAS 81-81-2

**PACKAGING:** Solid packaged under nitrogen in NENSURE with glass insert. Shipped in dry ice.**STABILITY AND STORAGE RECOMMENDATIONS:**

When warfarin, [<sup>14</sup>C]- is stored as a solid at -20°C, the rate of decomposition is initially less than 3% per year from the date of purification. The rate of decomposition can accelerate. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

**SPECIFIC ACTIVITY RANGE:** >40 mCi/mmol (>1480 MBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 96% when determined by the following method. It is advisable to check purity prior to use.

High performance liquid chromatography on a C18 column using the following mobile phase:

30 minute linear gradient from 60%A/40%B to 75%B

A) 10mM trifluoroacetic acid in water

B) acetonitrile

**QUALITY CONTROL:** The radiochemical purity of warfarin, [<sup>14</sup>C]- is checked at appropriate intervals using the above chromatography method.**HAZARD INFORMATION: WARNING:** This product contains a chemical known to the state of California to cause cancer.**Limited Warranty:** The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described

herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

NEC839-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.