

PRODUCT INFORMATION

Azide-Polysarcosine150

Item No. 39041

Formal Name: N-(3-azidopropyl)-2-(methyl((methylamino)methyl)amino)acetamide

Synonyms: Azide-pSar₁₅₀, N₃-pSar₁₅₀, N₃-Polysarcosine150

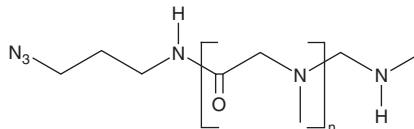
MF: (C₃H₅NO)_nC₅H₁₃N₅

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

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

Description

Azide-polysarcosine150 is an azide-containing form of the amine-functionalized hydrophilic polymer polysarcosine150 (Item No. 39051).¹ Azide-polysarcosines have been used in the formation of hydrophobicity-masking antibody-drug conjugates for the selective release of monomethyl auristatin E (Item No. 16267) to tumors in rats.

Reference

1. Viricel, W., Fournet, G., Beaumel, S., *et al.* Monodisperse polysarcosine-based highly-loaded antibody-drug conjugates. *Chem. Sci.* **10**(14), 4048-4053 (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, 02/26/2024

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