

PRODUCT INFORMATION



OVN73571

Item No. 36858

CAS Registry No.: 120173-57-1

Formal Name: N-[(9H-fluoren-9-ylmethoxy)carbonyl]-O-[3,4,6-tri-O-acetyl-2-(acetilamino)-2-deoxy- α -D-galactopyranosyl]-L-serine

Synonym: Fmoc-Ser(α -D-GalNAc(OAc)₃)-OH

MF: C₃₂H₃₆N₂O₁₃

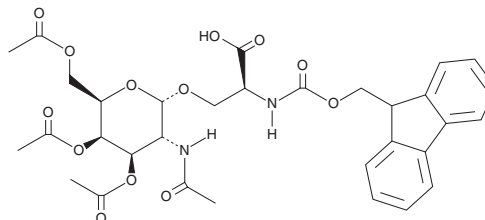
FW: 656.6

Purity: $\geq 98\%$

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

OVN73571 is supplied as a solid. A stock solution may be made by dissolving the OVN73571 in the solvent of choice, which should be purged with an inert gas. OVN73571 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of OVN73571 in DMF is approximately 16 mg/ml and approximately 14 mg/ml in ethanol and DMSO.

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 OVN73571 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of OVN73571 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

OVN73571 is a building block that has been used in the synthesis of the anticancer vaccine MAG-Tn3, a glycopeptide containing the tumor-associated Tn antigen.¹

Reference

1. Ganneau, C., Simenel, C., Emptas, E., *et al.* Large-scale synthesis and structural analysis of a synthetic glycopeptide dendrimer as an anti-cancer vaccine candidate. *Org. Biomol. Chem.* **15**(1), 114-123 (2017).

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, 12/21/2022

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