

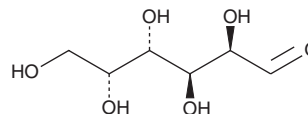
# PRODUCT INFORMATION



## D-Galactose

Item No. 20890

<b>CAS Registry No.:</b>	59-23-4
<b>Formal Name:</b>	(2R,3S,4S,5R)-2,3,4,5,6-pentahydroxyhexanal
<b>Synonym:</b>	D-(+)-Galactose
<b>MF:</b>	C <sub>6</sub> H <sub>12</sub> O <sub>6</sub>
<b>FW:</b>	180.2
<b>Purity:</b>	≥95%
<b>UV/Vis.:</b>	λ <sub>max</sub> : 260 nm
<b>Supplied as:</b>	A crystalline solid
<b>Storage:</b>	Room temperature
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

D-Galactose is supplied as a crystalline solid. A stock solution may be made by dissolving the D-galactose in the solvent of choice, which should be purged with an inert gas. D-Galactose is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of D-galactose in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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-galactose can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of D-galactose 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-Galactose is a natural aldohexose and C-4 epimer of glucose. D-galactose is converted enzymatically into D-glucose for metabolism or polysaccharides for storage. Chronic, systemic exposure to D-galactose accelerates senescence in invertebrates and mammals and has been used as a model for aging.<sup>1</sup> In bacteria, D-galactose is imported by a methyl-galactoside transport system to drive chemotaxis.<sup>2</sup>

### References

1. Cui, X., Zuo, P. Zhang, Q., *et al.* Chronic systemic D-galactose exposure induces memory loss, neurodegeneration, and oxidative damage in mice: Protective effects of R- $\alpha$ -lipoic acid. *J. Neurosci. Res.* **84**(3), 647-654 (2006).
2. Borrok, M.J., Kiessling, L.L., and Forest, K.T. Conformational changes of glucose/galactose-binding protein illuminated by open, unliganded, and ultra-high-resolution ligand-bound structures. *Protein Sci.* **16**(6), 1032-1041 (2007).

#### 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, 11/16/2022

#### 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