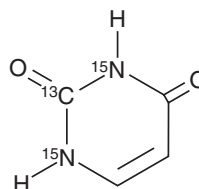


# PRODUCT INFORMATION



Uracil-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub>  
Item No. 33997

CAS Registry No.: 181517-11-3  
Formal Name: 2,4(1H,3H)-pyrimidinedione-2-<sup>13</sup>C-1,3-<sup>15</sup>N<sub>2</sub>  
MF: C<sub>3</sub>[<sup>13</sup>C]H<sub>4</sub>[<sup>15</sup>N]<sub>2</sub>O<sub>2</sub>  
FW: 115.1  
Purity: ≥95%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

Uracil-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> is supplied as a solid. A stock solution may be made by dissolving the uracil-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> in the solvent of choice, which should be purged with an inert gas. Uracil-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> is soluble in DMSO and a 1:1 solution of methanol:water. We do not recommend storing the aqueous solution for more than one day.

## Description

Uracil-<sup>13</sup>C,<sup>15</sup>N<sub>2</sub> is intended for use as an internal standard for the quantification of uracil (Item No. 26088). Uracil is a pyrimidine base and a fundamental component of RNA where it binds to adenine *via* hydrogen bonds.<sup>1</sup> It is converted into the nucleoside uridine through the addition of a ribose moiety, then to the nucleotide uridine monophosphate by the addition of a phosphate group.

## Reference

1. Voet, D. and Voet, J.G. *Biochemistry*. 3<sup>rd</sup> ed., John Wiley & Sons, Hoboken (2004).

### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/18/2024

## CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM