

# PRODUCT INFORMATION



## Sulfosuccinimidyl 3-[[2-(Biotinamido)ethyl]dithio]propionate

Item No. 39537

CAS Registry No.: 325143-98-4

Formal Name: 3-[[2-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]ethyl]dithio]-propanoic acid, monosodium salt, 2,5-dioxo-3-sulfo-1-pyrrolidinyl ester

Synonym: Sulfo-NHS-SS-biotin

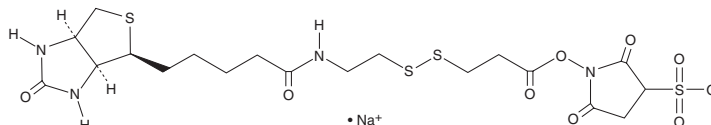
MF:  $C_{19}H_{27}N_4O_9S_4 \cdot Na$

FW: 606.7

Supplied as: A solid

Storage: -20°C

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Sulfosuccinimidyl 3-[[2-(biotinamido)ethyl]dithio]propionate (sulfo-NHS-SS-biotin) is supplied as a solid. A stock solution may be made by dissolving the sulfo-NHS-SS-biotin in the solvent of choice, which should be purged with an inert gas. Sulfo-NHS-SS-biotin is slightly soluble in DMSO and methanol.

### Description

Sulfo-NHS-SS-biotin is a membrane-impermeable amine-reactive reagent that contains biotin (Item No. 22582) conjugated to N-hydroxysulfosuccinimide (sulfo-NHS; Item No. 20680) via a disulfide bond.<sup>1,2</sup> It has been used as a derivatization reagent to identify post-translational modifications (PTMs) in lysine-rich proteins and novel cell surface proteins of *R. rickettsii*.

### References

1. Markoutsas, S., Bahr, U., Papasotiriou, D.G., *et al.* Sulfo-NHS-SS-biotin derivatization: A versatile tool for MALDI mass analysis of PTMs in lysine-rich proteins. *Proteomics* **14**(6), 659-667 (2014).
2. Gong, W., Xiong, X., Qi, Y., *et al.* Identification of novel surface-exposed proteins of *Rickettsia rickettsii* by affinity purification and proteomics. *PLoS One* **9**(6), e100253 (2014).

#### 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, 09/22/2023

#### 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