

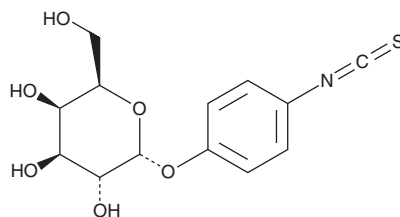
PRODUCT INFORMATION



α -D-Galactopyranosylphenyl isothiocyanate

Item No. 19544

CAS Registry No.: 120967-92-2
Formal Name: α -D-galactopyranoside, isothiocyanatophenyl
MF: $C_{13}H_{15}NO_6S$
FW: 313.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 228, 272, 283 nm
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

α -D-Galactopyranosylphenyl isothiocyanate is supplied as a crystalline solid. A stock solution may be made by dissolving the α -D-galactopyranosylphenyl isothiocyanate in the solvent of choice, which should be purged with an inert gas. α -D-Galactopyranosylphenyl isothiocyanate is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of α -D-galactopyranosylphenyl isothiocyanate in these solvents is approximately 20 and 30 mg/ml, respectively.

α -D-Galactopyranosylphenyl isothiocyanate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, α -D-galactopyranosylphenyl isothiocyanate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. α -D-Galactopyranosylphenyl isothiocyanate has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

α -D-Galactopyranosylphenyl isothiocyanate is a chemically activated form of galactose that has been used to prepare various neoglycoproteins, which consist of a glycosylated serum albumin substituted with either fluorescein or methotrexate.¹ Neoglycoproteins are used to identify the presence and sugar specificity of membrane lectins on the cell surface of mouse L1210 leukemia cells and to study membrane lectin-mediated endocytosis.¹

Reference

1. Monsigny, M., Roche, A. C., and Midoux, P. Uptake of neoglycoproteins via membrane lectin(s) of L1210 cells evidenced by quantitative flow cytometry and drug targeting. *Biol. Cell* **51**(2), 187-196 (1984).

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, 12/15/2022

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