

# PRODUCT INFORMATION



## Polysarcosine100

Item No. 39050

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

**Synonym:** pSar<sub>100</sub>

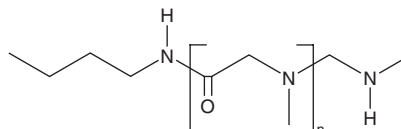
**MF:** (C<sub>3</sub>H<sub>5</sub>NO)<sub>n</sub>C<sub>6</sub>H<sub>16</sub>N<sub>2</sub>

**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

Polysarcosine100 is supplied as a solid. A stock solution may be made by dissolving the polysarcosine100 in the solvent of choice, which should be purged with an inert gas. Polysarcosine100 is soluble in chloroform and methanol.

### Description

Polysarcosine100 is an amine-functionalized hydrophilic polymer.<sup>1-3</sup> Surface modification of liposomes with polysarcosine polymers between 10 and 100 units long have been used to increase circulation half-life.<sup>2,3</sup>

### References

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).
2. Weber, B., Seidl, C., Schwiertz, D., *et al.* Polysarcosine-based lipids: From lipopolypeptoid micelles to stealth-like lipids in Langmuir Blodgett monolayers. *Polymers (Basel)* **8(12)**, 427 (2016).
3. Bleher, S., Buck, J., Muhl, C., *et al.* Poly(sarcosine) surface modification imparts stealth-like properties to liposomes. *Small* **15(50)**, e1904716 (2019).

#### 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