

Cyanine 5-dATP [Cy5-dATP]Catalog number: 17038
Unit size: 25 nmoles**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	1182.05
Soluble In	Water

Spectral Properties

Excitation Wavelength	651 nm
Emission Wavelength	670 nm

Applications

The dye-modified deoxyadenosine 5'-triphosphates are widely used for a variety of non-radioactive DNA labeling reactions including nick translation, random prime labeling, cDNA labeling and 3'-end labeling. Cy5-dATP can be used to produce Cy5-containing DNA by conventional enzymatic incorporation methods such as reverse transcription, nick translation, random primed labeling, or PCR. This enzymatic fluorescence labeling method is widely used for both FISH probes and microarray-based experiments. This Cy5-dATP conjugate can be used as a red fluorescence color with SpectrumFRed™ filter set (SpectrumFRed™ is the trademark of Vysis). It is functionally equivalent to Cyanine 5-dATP (Perkin Elmer's NEL593001EA) with significantly improved stability.