

EZ4U (NONRADIOACTIVE CELL PROLIFERATION AND CYTOTOXICITY)

ELISA KIT

NONRADIOACTIVE CELL PROLIFERATION AND CYTOTOXICITY ASSAY.
CAT. NO. BI-5000 10 X 96 DETERMINATIONS

FOR RESEARCH USE ONLY
NOT FOR USE IN DIAGNOSTIC PROCEDURES

rev.no. 301004 (replacing 020604)

Biomedica Medizinprodukte GmbH & Co KG, A-1210 Wien, Divischgasse 4
Tel. +43/1/291 07 45, Fax +43/1/291 07 85, E-mail export@bmgmp.com



CONTENT

1) ENGLISH 3

Additional information on our products is available on our website.

www.bmgrp.com

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES

1) INTRODUCTION

Proliferation assays are widely used in cell biology for the study of growth factors, cytokines, nutrients and for the screening of cytotoxic or chemotherapeutic agents. There are several ways to determine the number of cells either by microscopic inspection, or by the use of an electronic particle counter, indirectly by measuring the incorporation of radioactive precursors, quantitating total protein with chromogenic dyes, or by measuring metabolic activity of cellular enzymes. The most common assay for cell proliferation is the incorporation of ³H-thymidine into cellular DNA. The ³H-thymidine assay is, however, labour intensive as it requires the removal of excess, unincorporated label by using some method of cell harvesting before measurement. In 1956, the first paper was published on the use of tetrazolium salts as indicators of cell viability. The method was based on the finding that living cells are capable to reduce slightly or uncoloured tetrazolium salts into intensely coloured formazan derivatives. This reduction process requires functional mitochondria, which are inactivated within a few minutes after cell death.

This method therefore provides an excellent tool for the discrimination of living and death cells. However, the early tetrazolium salts did have some disadvantages, such as the insolubility of the resulting formazan products. Time and labour consuming resolubilisation procedures were necessary, including repipetting and mixing, or the application of hazardous solubilisers. This necessary post assay treatment, however, irreversibly terminated cell proliferation and thus made it impossible to prolong incubation in order to achieve an increase in sensitivity or continue cell culture. These inconveniences led to the development of non-toxic tetrazolium salts which yield soluble reduction products. Although the assay procedure was made easier by these soluble dyes, in practice the use was limited due to the instability of the formazan dye and a relatively low absorbance of the end product as compared to the classical MTT assay.

The BIOMEDICA research department has solved both problems and created an easy to use, rapid and reliable non-isotopic cell proliferation assay. For convenience, we have made it highly compatible with the standard thymidine incorporation assay. Therefore, no changes are required in the setup of the test and in the "labelling" procedure. Furthermore, there is no need for the removal of culture medium before or after the addition of the chromogenic substrate and neither solubilisation nor harvesting procedures are necessary. The work performed by BIOMEDICA resulted in an assay which combines the best of the thymidine and MTT methods, namely: accuracy, speed, reliability and ease of use. Also, according to our data achieved so far, the chromophore appears to be non-toxic. A double labelling with EZ4U and a radioactive nucleotide to obtain more information about cell viability and DNA content is now feasible.

2) CONTENTS OF THE KIT

CONT	KIT COMPONENTS	QUANTITY
SUB	Substrate, lyophilised	10 vials
ACT	Activator solution, ready to use	1 x 30 ml
	Package insert	1 piece

3) MATERIAL AND EQUIPMENT REQUIRED BUT NOT SUPPLIED

Microtiterplate reader equipped with a 450 nm or 492 nm filter (see Fig.1). We recommend the use of a 620 nm to 690 nm reference wavelength, which is beneficial but not absolutely necessary.

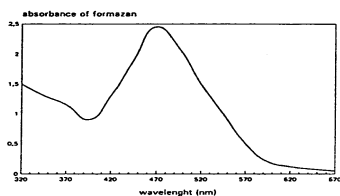


Fig. 1. Absorption spectrum of Formazan.

4) REAGENTS AND SAMPLE PREPARATION

One vial SUB in this EZ4U (Nonradioactive Cell Proliferation and Cytotoxicity) ELISA Assay kit contains the amount of substrate sufficient for one 96 well plate with 200 µl of cell culture medium/well. If more substrate is needed, combine the dissolved substrates prior to pulsing.

