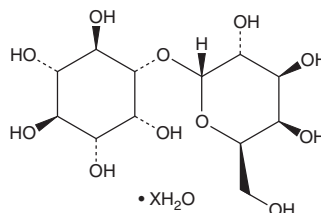


PRODUCT INFORMATION

Galactinol (hydrate)

Item No. 18426

Formal Name: 1-O- α -D-galactopyranosyl-myo-inositol, hydrate
MF: $C_{12}H_{22}O_{11} \cdot XH_2O$
FW: 342.3
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Galactinol (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the galactinol (hydrate) in the solvent of choice, which should be purged with an inert gas. Galactinol (hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 0.5 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 galactinol (hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of galactinol (hydrate) in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Galactinol, an α -galactoside of inositol that has been isolated from sugar beets, serves as a galactosyl donor in the transglycosidase-catalyzed biosynthesis of the raffinose family.¹ It is generated from myo-inositol and uridine diphosphate galactose in a galactosyltransferase-catalyzed reaction.¹

Reference

1. Mayer, T.G. and Schmidt, R.R. An efficient synthesis of galactinol and isogalactinol. *Liebigs Ann. Recl.* **859**, 859-863 (1997).

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, 12/02/2022

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