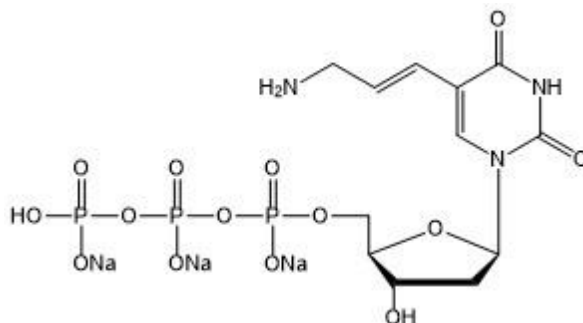


## Datasheet for 5-Aminoallyl-dUTP (AA-dUTP), 10 mM solution

**Catalog number:** C401A

**Product:** 5-Aminoallyl-dUTP (AA-dUTP), 10 mM solution



**Description:** 5-Aminoallyl-dUTP (5-[3-aminoallyl]-2'-deoxyuridine-5'-triphosphate) (cat. no. C401A) is supplied as 10 mM solutions in TE buffer. 5-Aminoallyl-dUTP is also available in lyophilized form (cat. no. C401B).

5-Aminoallyl-dUTP can be enzymatically incorporated into DNA with Reverse Transcriptases, Taq DNA polymerase, phi29 DNA Polymerase, Klenow Fragment, Klenow Fragment, exo- and DNA Polymerase I. The resulting amine-containing DNA can be subsequently labeled with any amine-reactive fluorescent dye, biotin or hapten. This two-step method for labeling nucleic acids is considerably more economical than the one-step method using a prelabeled dUTP.

**Concentration:** 10 mM solution in TE buffer

**Storage conditions:** -20°C

**Molecular formula:** C<sub>12</sub>H<sub>17</sub>N<sub>3</sub>Na<sub>3</sub>O<sub>14</sub>P<sub>3</sub>

**Molecular weight:** 589.17

**Citation of product:** If use of this item results in a publication, please use this information: 5-Aminoallyl-dUTP (AA-dUTP), 10 mM solution (C401A; GeneCopoeia, Inc., Rockville, MD).

**Limited Use License**

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Copyright ©2015 GeneCopoeia, Inc.

LNC401A -DS-031815

**GeneCopoeia, Inc.**

9620 Medical Center Drive, #101

Rockville, MD 20850 USA

Tel: 301-762-0888; Fax: 301-762-3888

Email: [support@genecopoeia.com](mailto:support@genecopoeia.com)

Web: [www.genecopoeia.com](http://www.genecopoeia.com)