

Cell Counting Kit-F (500 tests)

ALX-850-245-KI01: ~500 tests

ALX-850-245-KI02: 5x ~500 tests

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

TABLE OF CONTENTS

Introduction	2
Materials Supplied	2
Storage	2
Other Materials Needed	2
Working Solution	2
Cytotoxicity Assay Procedure	3
Determination of IC ₅₀	3
Caution.....	3
Typical Results.....	4
Contact Information.....	6

INTRODUCTION

Cell Counting Kit-F (CCK-F) offers simplified and highly sensitive cell proliferation and cytotoxicity assay method by employing Calcein-AM (3',6'-Di(O-acetyl)-4',5'-bis[N,N-bis(carboxymethyl)amino-methyl]fluorescein, tetraacetoxymethyl ester) that produces a highly sensitive fluorescent dye ($\lambda_{\text{ex}}=490 \text{ nm}$, $\lambda_{\text{em}}=515 \text{ nm}$) upon enzymatic hydrolysis by esterases in living cells. The amount of the fluorescent dye, calcein, produced by hydrolysis by esterases in cells is directly proportional to the number of viable cells in a culture medium.

The CCK-F assay has a detection range from 50 or fewer to at least 25,000 cells. Furthermore, CCK-F assay does not require any radioisotopes (such as in [^3H]-thymidine incorporation assay) nor solubilization procedure (such as in MTT assay), and no special skills are necessary for use. Therefore, it allows you to obtain highly reproducible and accurate results.

MATERIALS SUPPLIED

Calcein-AM DMSO solution, 110 $\mu\text{l} \times 1$

STORAGE

Store at -20°C . This product is stable for six months at -20°C before opening.

If the solution remains, close the bottle cap tightly and store in a freezer at -20°C .

OTHER MATERIALS NEEDED

1. Fluorescence microplate reader (excitation filter: 480-500 nm, emission filter: 500-535 nm)
2. 96-well microplate for fluorescent measurements (black plate or white plate)
3. Phosphate Buffered Saline (PBS)

WORKING SOLUTION

Required amount of Calcein-AM DMSO solution is diluted 50 times with PBS. Since Calcein-AM is not stable in PBS, prepare the working solution immediately before use.



Reagents require separate storage conditions.

CYTOTOXICITY ASSAY PROCEDURE

1. Dispense 100 µl of cell suspension (5,000 cells/well) onto a 96-well microtiter plate.
2. Preincubate the plate for 24 hours in an incubator (humidified atmosphere, e.g., at 37°C, 5% CO₂).
3. Add 10 µl of various concentrations of the test substance into the culture medium of the plate.
4. Incubate cell cultures for 48 hours in the incubator.
5. Replace the medium with 100 µl of PBS. For nonadherent cells, centrifuge the microplate before removing the culture medium.
6. Add 10 µl of the working solution and incubate cell cultures for 15-30 min in the incubator.
7. Measure the fluorescence ($\lambda_{ex}=490$ nm, $\lambda_{em}=515$ nm) using a fluorescence microplate reader.

DETERMINATION OF IC₅₀

Cell viability is calculated using the following equation. IC₅₀, concentration killing 50% of the cells, is determined from plot of viability versus concentration of the test substance.

$$\text{Viability (\%)} = [(A_s - A_b) / (A_c - A_b)] \times 100$$

A_s: Fluorescence intensity of sample (cell + test substance + CCK-F)

A_c: Fluorescence intensity of control (cell + CCK-F, no test substance)

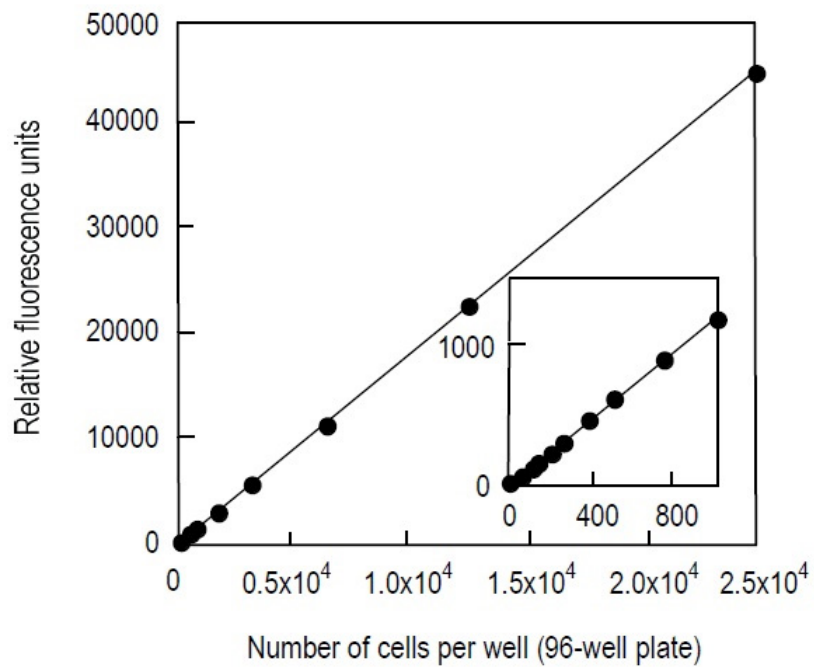
A_b: Fluorescence intensity of blank (medium + CCK-F, no cell)

Example of toxicological test using CCK-F is shown in Fig. 3.

