

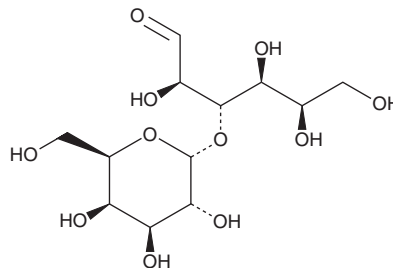
PRODUCT INFORMATION



Galactose- α -1,3-galactose

Item No. 45639

CAS Registry No.: 13168-24-6
Formal Name: 3-O- α -D-galactopyranosyl-D-galactose
Synonyms: α -Gal, 3 α -Galactobiose
MF: C₁₂H₂₂O₁₁
FW: 342.3
Purity: $\geq 95\%$
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

Galactose- α -1,3-galactose (α -Gal) is supplied as a solid. A stock solution may be made by dissolving the α -Gal in the solvent of choice, which should be purged with an inert gas. α -Gal is soluble (≥ 10 mg/ml) in DMSO.

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 α -Gal can be prepared by directly dissolving the solid in aqueous buffers. α -Gal is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

α -Gal is a carbohydrate allergen found in non-catarrhine mammals.^{1,2} It causes alpha-Gal syndrome (AGS), which is mediated by anti- α -Gal IgE antibodies primarily arising from tick bites, which results in red meat allergies. AGS can also originate from cetuximab exposure or xenotransplantation.

References

1. Wilson, J.M., Schuyler, A.J., Schroeder, N., *et al.* Galactose- α -1,3-galactose: Atypical food allergen or model IgE hypersensitivity? *Curr. Allergy Asthma Rep.* **17**(1), 8 (2017).
2. Cabezas-Cruz, A., Hodžić, A., Román-Carrasco, P., *et al.* Environmental and molecular drivers of the α -Gal syndrome. *Front. Immunol.* **10**, 1210 (2019).

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, 07/20/2026

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