

PRODUCT INFORMATION



Sulfo-SMCC (sodium salt)

Item No. 35692

CAS Registry No.: 92921-24-9
Formal Name: 4-[(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)methyl]-cyclohexanecarboxylic acid, 2,5-dioxo-3-sulfo-1-pyrrolidinyl ester, monosodium salt

Synonym: Sulfosuccinimidyl 4-(N-maleimidomethyl) cyclohexane-1-carboxylate

MF: C₁₆H₁₇N₂O₉S • Na

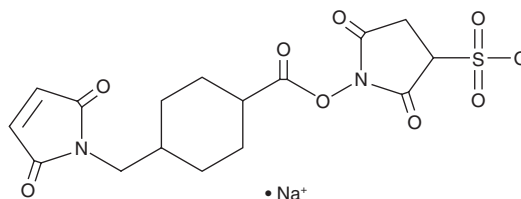
FW: 436.4

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

Sulfo-SMCC (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the sulfo-SMCC (sodium salt) in the solvent of choice, which should be purged with an inert gas. Sulfo-SMCC (sodium salt) is soluble in acetonitrile. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfo-SMCC is a water-soluble protein cross-linking reagent that contains an anime-reactive sulfo-N-hydroxysuccinimide (sulfo-NHS) ester and a sulfhydryl-reactive maleimide group that can be used in the conjugation of proteins.^{1,2} It has been used to conjugate anti-programmed cell death ligand 1 (PD-L1) antibodies to the surface of platelets to induce antitumor immunity in xenograft models, as well as to conjugate Golgi-targeting peptides to platelet microparticle-mimetic nanoplatfoms for the delivery of all-trans retinoic acid (Item No. 11017) to isolated human synovial fibroblasts from patients with rheumatoid arthritis.

References

1. Gao, Y., Chen, X., Wang, B., *et al.* Engineering platelets with PDL1 antibodies and iron oxide nanoparticles for postsurgical cancer immunotherapy. *ACS Appl. Bio Mater.* **6**(1), 257-266 (2023).
2. Deng, C., Zhao, X., Chen, Y., *et al.* Engineered platelet microparticle-membrane camouflaged nanoparticles for targeting the golgi apparatus of synovial fibroblasts to attenuate rheumatoid arthritis. *ACS Nano* **16**(11), 18430-18447 (2022).

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, 03/27/2024

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