CAUTION

1. Phenol red and serum in a culture medium interfere with the fluorescence measurement. Replace a culture medium with PBS or phenol red and serum free medium prior to adding the working solution.
2. Dilute the test substance with non-toxic solution, such as culture medium or PBS or saline. And use same solution for control and blank wells instead of sample.
3. If a 24-well or 6-well plate is used for this assay, calculate the number of cells per well accordingly, and adjust the volume of the working solution in a well to 10% of the total volume.

TYPICAL RESULTS



Cell line: HL60

Incubation time: 30 min

Detection: $\lambda_{ex}=485$ nm, $\lambda_{em}=535$ nm

Fig.1 A relationship between fluorescence intensity and number

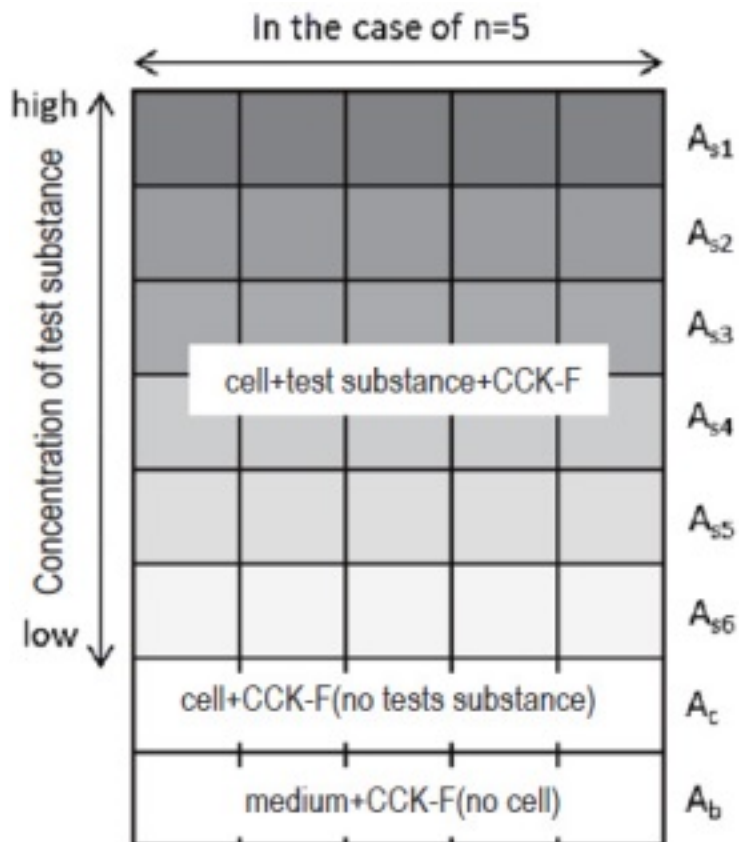
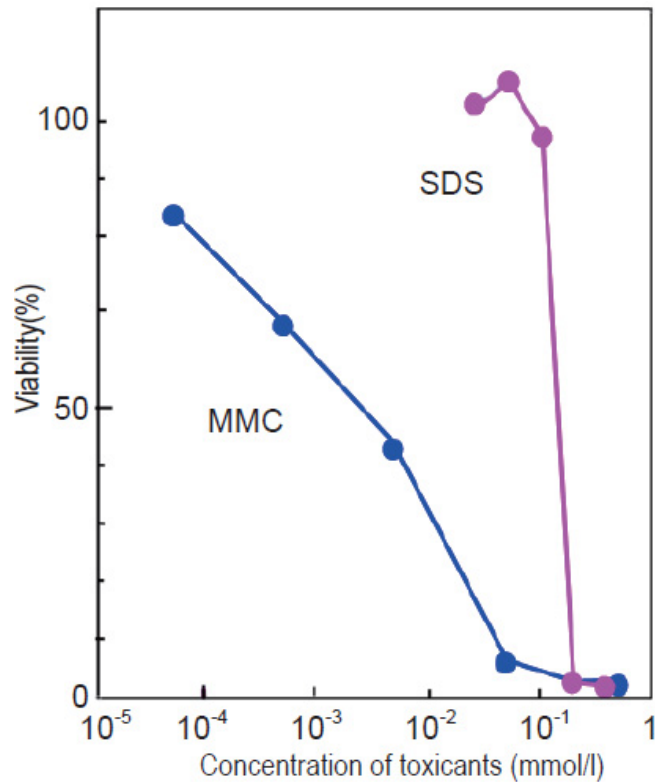


Fig. 2 An example of sample arrangement



Cell line: HL60

Toxicants: Mitomycin C (MMC), Sodium dodecylsulfate (SDS)

Condition: 37°C, 5% CO₂, 48 hrs

Incubation time: 30 min

Detection: λ_{ex} =485 nm, λ_{em} =535 nm

Fig.3 Toxicological test for MMC and SDS by CCK-F.



Product Manual

www.enzolifesciences.com

Enabling Discovery in Life Science®

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

www.enzolifesciences.com