

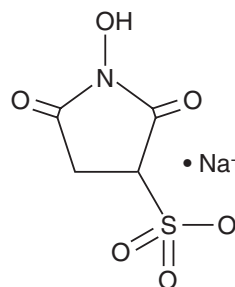
PRODUCT INFORMATION



N-Hydroxysulfosuccinimide (sodium salt)

Item No. 20680

CAS Registry No.: 106627-54-7
Formal Name: 1-hydroxy-2,5-dioxo-3-pyrrolidinesulfonic acid, monosodium salt
Synonym: Sulfo-NHS
MF: $C_4H_4NO_6S \cdot Na$
FW: 217.1
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 210 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

N-Hydroxysulfosuccinimide (sodium salt) (sulfo-NHS) is supplied as a crystalline solid. Aqueous solutions of sulfo-NHS can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of sulfo-NHS in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfo-NHS is a modifying reagent that, in the presence of a cross-linking reagent like carbodiimide or EDAC, converts carboxyl groups to amine-reactive sulfo-NHS esters.¹⁻³ The cross-linker converts the carboxylic acid to an unstable *o*-acylisourea intermediate, which then reacts with the primary amine of sulfo-NHS to produce the stable product.

References

1. Bogdanov, Jr., A.A., Klibanov, A.L and Torchilin, V.P. Protein immobilization on the surface of liposomes via carbodiimide activation in the presence of N-hydroxysulfosuccinimide. *FEBS Lett.* **231(2)**, 381-384 (1988).
2. Lelong, C., Sétif, P., Lagoutte, B., *et al.* Identification of the amino acids involved in the functional interaction between photosystem I and ferredoxin from *Synechocystis* sp. PCC 6803 by chemical cross-linking. *J. Biol. Chem.* **269(13)**, 10034-10039 (1994).
3. Klyachko, N.L., Manickam, D.S., Brynskikh, A.M., *et al.* Cross-linked antioxidant nanozymes for improved delivery to CNS. *Nanomedicine* **8(1)**, 119/129 (2012).

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, 11/29/2022

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