

PRODUCT INFORMATION

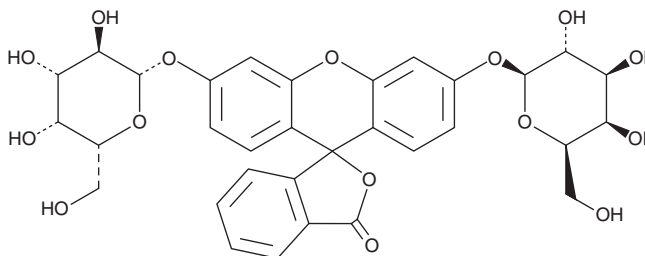


Fluorescein di-β-Galactopyranoside

Item No. 26860

CAS Registry No.: 17817-20-8
Formal Name: 3',6'-bis(β-D-galactopyranosyloxy)-
spiro[isobenzofuran-1(3H),9'-[9H]
xanthen]-3-one

Synonym: FDG
MF: C₃₂H₃₂O₁₅
FW: 656.6
Purity: ≥95%
Ex./Em. Max: 490/514 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Protect from light and moisture

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

Laboratory Procedures

Fluorescein di-β-galactopyranoside is supplied as a solid. A stock solution may be made by dissolving the fluorescein di-β-galactopyranoside in the solvent of choice, which should be purged with an inert gas. Fluorescein di-β-galactopyranoside is soluble in the organic solvent DMSO.

Description

Fluorescein di-β-galactopyranoside is a fluorogenic substrate for β-galactosidase.¹ Upon enzymatic cleavage by β-galactosidase, fluorescein is released and its fluorescence can be used to quantify β-galactosidase activity. Fluorescein displays excitation/emission maxima of 490/514 nm, respectively.^{1,2}

References

1. Plovins, A., Alvarez, A.M., Ibañez, M., *et al.* Use of fluorescein-di-β-D-galactopyranoside (FDG) and C₁₂-FDG as substrates for β-galactosidase detection by flow cytometry in animal, bacterial, and yeast cells. *Appl. Environ. Microbiol.* **60(12)**, 4638-4641 (1994).
2. Hermanson, G.T. Fluorescent probes. *Bioconjugate Techniques*. Audet, J., editor, 3rd edition, Elsevier (2013).

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, 11/03/2022

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