

PRODUCT INFORMATION



PAMAM Dendrimer G6.0 Amidoethanol (water solution)

Item No. 39115

Synonyms:	PAMAM 6.0 Amidoethanol, Polyamidoamine Dendrimer G6.0 Amidoethanol
MF:	$[\text{NH}_2(\text{CH}_2)_2\text{NH}_2]:(\text{G}=6); \text{dendri}$ $[\text{NH}_2(\text{CH}_2)_2\text{NH}_2]:(\text{G}=6); \text{dendri PAMAM}(\text{NHCH}_2\text{CH}_2\text{OH})_{256}$ PAMAM(NHCH ₂ CH ₂ OH) ₂₅₆
FW:	58,298.2
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 amidoethanol is a polyamidoamine (PAMAM G6.0-OH) dendrimer with hydroxyl termini. It is approximately 65 Å in diameter in PBS.¹ PAMAM G6.0-OH induces mortality in zebrafish embryos to a lesser extent than PAMAM dendrimer G6.0 amine (Item No. 39076) but to a greater extent than PAMAM dendrimer G6.0 succinamic acid (Item No. 39123) when used at a concentration of 250 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, 12/12/2023

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