

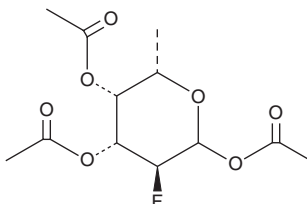
PRODUCT INFORMATION



2F-Peracetyl-Fucose

Item No. 30865

CAS Registry No.: 188783-78-0
Formal Name: 2,6-dideoxy-2-fluoro-L-galactopyranose 1,3,4-triacetate
Synonyms: 2F-PerAcFuc, 1,3,4-Tri-O-acetyl-2-deoxy-2-fluoro-L-fucose
MF: $C_{12}H_{17}FO_7$
FW: 292.3
Purity: $\geq 95\%$
Supplied as: A crystalline 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

2F-Peracetyl-fucose is supplied as an oil. A stock solution may be made by dissolving the 2F-peracetyl-fucose in the solvent of choice, which should be purged with an inert gas. 2F-Peracetyl-fucose is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml. 2F-Peracetyl-fucose is soluble in chloroform and methanol.

Description

2F-Peracetyl-fucose is an inhibitor of protein fucosylation.¹ It inhibits fucosylation of IgG1 monoclonal antibodies produced by CHO-DG44 cells when used at a concentration of 50 μM . It reduces intracellular GDP-fucose levels in CHO-DG44 cells in a concentration-dependent manner. 2F-Peracetyl-fucose (100 μM) also decreases intracellular fucose levels in *A. thaliana* seedlings.² 2F-Peracetyl-fucose decreases root length in *A. thaliana* seedlings in a concentration-dependent manner.

References

1. Okeley, N.M., Alley, S.C., Anderson, M.E., *et al.* Development of orally active inhibitors of protein and cellular fucosylation. *Proc. Natl. Acad. Sci. USA* **110**(14), 5404-5409 (2013).
2. Villalobos, J.A., Yi, B.R., and Wallace, I.S. 2-Fluoro-L-fucose is a metabolically incorporated inhibitor of plant cell wall polysaccharide fucosylation. *PLoS One* **10**(9), e0139091 (2015).

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, 03/21/2024

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