

SEAP Reporter Gene Assay Kit (Luminescence) ab133077

[2 References](#) [1 图像](#)

概述

产品名称	SEAP Reporter Gene Assay试剂盒(Luminescence)
检测方法	Luminescent
样品类型	Cell culture media
检测类型	Quantitative
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	SEAP Reporter Gene Assay Kit (Luminescence) (ab133077) provides a simple chemiluminescence method for the sensitive quantitation of SEAP (secreted alkaline phosphatase) in conditioned cell culture medium from transfected cells. The assay can detect SEAP activity in the microunits/well (mU/ml) range. The kit includes enough reagents to run three 96-well plates. The assay is easy to perform and can be completed within one hour.

说明

Secreted alkaline phosphatase (SEAP) is commonly used as a reporter of gene expression. Compared to other conventional intracellular reporters such as chloramphenicol acetyltransferase (CAT) and firefly luciferase, SEAP has the advantage of being secreted from transfected cells into the culture medium. SEAP activity in the culture medium is directly proportional to changes in intracellular concentrations of SEAP mRNA and protein. In addition, the kinetics of gene expression can be studied using the same cultures by repeatedly collecting culture medium at different time points. The intact cells can be used for further analysis of RNA or protein expression.

SEAP activity was first measured using the chromogenic alkaline phosphatase substrate *p*-nitrophenyl phosphate (*p*NPP). Today, the most sensitive SEAP assays employ chemiluminescent alkaline phosphatase substrates such as the 1,2-dioxetane CSPD. Chemiluminescent detection of SEAP is fast, easy to perform, and sensitive.

平台 Microplate reader

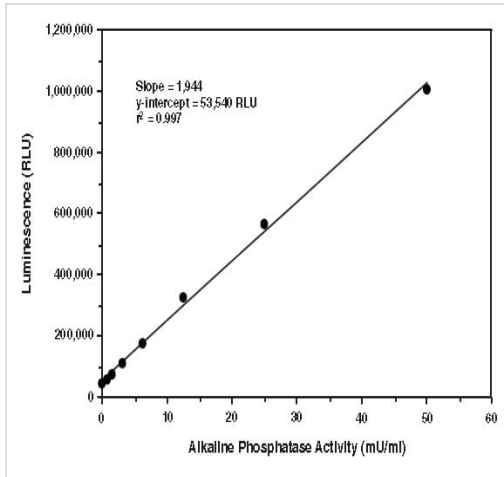
性能

存放说明 Store at +4°C. Please refer to protocols.

组件	300 tests	300 tests
96-Well Solid Plate (white) with lid	3 units	3 units

组件	300 tests	300 tests
Cell-Based Alkaline Phosphatase Standard	1 x 200µl	1 x 200µl
SEAP Substrate (Luminescence)	1 x 15ml	1 x 15ml

图片



Cell-Based Alkaline Phosphatase Standard curve

Functional Studies - SEAP Reporter Gene Assay
Kit (Luminescence) (ab133077)

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