

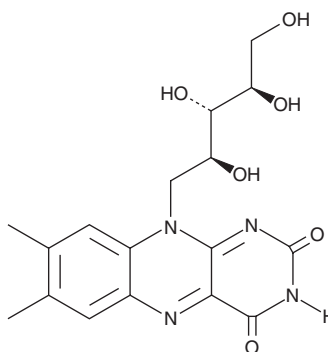
# PRODUCT INFORMATION



## Riboflavin

Item No. 30351

**CAS Registry No.:** 83-88-5  
**Formal Name:** 7,8-dimethyl-10-((2S,3S,4R)-2,3,4,5-tetrahydroxy-pentyl)benzo[g]pteridine-2,4(3H,10H)-dione  
**Synonyms:** Lactochrome, NSC 33298, Vitamin B<sub>2</sub>  
**MF:** C<sub>17</sub>H<sub>20</sub>N<sub>4</sub>O<sub>6</sub>  
**FW:** 376.4  
**Purity:** ≥98%  
**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

Riboflavin is supplied as a solid. A stock solution may be made by dissolving the riboflavin in the solvent of choice, which should be purged with an inert gas. Riboflavin is slightly soluble (0.1-1 mg/ml) in chloroform and methanol.

### Description

Riboflavin is an essential B vitamin.<sup>1</sup> It is a biosynthetic precursor to riboflavin 5'-monophosphate (Item No. 18167) and flavin adenine dinucleotide (FAD; Item No. 23386), which are necessary coenzymes in oxidation-reduction reactions involved in diverse biological processes, including electron transport and lipid and drug metabolism.<sup>1-3</sup> Formulations containing riboflavin have been used as dietary supplements.

### References

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. Section 25.5 NAD<sup>+</sup>, FAD, and coenzyme A are formed from ATP. *Biochemistry*. 5<sup>th</sup> edition, W.H. Freeman (2002).
2. Walsh, C.T. and Wencewicz, T.A. Flavoenzymes: Versatile catalysts in biosynthetic pathways. *Nat. Prod. Rep.* **30(1)**, 175-200 (2013).
3. Pinto, J.T. and Zemleni, J. Riboflavin. *Adv. Nutr.* **7(5)**, 973-975 (2016).

#### 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, 08/21/2025

#### 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