

PRODUCT INFORMATION



Polysarcosine150

Item No. 39051

Formal Name: N-butyl-2-(methyl((methylamino)methyl)amino)acetamide

Synonym: pSar₁₅₀

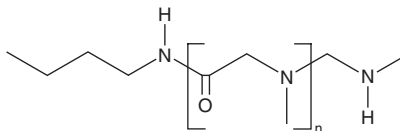
MF: (C₃H₅NO)_nC₆H₁₆N₂

Purity: ≥90%

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

Polysarcosine150 is supplied as a solid. A stock solution may be made by dissolving the polysarcosine150 in the solvent of choice, which should be purged with an inert gas. Polysarcosine150 is soluble in chloroform.

Description

Polysarcosine150 is an amine-functionalized hydrophilic polymer.¹ Surface modification of liposomes with polysarcosine polymers between 10 and 400 units long have been used to increase circulation half-life.

Reference

1. Weber, B., Birke, A., Fischer, K., *et al.* Solution properties of polysarcosine: From absolute and relative molar mass determinations to complement activation. *Macromolecules* **51**(7), 2653–2661 (2018).

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, 02/08/2024

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