

PRODUCT INFORMATION



Polysarcosine20

Item No. 39048

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

Synonym: pSar₂₀

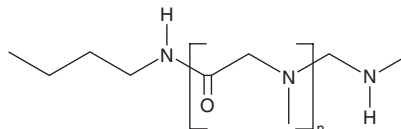
MF: (C₉H₁₆NO)_nC₆H₁₆N₂

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

Polysarcosine20 is supplied as a solid. A stock solution may be made by dissolving the polysarcosine20 in the solvent of choice, which should be purged with an inert gas. Polysarcosine20 is soluble in chloroform and methanol.

Description

Polysarcosine20 is an amine-functionalized hydrophilic polymer.¹ Addition of polysarcosine20 to surface-grafted polymer brushes reduces protein absorption and mammalian and bacterial cell attachment, indicating potential use as an antifouling agent. Surface modification of liposomes with polysarcosine polymers between 10 and 100 units long have been used to increase circulation half-life.^{2,3}

References

1. Lau, K.H.A., Ren, C., Sileika, T.S., *et al.* Surface-grafted polysarcosine as a peptoid antifouling polymer brush. *Langmuir* **28**(46), 16099-16107 (2012).
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, 02/20/2024

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