

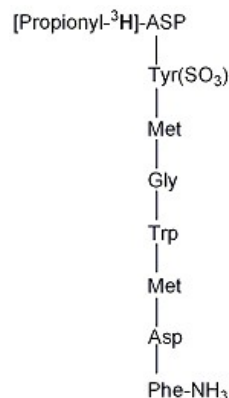
CCK-8 (sulphated) Propionylated, [Propionyl- $^3\text{H}(\text{N})$]-

Product Number: NET1162

LOT SPECIFIC INFORMATION

Lot Number:	3437894
Specific Activity:	110.2 Ci/mmol
	4077.4 GBq/mmol
Production Date:	2-Oct-25

M.W. 1199



C₅₂H₆₆N₁₀O₁₇S₃
CAS 177564-74-8

PACKAGING: 0.2 mCi/ml (7.4 MBq/ml) in triethylammonium phosphate (TEAP), pH 3.4 : acetonitrile (1:1), containing 0.2% 2-mercaptoethanol. Shipped in blue ice.

STABILITY AND STORAGE RECOMMENDATIONS: When CCK-8 (sulphated) propionylated, [propionyl- $^3\text{H}(\text{N})$]- is stored at 4°C in its original solvent and at its original concentration, the rate of decomposition is initially 4-8% per month 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: 60-145 Ci/mmol (2220-5365 GBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 90% when determined by the following method. It is advisable to check purity prior to use.

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

Isocratic 0.1% trifluoroacetic acid: acetonitrile (70:30)

QUALITY CONTROL: The radiochemical purity of CCK-8 (sulphated) propionylated, [propionyl- $^3\text{H}(\text{N})$]- 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.

NET1162-R Rev01