Dissolve the SUB (substrate) in 2.5 ml ACT (activator). Prewarm this solution to 37°C prior to addition. If necessary, warm up the substrate vial in your hand while mixing with activator.

This procedure yields a straw-coloured solution.

The mixed substrate is designed for immediate use only and should not be stored.

5) GENERAL CONSIDERATIONS FOR SETTING UP ASSAYS WITH EZ4U

The assay set-up is performed in a manner similar to the standard 3H-thymidine incorporation method. Instead of pulsing with tritiated nucleotide, 20 µl of dye solution is added to 200 µl sample. Incubation time is dependent on the metabolic capacity of the cells. Usually 2 to 5 hours of incubation at 37°C are sufficient to yield a significant increase in color intensity. As different cells vary in their ability to convert the yellow colored tetrazolium compound to its red formazan derivative, we recommend testing every new cell-line's metabolic capacity as described in Fig.2. After incubation, the plate is removed from the incubator and gently mixed by tipping the plate at all four sides. To avoid increased standard deviations, the plate must be shaken before reading the optical density.

The absorbance is measured by a microplate-reader, set at 450 nm or 492 nm with 620 nm as a reference. The reference absorbance at 620 nm (or any wavelength between 620-690 nm) is used to correct for nonspecific background values, caused by cell debris, fingerprints, or other potential interferences. However, the reference may be omitted without significant changes in the accuracy of the assay.

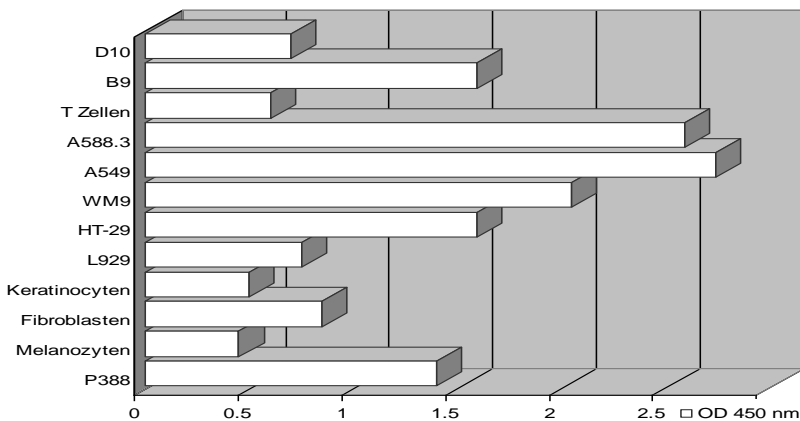


Fig. 2. Different metabolic capacity of various cell lines.

3×10^3 cells/well were cultivated in 200 µl RPMI 1640. Following a cultivation period of 3 days, 25 µl of the dye substrate were added to each well. Optical density was recorded after 4 hours, showing significant differences in the metabolic capacity of the various cell lines.

6) IMPORTANT CONSIDERATIONS

- The substrate is not sterile. If sterile conditions are demanded, the solubilised ready to use dye substrate must be sterile filtered. (A minor turbidity prior filtration interferes neither with the filtration, nor with the assay performance).
- Due to the high sensitivity of this test, it is advisable to use as little cells as possible. Otherwise the occurrence of a non linear titration-curve may be possible.
- To achieve reproducible time kinetics in color development, equilibrate cell cultures at 37°C.
- Do not prolong incubation times without pretesting, this might result in an increased background without improved sensitivity.
- The use of a reference wavelength of 620 nm (which is subtracted from the values obtained at 450 or 492 nm) is not absolutely necessary, but increases the performance of the test.
- The chromophore appears to be non-toxic and therefore prolongation of cell culture is possible after removal of the formazane derivative.

7) ASSAY PROTOCOL

All reagents and samples in this EZ4U, (Nonradioactive Cell Proliferation and Cytotoxicity) ELISA Assay kit must be at room temperature (18-26°C) before use in the assay.

Add 200 µl cell culture into respective wells.

Add 20 µl SUB (substrate) into each well, swirl gently.

