

Product datasheet

PEG Virus Precipitation Kit ab102538

★★★★★ 1 Abreviews 6 References 1 图像

概述

产品名称	PEG Virus Precipitation试剂盒
样品类型	Other biological fluids, Cell culture media
检测类型	Direct
灵敏度	> 40 %
范围	40 % - 100 %
产品概述	<p>Abcam's PEG Virus Precipitation Kit provides an easy, convenient and time-saving method to concentrate viruses without ultra-centrifugation. The kit can be used for small lab samples or large scale virus preparation with high yield and high viral titer. The kit can be used to concentrate retroviruses, baculoviruses, lentiviruses, and phages etc. in cell culture medium or environmental samples. Viruses can be concentrated over 100 fold. An optimized Virus Re-suspension Solution is provided to maximize viral recovery by 40-100% depending on the virus type and sources. The whole process uses non-toxic reagents. The concentrated virus can be used for infection, viral DNA or RNA purification, etc.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<p>Viruses are usually produced at a low titer. They often need to be concentrated for storage or further applications. A quick, easy and inexpensive method is desired to concentrate viruses and remove impurities.</p>

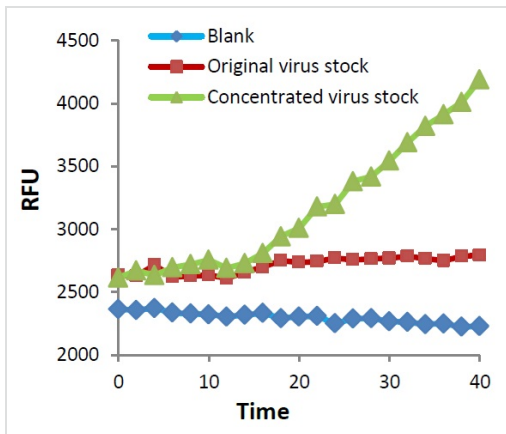
性能

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

组件	50 tests
PEG Solution	1 x 125ml
Virus Re-suspension Solution	1 x 10ml

相关性 Virus is usually produced at a low titer. It often needs to be concentrated for storage or further applications. A quick, easy and inexpensive method is desired to concentrate virus and remove impurities. PEG (Polyethylene glycol) Virus Precipitation provides an easy, convenient and time-saving method to concentrate virus without ultra-centrifugation.

图片



Functional Studies - PEG Virus Precipitation Kit
(ab102538)

Concentration of baculovirus: Low titer baculovirus (10 ml) expressing human Granzyme B was precipitated following the kit protocol and the precipitate was suspended in 1 ml of Virus Re-suspension Solution. Both low titer and precipitated baculovirus was subsequently used to infect SF9 Insect cells. The activity of recombinant Granzyme B secreted into the culture medium by low titer and precipitated baculovirus infected insect cells was monitored using [ab157403](#) - Granzyme B Activity Assay Kit (Fluorometric).

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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