

# PEG Virus Precipitation Kit ab102538

★★★★★ [2 Abreviews](#) [15 References](#) [画像数 1](#)

### 製品の概要

製品名	PEG Virus Precipitation Kit
サンプルの種類	Other biological fluids, Cell culture media
アッセイタイプ	Direct
検出感度	> 40 %
検出範囲	40 % - 100 %
製品の概要	<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 <a href="#">FAQs page</a> for tips and troubleshooting.</p>
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K904 PEG Virus Precipitation Kit. K904-50 is the same size as the 50 test size of ab102538.</p> <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>

### 製品の特性

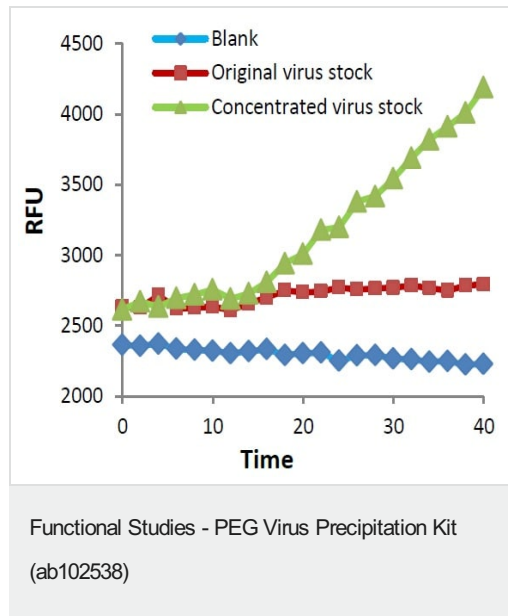
**保存方法** Store at -20°C. Please refer to protocols.

内容	50 tests
5X PEG Solution	1 x 125ml
Virus Resuspension Solution	1 x 10ml

**関連性** 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.

## 画像



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. 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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

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