

**ALDRICH®**

CHEMISTS HELPING CHEMISTS

Aldrich is a member of the Sigma-Aldrich family

1001 West Saint Paul Ave., Milwaukee, WI 53233  
Telephone: 414-273-3850 • 800-231-8327  
Fax: 414-273-4979 • 800-962-9591  
Website: <http://www.sigma-aldrich.com>  
E-mail: [aldrich@sial.com](mailto:aldrich@sial.com)

## Technical Bulletin

# PATENT NOTICE FOR TFE/PDD COPOLYMER (TEFLON® AF)

**46,961-0 Poly[4,5-difluoro-2,2-bis(trifluoromethyl)-1,3-dioxole-co-tetrafluoroethylene],**  
65 mole % dioxole (TEFLON® AF 1600)

**46,962-9 Poly[4,5-difluoro-2,2-bis(trifluoromethyl)-1,3-dioxole-co-tetrafluoroethylene],**  
87 mole % dioxole (TEFLON® AF 2400)

### Notice of License

Teflon® AF fluoropolymer resins are products of E.I DuPont de Nemours & Company and are protected under one or more of the following U.S. Patents and/or their foreign equivalents: 4,399,264; 4,485,250; 4,754,009; 4,935,477; 5,276,121; 5,326,839; 5,353,368; 5,324,889; 5,338,608; 5,310,838; 5,354,910; 5,408,020; 4,530,569; 4,948,851; 4,975,505; 4,982,056; 5,000,547; 4,977,025; 4,977,026; 4,977,008; 4,977,297; 4,973,142; 4,999,248; 4,985,308; 5,006,382; 5,296,283; 5,308,685; 5,076,659 and 5,051,114. DEPENDING UPON THE APPLICATION, IT MAY BE NECESSARY FOR THE PURCHASER OF THIS RESIN TO OBTAIN A SEPARATE LICENSE FROM DUPONT.

### Restricted Use

The purchase price of Teflon® AF fluoropolymer resin DOES NOT include a royalty for use of the purchased quantity of resin under DuPont's Patents, U.S. or foreign. NO RIGHT OR LICENSE IS GRANTED OR IS TO BE IMPLIED BY THE MERE PURCHASE OF TEFLON® AF FLUOROPOLYMER RESIN ACCORDING TO THIS INVOICE. Purchaser may request a separate, written, non-exclusive license to permit practice under one or more of the above-cited DuPont patents and the payment of royalties separately, such request being directed to DuPont Fluoroproducts, Sales and Marketing Manager--Teflon® AF, P.O. Box 80702, Wilmington, Delaware 19880-0702, U.S.A.