Incubate at 37°C for 2-5 hours, depending on the metabolic capacity of the cells.

If no microplate reader with shaking plate carrier is available, mix the plate on a vibrating platform or by tipping with the fingers.

Read absorbance at 450 nm or 492 nm, with 620 nm as reference.

For the most accurate results, absorbance from a substrate blank in assay medium without cells should be subtracted from all other values.

8) TECHNICAL HINTS

- Do not mix or substitute reagents with those from other lots or sources.
- Do not mix stoppers and caps from different reagents or use reagents between lots.
- Do not use reagents beyond expiration date.
- Protect reagents from direct sunlight.
- Avoid foaming when mixing reagents.

9) PRECAUTIONS

- Do not pipette by mouth.
- Do not eat, drink, smoke or apply cosmetics where reagents are used.
- Avoid all contact with the reagents by using gloves.

10) LITERATURE

- Hansen M. B., Nielsen S.E. and Berg K., J. Immunol. Meth. 119, 203-210, (1989)
- Denizoit F. and Lang R., J. Immunol. Meth. 89, 271-277, (1986)
- Carmichael J. et al., Cancer Res., 47, 943-946, (1987)
- Karasek M; Gruszka. Et al. Department of Electron Microscopy, J Pineal Res 2003 May;34(4):294-6
- Klocking R. et al; Institute for Antiviral Chemotherapy; Antivir Chem Chemother 2002 13(4):241-9;
- Wutzler P. et al.; Institute for Antiviral Chemotherapy; Antiviral Res 2002 May;54(2):89-97
- Hutter R. et al., Circulation, Apr 2003; 107: 1658-1663.
- Sturlan S. et al., Blood, Jun 2003; 101: 4990-4997.
- Völlenkle Ch. et al., Appl Envir Microbiol., Mar 2004; 70: 1514-1521.

- Bauer S. et al., J Immunol., Mar 2004; 172: 3930-3939.
- Thurow K. et al., Journal of the Association for Laboratory Automation, Jun 2004; 9: 159-162.
- Heffeter P. et al., J Pharmacol Exp Ther., Jan 2005; 312: 281-289.
- Kuhl A. et al., Antimicrob Agents Chemother., Mar 2005; 49: 987-995.
- Sulochana KN. et al., J Biol Chem, Jul 2005; 280: 27935-27948.
- Ank N. et al., J. Virol., Aug 2005; 79: 9831-9841.
- Wegner B. et al., Nephrol Dial Transplant., Oct 2005; 20: 2071-2079.
- Fineschi S. et al., FASEB J, Jan 2006; 10.1096/fj.05-4870fje.
- Perabo F. et al., Anticancer Res, May 2006; 26: 2129-2135.
- Bauer S. et al., J. Immunol., Aug 2006; 177: 2423-2430.
- Wolf AM et al., Haematologica, Sep 2006; 91: 1165-1171.
- Yoshida A. et al., Anticancer Res, Nov 2006; 26: 4003-4007.
- Molnarfi F. et al., J Immunol., Jan 2007; 178: 446-454.
- Demyanets S. et al., Am J Physiol Heart Circ Physiol, Sep 2007; 293: H1962-H1968.
- Tonack S. et al., Endocrinology, Dec 2007; 148: 5902-5912.
- Sonvilla G. et al., Carcinogenesis, Jan 2008; 29: 15-24.
- Hohegger K. et al., J Am Soc Nephrol., Aug 2008; 19: 1520-1529.
- Stewart-Jones G. et al., PNAS, Apr 2009; 106: 5784-5788.
- Walochnik J. et al., J Antimicrob Chemother., Sep 2009; 64: 539-545.
- Stuhlmeier K. et al., Exp Biol Med., Nov 2009; 234: 1327-1338.
- Ang ChW. et al., Carcinogenesis, Sep 2010; 31: 1541-1551.
- Shehata M. et al., Blood, Oct 2010; 116: 2513-2521.
- Cindric M. et al., Anticancer Res., Oct 2010; 30: 4063-4069.
- Gomez A. et al., Mol Pharmacol., Dec 2010; 78: 1004-1011.
- Breinig M. et al., Clin Cancer Res., Apr 2011; 17: 2237-2249.
- Jungwirth U. et al., Mol Pharmacol., May 2012; 81: 719-728.
- Swoboda A. et al., Hum Mol Genet., Aug 2012; 21: 3387-3396.
- Cao R. et al., PNAS, Sep 2012; 109: 15894-15899.

