

# PRODUCT INFORMATION



**C<sub>12</sub>FDG**

Item No. 25583

**CAS Registry No.:** 138777-25-0  
**Formal Name:** N-[3',6'-bis(β-D-galactopyranosyloxy)-3-oxospiro [isobenzofuran-1(3H),9'-[9H] xanthen]-5-yl]-dodecanamide  
**Synonym:** 5-Dodecanoylaminofluorescein di-β-D-Galactopyranoside

**MF:** C<sub>44</sub>H<sub>55</sub>NO<sub>16</sub>

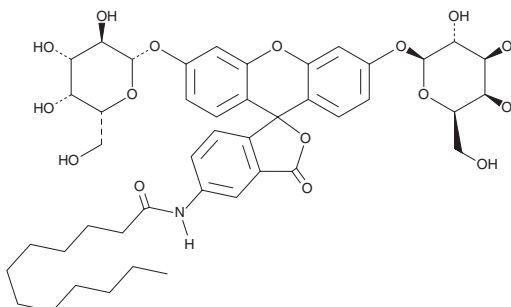
**FW:** 853.9

**Purity:** ≥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

C<sub>12</sub>FDG is supplied as a solid. A stock solution may be made by dissolving the C<sub>12</sub>FDG in the solvent of choice. C<sub>12</sub>FDG is soluble in the organic solvent DMSO, which should be purged with an inert gas.

## Description

C<sub>12</sub>FDG is a fluorogenic substrate for β-galactosidase.<sup>1</sup> 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 485/530 nm, respectively.

## Reference

1. Zhang, Y.-Z., Naleway, J.J., Larison, K.D., *et al.* Detecting lacZ gene expression in living cells with new lipophilic, fluorogenic β-galactosidase substrates. *FASEB J.* **5(15)**, 3108-3113 (1991).

### 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, 09/28/2022

## 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