

# Whole Blood Polymorphonuclear Cell Isolation Kit ab241001

[2 Images](#)

### Overview

<b>Product name</b>	Whole Blood Polymorphonuclear Cell Isolation Kit
<b>Sample type</b>	Whole Blood
<b>Product overview</b>	<p>Whole Blood Polymorphonuclear Cell Isolation Kit (ab241001) is unique in the market because it enables the user to isolate Polymorphonuclear cells (PMNs), assess viability of the obtained cell fractions and also determine purity with three simple, easy-to-use reagents. High yields, or 60-80% of total PMNs have been obtained yielding <math>\geq 5 \times 10^6</math> cells per mL of human whole blood. Purity of the cell fraction is assessed with giemsa stain (included). The kit's use results in highly enriched (&gt;95%) cell fraction of polymorphonuclear cells with less than 3% red blood cell counts. More than 99% of the cells are viable, as determined by the Viability Stain.</p> <p>30 test refers to 30 mL isolation.</p>

<b>Notes</b>	This product is manufactured by BioVision, an Abcam company and was previously called K483 Human Whole Blood Polymorphonuclear Cell Isolation Kit. K483-30 is the same size as the 30 test size of ab241001.
--------------	--

### Properties

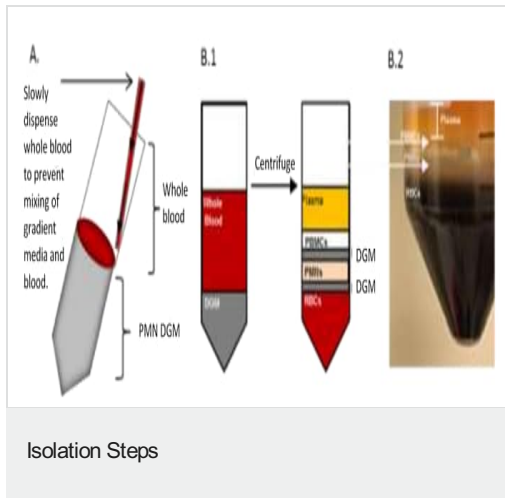
<b>Storage instructions</b>	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	30 tests
Blunt-End Needle	15 units
Giemsa Stain	1 x 7ml
PMN Density Gradient Solution	2 x 20ml
PMN Isolation Buffer	2 x 30ml
Viability Stain Mix	1 x 200µl

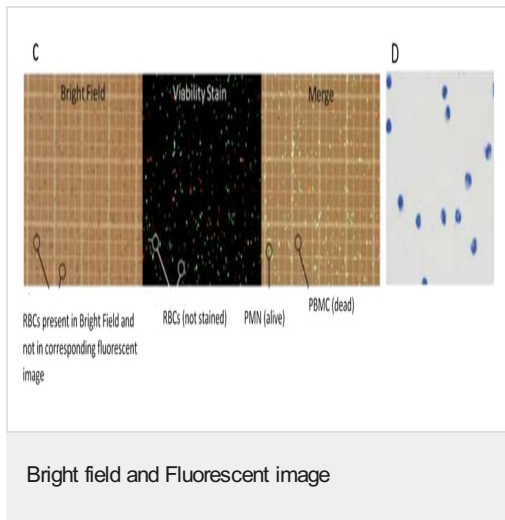
<b>Relevance</b>	Polymorphonuclear cells (PMN) have a multilobed nucleus and include the following cells types: mast cells, neutrophils, eosinophils and basophils. Upon isolation, these cells can be used for
------------------	--

migration assays and determination of cytokine/chemokine production in response to pharmacological compounds.

## Images



A. Illustration of conical tube held at 45 degrees while blood layered on top of PMN DGM. B.1. Layers of PMN DGM and whole blood prior to and after centrifugation showing the separation of layers in the conical tube. B.2 Illustrates separation of six layers (plasma, PBMCs, PMN DGM, PMNs, PMN DGM and RBCs) obtained after spinning whole blood with PMN DGM.



Bright-Field image of hemocytometer showing total cells(left); image from Fluorescent microscope with Rhodamine/FITC filters of same ROI showing live (green) and dead (red) cells (middle); merge of two panels (right). D. Polymorphonuclear Cells stained with Giemsa stain indicating purity of cell suspension obtained.

**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.com/abpromise> or contact our technical team.

## Terms and conditions

---

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