

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 [³H]-thymidine incorporation assay.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-2450	500 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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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