

PRODUCT INFORMATION



Semaglutide (FITC labeled)

Item No. 44292

Synonyms:	FITC-Semaglutide, FITC-SET
Purity:	≥95%
Ex./Em. Max.:	485/535 nm
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

Semaglutide (FITC labeled) is supplied as a solid. A stock solution may be made by dissolving the semaglutide (FITC labeled) in the solvent of choice, which should be purged with an inert gas. Semaglutide (FITC labeled) is soluble (≥10 mg/ml) in 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 semaglutide (FITC labeled) can be prepared by directly dissolving the solid in aqueous buffers. Semaglutide (FITC labeled) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Semaglutide (FITC labeled) is a FITC-conjugated form of the glucagon-like peptide 1 (GLP-1) agonist semaglutide (Item Nos. 40231 | 29969 | 40170). Semaglutide (FITC labeled) has been used in the study of milk exosome-based hybrid vesicle movement and penetration capacity in porcine intestinal mucus *ex vivo*.¹ It has also been used in the study of hybrid vesicle permeability in Caco-2/E12 co-cultured cells as a model of mucosal tissue and in the study of intestinal epithelial uptake in rats. FITC displays excitation/emission maxima of 485/535 nm, respectively.

Reference

1. Xiao, P., Yuan, H., Liu, H., *et al.* Modulating the elasticity of milk exosome-based hybrid vesicles to optimize transepithelial transport and enhance oral peptide delivery. *J. Control Release* **380**, 36-51 (2025).

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, 12/12/2025

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