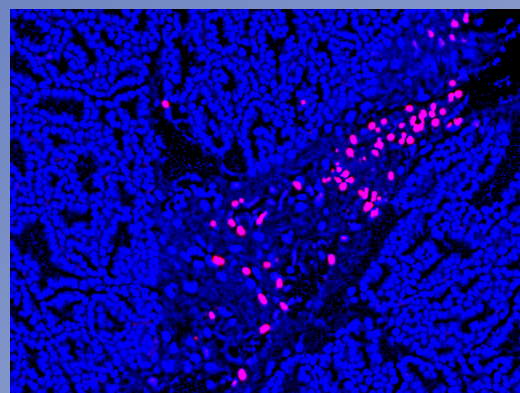
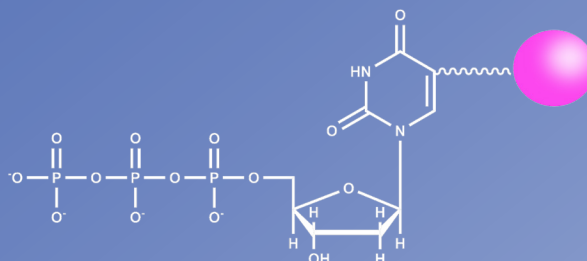


Fluorescent Nucleotide Conjugates and Related Probes

- Labeled dUTP, UTP, dCTP, CTP and ddCTP for nucleic acid probe labeling or TUNEL assay
- Choose from a selection of super bright and exceptionally photostable CF™ dyes, or biotin
- Aminoallyl dUTP and aminoallyl UTP for economical two-step labeling of nucleotides
- BrdU, 5-Bromo-dUTP, and BrUTP for antibody-based detection and cell proliferation studies
- TUNEL kits for labeling apoptotic cells with green fluorescent CF™488A-dUTP, deep red fluorescent CF™594-dUTP, or far-red fluorescent CF™640R-dUTP



TUNEL detection of apoptotic cells in involuting mammary gland using CF594 dUTP (red). Nuclei are counterstained with DAPI (blue).

Catalog No	Product Name	Unit size
40001	5-Tetramethylrhodamine-dUTP, 1 mM in Tris buffer	25 uL
40004	CF405S-dUTP, lyophilized powder	25 nmol
40008	CF488A-dUTP, lyophilized powder	25 nmol
40002	CF543-dUTP, lyophilized powder	25 nmol
40005	CF568-dUTP, lyophilized powder	25 nmol
40006	CF594-dUTP, lyophilized powder	25 nmol
40007	CF640R-dUTP, lyophilized powder	25 nmol
40003	CF680R-dUTP, lyophilized powder	25 nmol
40057	CF532-dCTP, lyophilized powder	25 nmol
40058	CF543-dCTP, lyophilized powder	25 nmol
40027	CF555-dCTP, lyophilized powder	25 nmol
40055	CF568-dCTP, lyophilized powder	25 nmol
40056	CF594-dCTP, lyophilized powder	25 nmol
40028	CF647-dCTP, lyophilized powder	25 nmol
40031	CF555-ddCTP, lyophilized powder	25 nmol
40032	CF640R-UTP, lyophilized powder	25 nmol
30063	CF488A TUNEL Assay Apoptosis Detection Kit	50 reactions
30064	CF594 TUNEL Assay Apoptosis Detection Kit	50 reactions
30074	CF640R TUNEL Assay Apoptosis Detection Kit	50 reactions

Catalog No	Product Name	Unit size
40020	5-Aminoallyl-dUTP, 10 mM in TE buffer	100 uL
40020-1	5-Aminoallyl-dUTP, lyophilized powder	1 mg
40021	5-Aminoallyl-UTP, 10 mM in TE buffer	100 uL
40021-1	5-Aminoallyl-UTP, lyophilized powder	1 mg
40022	Biotin-16-dUTP, 1 mM in Tris buffer	50 uL
40022-1	Biotin-16-dUTP, lyophilized powder	50 ug
40023	Biotin-16-UTP, 10 mM in Tris buffer	25 uL
40024	5-Bromo-2'-deoxyuridine (BrdU)	100 mg
40025	5-Bromo-dUTP, 10 mM in TE buffer	25 uL
40026	5-Bromouridine-5'-triphosphate (BrUTP), 10 mM in TE buffer	25 uL
40029	Biotin-11-dUTP, 1 mM in Tris buffer	50 uL
40029-1	Biotin-11-dUTP, lyophilized powder	50 ug
40030	Biotin-20-dUTP, 1 mM in Tris buffer	50 uL
40030-1	Biotin-20-dUTP, lyophilized powder	50 ug
40033	Biotin-11-UTP, 10 mM in Tris buffer	25 uL
40034	Biotin-20-UTP, 10 mM in Tris buffer	25 uL
40035	Biotin-11-CTP, 10 mM in Tris buffer	25 uL
40036	Biotin-11-dCTP, 1 mM in Tris buffer	50 uL

Also see our selection of CF™ dye-labeled monoclonal anti-biotin antibodies and streptavidin conjugates for detection of biotinylated probes.