

PRODUCT INFORMATION

PAMAM Dendrimer G6.0 Succinamic Acid (water solution)

Item No. 39123

Synonyms: PAMAM G6.0 Succinamic Acid, Polyamidoamine Dendrimer G6.0 Succinamic Acid
FW: 83,648.0
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 G6.0 succinamic acid (PAMAM G6.0 succinamic acid) is a succinamic acid-terminated PAMAM dendrimer. It is approximately 76 Å in diameter in PBS and has 256 surface groups.^{1,2} PAMAM G6.0 succinamic acid does not induce mortality in zebrafish embryos when used at a concentration of 100 ppm in the tank water.²

References

1. Dobrovolskaia, M.A., Patri, A.K., Simak, J., *et al.* Nanoparticle size and surface charge determine effects of PAMAM dendrimers on human platelets *in vitro*. *Mol. Pharm.* **9(3)**, 382-393 (2012).
2. Pryor, J.B., Harper, B.J., and Harper, S.L. Comparative toxicological assessment of PAMAM and thiophosphoryl dendrimers using embryonic zebrafish. *Int. J. Nanomedicine* **9**, 1947-1956 (2014).

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