

# Cell Counting Kit-F

Fluorometric kit for simple and accurate cell proliferation and cytotoxicity assays. **Advantages:** No radioisotopes, more sensitive than the colorimetric assay, short staining reaction time, no solubilization step, correlates with the [<sup>3</sup>H]-thymidine incorporation assay.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |             |
|------------------|-------------|
| ALX-850-245-KI01 | 500 tests   |
| ALX-850-245-KI02 | 5x500 tests |

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Protect from light.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CCKi-F

**Application** Fluorescence microscopy

**Technical Info / Product Notes** **Principle:** The amount of a fluorescent dye, calcein, produced from calcein-AM (3',6'-Di(O-acetyl)-2',7'-bis[N,N-bis-(carboxymethyl)aminomethyl]fluorescein . tetraacetoxymethyl ester) by esterases in cells is directly proportional to the number of viable cells in a culture medium. Since calcein-AM is highly lipophilic due to the acetoxymethyl groups, it can rapidly permeate into the cytoplasm through the cell membrane. The detection range of the number of viable cells by CCKi-F is from <50 to at least 25,000 cells.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)