

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352 www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:

APTS (Full chemical name: 8-Aminopyrene-1,3,6-trisulfonic acid, trisodium salt)

CATALOG # 90028

MOLECULAR INFORMATION: C₁₆H₈NNa₃O₉S₃ MWt: 523.39 [196504-57-1]



PROPERTIES:

Color & Form Solubility Absorption/Emission Extinction Coefficient STORAGE AND HANDLING:	Yellow solid Soluble in water 424nm/505nm (before conjugation) 17,000 Store at 4 °C. Protect from light, especially when in solution.
APPLICATION:	APTS is a very useful green fluorescent dye for labeling glycoproteins or sugar molecules in general. The labeling occurs via reductive amination which involves the formation of a Schiff base between the amine of APTS and the aldehyde or ketone of the sugar, followed by the reduction of the Schiff base linkage to a stable carbon- nitrogen bond. The multi-anionic nature of the dye makes it ideal for the studies of carbohydrate molecules by high resolution capillary electrophoresis. Ref.: 1) <i>Anal. Chem.</i> 70 , 568(1998); 2) <i>Electrophoresis</i> , 17 , 347(1996); 3) <i>Electrophoresis</i> , 17 , 412(1996); 4) <i>Anal. Chem.</i> 69 , 4554(1997); 5) <i>Anal. Chem.</i> 69 , 3846(1997).
TOXICITY:	Not established. Not listed by NTP, IARC or OSHA.
FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: *Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.*