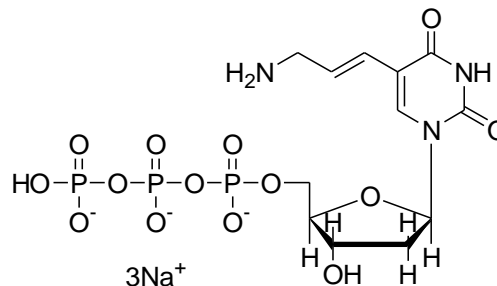


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Aminoallyl-dUTP, sodium salt, lyophilized powder

CATALOG # 40020-1

MOLECULAR INFORMATION: C₁₂H₁₇N₃Na₃O₁₄P₃
MWt: 589.17



PROPERTIES:

Color & Form White solid
Purity ≥ 95% by HPLC
Solubility Soluble in water

STORAGE AND HANDLING: Stored desiccated at -20 °C.

APPLICATION: 5-Aminoallyl-dUTP can be enzymatically incorporated into DNA. The resulting amine-containing DNA can be subsequently labeled with a fluorescent dye, biotin or other haptens via conventional peptide coupling method.^{1,2} This two-step method for labeling nucleic acids is considerably more economical than the one-step method using a pre-labeled dUTP.³ Note: for PCR applications, Taq polymerase should be used with dUTP conjugates, because dUTP inhibits archaeal polymerases such as Pfu and Vent.^{4,5}

Ref.: 1) *Biol Chem Hoppe Seyler* **371**, 953(1990) ; 2) *Biol Chem Hoppe Seyler* **371**, 953(1990) ; 3) *Biotechniques* **28**, 518 (2000); 4) *Anal Biochem.* **211**,164(1993); 5) *PNAS* **99**,596(2002).

TOXICITY: Unknown

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: <i>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.</i>
--