Glowing products for science

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	30063
CF®594	593/614 nm	50 reactions	<u>30064</u>
CF®640R	642/662 nm	50 reactions	<u>30074</u>

Kit sized based on 50 uL reaction volume.

This datasheet was generated on November 4, 2025 at 04:27:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cf-dye-tunel-assay-apoptosis-detection-kit/

Product attributes

Call us: 800-304-5357

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

Email: btinfo@biotium.com