



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## TUNEL Assay Apoptosis Detection Kit

The CF® dye TUNEL Assay Apoptosis Detection Kit label apoptotic cells in fixed tissue sections or cellular samples for fluorescence microscopy or flow cytometry.



### Product Description

These kits use enzymatic labeling of DNA double-strand breaks for fluorescence staining of apoptotic cells in fixed samples or tissue sections by TUNEL.

In TUNEL (terminal deoxynucleotidyl transferase-mediated dUTP nick-end labeling), TdT enzyme catalyzes the addition of labeled dUTP to the 3'- ends of the internucleosomal DNA strand breaks that are a hallmark of apoptosis. The CF® dye TUNEL Assay Apoptosis Detection Kits employs dUTP conjugated to the exceptionally bright and photostable fluorescent dyes, for single step fluorescent TUNEL labeling of fixed cells or tissues for analysis by fluorescence microscopy or flow cytometry. Learn more about Biotium's next-generation [CF® Dyes](#).

See our full selection of [Cell Viability & Apoptosis Kits](#).

### CF® Dye TUNEL Assay Apoptosis Detection Kits

Dye	Ex/Em	Size	Catalog No.
CF®488A	490/515 nm	50 reactions	<a href="#">30063</a>
CF®594	593/614 nm	50 reactions	<a href="#">30064</a>
CF®640R	642/662 nm	50 reactions	<a href="#">30074</a>

Kit sized based on 50 uL reaction volume.

### Product attributes

<b>Apoptosis/viability marker</b>	DNA fragmentation
<b>For live or fixed cells</b>	For fixed cells/tissue sections
<b>Detection method/readout</b>	Fluorescence microscopy, Flow cytometry
<b>Assay type/options</b>	Endpoint assay
<b>Fixation options</b>	Fix before staining (formaldehyde), Fix before staining (methanol), Permeabilize before staining
<b>Colors</b>	Green, Red, Far-red
<b>Product origin</b>	TdT enzyme (calf thymus); recombinant, produced in E. coli
<b>Storage Conditions</b>	Store at -10 to -35 °C, Protect from light

This datasheet was generated on January 6, 2026 at 11:53:15 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cf-dye-tunel-assay-apoptosis-detection-kit/>