

Technical Bulletin

Polyester-8-hydroxyl-1-carboxyl bis-MPA dendron, generation 3

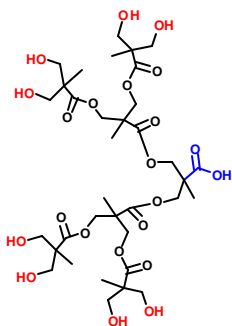
Catalog Number **686670**

Storage Temperature 2–8 °C

Technical Bulletin AL-240

Synonym: Dendron-G3-Carboxyl-OH

Product Description



Theoretical MW: 830.82 g/mol

Average Molecular Formula: $C_{35}H_{58}O_{22}$

Number of hydroxyl groups: 8

Number of carboxylic acid groups: 1

This monodisperse compound has multiple surface hydroxyl groups and a single carboxylic acid focal point function. The carboxylic acid function is well suited for EDC-mediated coupling to amine-containing reagents. The hydroxyl groups are selectively reactive with anhydrides or acyl halides (see Figure 1). It is transparent to light in the UV/Vis wavelength region.

The product is a freeze-dried, white powder, which may contain traces of methanol or ether.

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

The product is soluble in polar organic solvents (THF, DMSO, DMF, and methanol) and water.

Storage/Stability

Polyester dendrons are susceptible to hydrolysis, which is accelerated when material is dissolved in water, at acidic/basic pH, or at elevated temperatures. For extended storage, store the dry powder at 2–8 °C.

References

EDC-mediated couplings:

1. Desai, M.C., and Stramiello, L.M.S., Polymer Bound EDC (P-EDC): A Convenient Reagent for Formation of an Amide Bond. *Tetrahedron Letters*, **34**, 7685-7866 (1993).
2. Xu, Y., and Miller, M.J., Total Syntheses of Myobactin Analogues as Potent Antimycobacterial Agents Using a Minimal Protecting Group Strategy. *J. Org. Chem.*, **63**, 4314-4322 (1998).
3. Futami, J. et al., Preparation of Potent Cytotoxic Ribonucleases by Cationization. Enhanced Cellular Uptake and Decreased Interaction with Ribonuclease Inhibitor by Chemical Modification of Carboxyl Group. *Biochemistry*, **40**, 7518-7524 (2001).

Related Products:

Polyester-16-hydroxyl-1-carboxyl bis-MPA dendron, generation 4 (Catalog Number 686662)

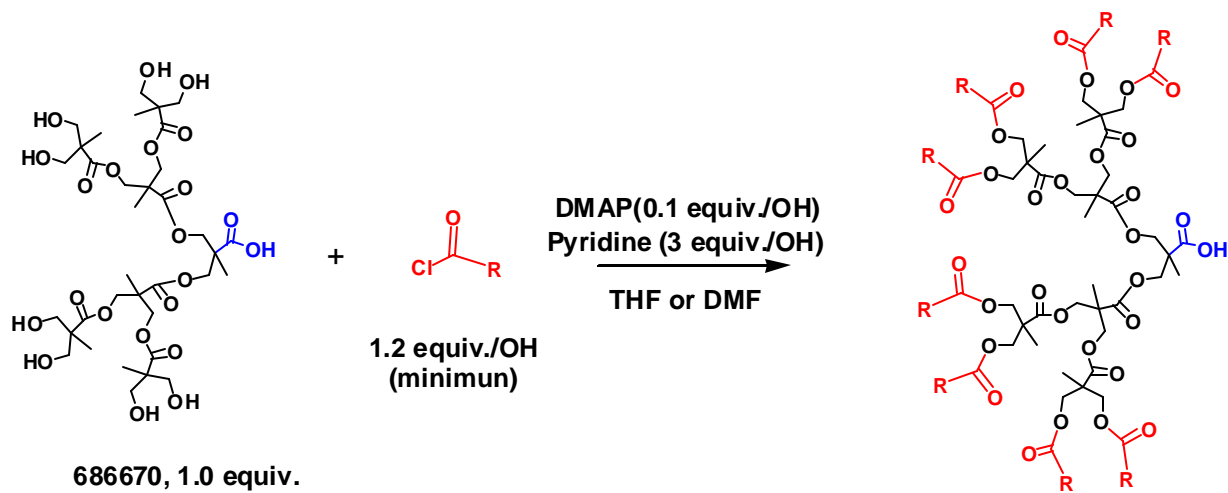
Polyester-32-hydroxyl-1-carboxyl bis-MPA dendron, generation 5 (Catalog Number 686654)

EDC, *N*-(3-Dimethylaminopropyl)-*N*'-ethylcarbodiimide hydrochloride (Catalog Number E1769)

Manufactured by Polymer Factory Sweden AB.

IK,MAM 09/07-1

Figure 1.
Reaction Scheme



Aldrich brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.