SYMBOLS



Expiry date / Verfallsdatum / Date de péremption / Data di scadenza / Fecha de caducidad / Data de validade / Uiterste gebruiksdatum / Udløbsdato / Utgångsdatum / Termin Ważności / Lejárati idő / Doba expirácie / Doba expirace



Consider instructions for use / Bitte Gebrauchsanweisung beachten / Consultez la notice d'utilisation / Consultare le istruzioni per l'uso / Consulte las instrucciones de utilización / Consulte as instruções de utilização / Raadpleeg de gebruiksaanwijzing / Se brugsanvisningen / Läs anvisningarna före användning / Proszę przeczytać instrukcję wykonania / Vegyük figyelembe a használati utasításban foglaltakat / Postupujte podľa pokynov na použitie / Postupujte dle návodu k použití



In vitro Diagnostic Medical Device (for in Vitro Diagnostic Use) / In vitro Diagnostikum (zur In-vitro-Diagnostik) / Dispositif médical de diagnostic in vitro (Pour usage diagnostique in vitro) / Dispositivo medico per diagnostica in vitro (per uso diagnostico in vitro) / Dispositivo médico de diagnóstico in vitro (para uso diagnóstico in vitro) / Dispositivo médico para diagnóstico in vitro (Para utilização de diagnóstico "in vitro") / Medisch hulpmiddel voor diagnostiek in vitro (Voor diagnostisch gebruik in vitro) / Medicinsk udstyr til in vitro-diagnostik (Udelukkende til in vitro diagnostisk anvendelse) / Medicinteknisk produkt avsedd för in vitro-diagnostik (För in vitro-diagnostiskt bruk) / Wyrób medyczny do Diagnostyki In Vitro / In vitro orvosdiagnosztikai termék / In vitro diagnostický zdravotnícký materiál (určené pre diagnostiku „in vitro“) / In vitro diagnostický zdravotnícký materiál (určeno pro diagnostiku „in vitro“)



Lot-Batch Number / Charge-Chargennummer / Lot-Code du lot / Lotto-Numero di lotto / Lote-Código de lote / Lote-Código do lote / Lot-Partijnummer / Lot-Batchkode / Lot-Satskod / Numer serii / Lot-Batch szám / Číslo šarže / Číslo šarže



Manufactured by / Hergestellt von / Fabriqué par / Prodotto da / Fabricado por / Fabricado por / Vervaardigd door / Fabrikation af / Tillverkad av / Wyprodukowane pr / Gyártotta / Vyrobené / Vyrobeno



Catalogue Number / Bestellnummer / Numéro de référence / Numero di riferimento / Número de referencia / Número de referência / Referentienummer / Referencenummer / Katalognummer / Numer katalogowy / Katalógusszám / Katalógové číslo / Katalogové číslo



Store at between / Lagerung bei zwischen / Conserver à entre / Conservare a tra / Conservar a temp. entre / Armazene a entre / Bewaar bij tussen / Opbevares mellem / Förvaras vid / Przechowywać w / Táróljuk között / Skladujte v rozsahu / Skladujte v rozmezí



Contains sufficient for x tests / Inhalt ausreichend für x Tests / Contient suffisant pour x tests / Contenuto sufficiente per x test / Contiene suficiente para x pruebas / Contém suficiente para x testes / Bevat voldoende voor x bepalingen / Indeholder tilstrækkeligt til x prøver / Innehållet räcker till x analyser / Zawartość na x testów / Tartalma X teszt elvégzésére elegendő / Obsahuje materiál pre x testov / Obsahuje materiál pro x testů

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.



Manufactured by:

Biomedica Medizinprodukte GmbH & Co KG
A-1210 Wien, Divischgasse 4
Tel.+43/1/291 0745 • Fax +43/1/291 0785
www.bmgrp.com • export@bmgrp.com

distributed in the US/Canada by:



EAGLE
BIOSCIENCES

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com • info@eaglebio.com