

PRODUCT INFORMATION



GPMA

Item No. 9004846

CAS Registry No.: 231934-29-5

Formal Name: N-[3-[(aminoiminomethyl)amino]propyl]-2-methyl-2-propenamide

Synonyms: GuaMA, 3-Guanidinopropyl Methacrylamide, N-(3-Methacrylamidopropyl)-guanidinium, 2-Methacrylic Acid 3-guanidino-propyl ester, MGUA

MF: $C_8H_{16}N_4O$

FW: 184.2

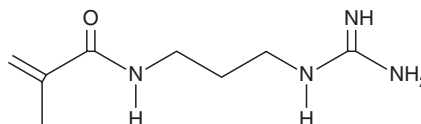
Purity: $\geq 90\%$

Supplied as: A 10 mg/ml solution in methanol

Storage: -20°C

Stability: ≥ 2 years

Special Conditions: Store under nitrogen



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

Description

GPMA is a cationic copolymerization agent.^{1,2} It has been used in the generation of polymer complexes for the delivery of plasmid DNA *in vitro*.

References

1. Funhoff, A.M., van Nostrum, C.F., Lok, M.C., *et al.* Poly(3-guanidinopropyl methacrylate): A novel cationic polymer for gene delivery. *Bioconjug. Chem.* **15**(6), 1212-1220 (2004).
2. Li, X., Qin, Z., Wu, Y., *et al.* Improvement of transfection efficiency by galactosylated N-3-guanidinopropyl methacrylamide-co-poly (ethylene glycol) methacrylate copolymers. *Eur. Polym. J.* **49**(10), 2884-2894 (2013).

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, 09/19/2025

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