

PRODUCT INFORMATION



PAMAM Dendrimer G2.0 Succinamic Acid (water solution)

Item No. 39119

Synonyms: PAMAM G2.0 Succinamic Acid, Polyamidoamine Dendrimer G2.0 Succinamic Acid
FW: 4,856.0
Supplied as: A solution in water
Storage: -20°C
Stability: ≥2 years

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

Description

PAMAM dendrimer G2.0 succinamic acid (PAMAM G2.0 succinamic acid) is a polyamidoamine (PAMAM) dendrimer with succinamic acid termini.¹ It has 16 surface groups. PAMAM G2.0 succinamic acid is an antagonist for the pore-forming channels anthrax toxin protective antigen 63 (PA63; IC₅₀ = 879 nM) in lipid bilayers.² It is active against *S. aureus*.³ PAMAM G2.0 succinamic acid inhibits plaque formation induced by Middle East respiratory syndrome coronavirus (MERS-CoV) in Vero cells when used at a concentration of 10 μM.¹

References

1. Kandeel, M., Al-Taher, A., Park, B.K., et al. A pilot study of the antiviral activity of anionic and cationic polyamidoamine dendrimers against the Middle East respiratory syndrome coronavirus. *J. Med. Virol.* **92(9)**, 1665-1670 (2020).
2. Yamini, G., Kalu, N., and Nestorovich, E.M. Impact of dendrimer terminal group chemistry on blockage of the anthrax toxin channel: A single molecule study. *Toxins (Basel)* **8(11)**, 337 (2016).
3. Altaher, Y. and Kandeel, M. Structure-activity relationship of anionic and cationic polyamidoamine (PAMAM) dendrimers against *Staphylococcus aureus*. *J. Nanomater.* 4013016 (2022).

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, 12/08/2023

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