

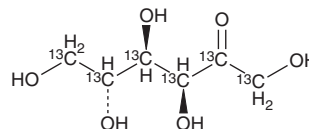
PRODUCT INFORMATION



D-Fructose-¹³C₆

Item No. 30071

CAS Registry No.: 201595-65-5
Formal Name: D-fructose-1,2,3,4,5,6-¹³C₆
Synonyms: D-(-)-Fructose-¹³C₆, D-(-)-Levulose-¹³C₆
MF: [¹³C]₆H₁₂O₆
FW: 186.1
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

D-Fructose-¹³C₆ is supplied as a crystalline solid. A stock solution may be made by dissolving the D-fructose-¹³C₆ in the solvent of choice, which should be purged with an inert gas. D-Fructose-¹³C₆ is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of D-fructose-¹³C₆ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of D-fructose-¹³C₆ in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

D-Fructose-¹³C₆ is intended for use as an internal standard for the quantification of D-fructose by GC- or LC-MS. D-Fructose is a ubiquitous monosaccharide and is derived, in addition to glucose, from the breakdown of sucrose by sucrase in the intestine.¹ It is a precursor in the biosynthesis of D-fructose-1,6-bisphosphate (Item No. 20516), which is an intermediate in the production of D-glucose *via* gluconeogenesis. Deficiencies in the enzymes that metabolize D-fructose are inborn errors of metabolism that range from benign, for fructokinase deficiency, to severe, for hereditary fructose intolerance, if D-fructose, sucrose, and sorbitol are not eliminated from the diet.² Increased consumption of D-fructose is associated with obesity, dyslipidemia, and impaired insulin sensitivity.³

References

1. Chen, M. and Whistler, R.L. Metabolism of D-fructose. *Adv. Carbohydr. Chem. Biochem.* **34**, 265-343 (1977).
2. Tran, C. Inborn errors of fructose metabolism. What can we learn from them? *Nutrients* **9**(4), E356 (2017).
3. Tappy, L. and Lê, K.-A. Metabolic effects of fructose and the worldwide increase in obesity. *Physiol. Rev.* **90**(1), 23-46 (2010).

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, 05/04/2020

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