

Cy5-11-dCTP *1 mM in Tris Buffer (pH 7.5)*

Catalog Number: 17121

Unit Size: 25 nmoles

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	24 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Liquid
------------	--------

Molecular Weight	1159.01
------------------	---------

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	651 nm
-----------------------	--------

Emission Wavelength	670 nm
---------------------	--------

Applications

Cy5-11-dCTP is a red-fluorescent deoxycytidine analog designed for direct incorporation into DNA using standard enzymatic methods. It can be used as a substitute for dCTP in applications such as PCR, nick translation, and random priming to generate Cy5-labeled DNA or cDNA probes.

The Cy5 dye is linked to the cytidine base and optimized for efficient enzymatic incorporation while preserving polymerase activity. Cy5-11-dCTP is ideal for fluorescence in situ hybridization (FISH), microarray analysis, and other hybridization-based detection platforms requiring far-red readouts.