

MDR Assay Kit (Fluorometric) ab112142

6 References 1 图像

概述

产品名称	MDR Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
检测时间	1h 0m
种属反应性	与反应: Mammals, Other species
产品概述	<p>Multi drug resistance (MDR) is a major factor in the failure of many forms of chemotherapy. In the past few years it has become widely accepted that the resistance to chemotherapy correlates with the overexpression of at least two ATP dependent drug efflux pumps. These cell membrane proteins, called P-glycoprotein (Pgp, MDR1), and multidrug resistance associated protein (MRP1) are members of the ABC transporter family. Abcam's Fluorimetric MDR Assay Kits use a fluorescent MDR indicator for assaying these two MDR pump activities. This hydrophobic fluorescent dye molecule rapidly penetrates cell membranes and becomes trapped in cells. Following a short incubation, the intracellular free dye concentration can increase significantly. In the MDR1 and/or MRP1-expressing cells this dye is extruded by the MDR transporters, thus decreasing the cellular fluorescence intensity. However, when its extrusion is blocked by an agent that interferes with the MDR1 and/or MRP1 pump-activity, its cellular fluorescence intensity increases significantly. Abcam's Fluorimetric MDR Assay Kits are ideal for high throughput screening of MDR pump inhibitors or identifying the cells that have high level of MDR pump activities.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	Store ab112142 dessicated.
平台	Microplate reader

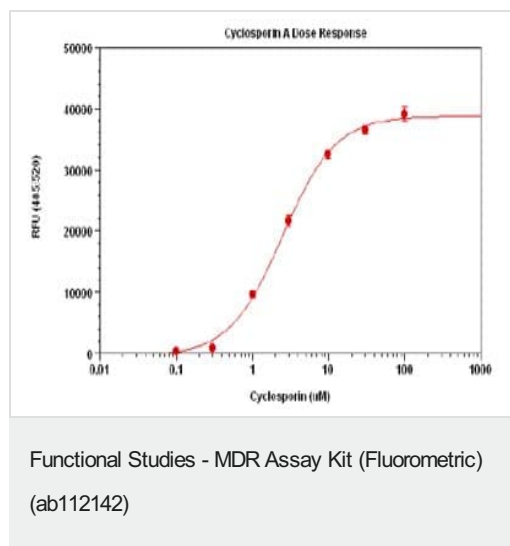
性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests	10 x 96 tests
Assay Buffer	1 x 10ml	1 x 100ml

组件	1 x 96 tests	10 x 96 tests
DMSO	1 x 100μl	1 x 300μl
MDR Sensor	1 vial	1 vial

图片



Effect of Cyclosporin A on the inhibition of P-gp pump in MCF7/ADR cells. The increased concentration of Cyclosporin A resulted in an increase in fluorescence signal caused by the inhibition of P-gp pump which enhanced the intracellular accumulation of MDR indicator dye. The $EC_{50} = 2.4 \mu M$ (measured with the kit) is similar to the value reported in the literature.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors