

Berthold Detection Systems GmbH  
Bleichstrasse 56–68  
D-75173 Pforzheim/Germany

Phone: +49(0)7231/9206-0  
Fax: +49(0)7231/9206-50  
E-Mail: [contact@berthold-ds.com](mailto:contact@berthold-ds.com)  
Internet: [www.berthold-ds.com](http://www.berthold-ds.com)

## Sirius L Tube Luminometer FB12/Sirius Software V2.0

### ABEL® Cell Activation Assay for Whole Blood or Isolated Cells with Pholasin® and Adjuvant-K™

The ABEL® (Analysis By Emitted Light) Cell Activation Assay is a chemiluminescent test that measures degranulation and the real-time production of free radicals during the activation of the NADPH oxidase system, the so-called 'respiratory burst'. The NADPH oxidase is most developed in granulocytic leucocytes, in particular neutrophils and monocytes, in which electrons are transported through the cell membrane to reduce oxygen to the free radical, superoxide anion  $O_2^-$ . Superoxide is released to the outside of the cell on initial activation and, during phagocytosis, into a phagosomal vacuole. The assay uses both the light-emitting protein Pholasin®, which reacts with free radicals to emit light, and an enhancer, Adjuvant-K™.

The assay can be used to:

- measure the respiratory burst in leucocytes
- study abnormalities in NADPH oxidase such as occur in different types of chronic granulomatous disease and various polymorphisms
- distinguish between receptor activation and intracellular activation of NADPH oxidase
- monitor changes in the activation of the NADPH oxidase in response to complement activation, infection, inflammation and medical intervention, thus having applications in:
  - biocompatibility monitoring
  - drug evaluation and testing
  - QC of vaccines
  - surgery
  - disease management
  - assessing neutrophil activity following chemotherapy, transplantation etc.
- monitor degranulation of myeloperoxidase
- measure free radicals and oxidants produced by a range of cell types including brain cells.

The test is simple, rapid and ultra-sensitive and capable of measuring fmols of superoxide from a small number of cells. The assay works on venous or capillary blood as well as on isolated cells and has been used on blood from humans, horses, cattle, dogs, cats, rats, mice, birds, fish and reptiles. The amount of blood required for a 1mL assay is 1µL but it is recommended to do an initial 1:100 dilution of 20µL.

The assay is available as a test kit that includes all the reagents, tubes to dilute the blood and two stimulants to initiate the respiratory burst: the receptor stimulant fMLP (f-Met-Leu-Phe), which activates NADPH oxidase via receptors on the surface, and the phorbol ester PMA that enters the cell and activates NADPH oxidase via the activation of protein kinase C. Other activators or receptor primers such as platelet activating factor or lipopolysaccharide can be used in place of, or together with, fMLP. Many non-human leucocytes do not have receptors for fMLP. In the presence of Pholasin® and Adjuvant-K™ the kinetics of this response can be determined by measurement of the light emitted during the few minutes of the analysis.

## Sirius L Tube Luminometer FB12/Sirius Software V2.0

### Materials:

Luminometer: Sirius L Tube Luminometer

Software: FB12/Sirius Software V2.0

Assay: ABEL® Cell Activation Test Kit with Pholasin® and Adjuvant-K™, Knight Scientific Ltd, Plymouth, UK (Product No. KSL-ABEL-04)

Cuvettes: Transparent polystyrene tubes (5 mL, 75 mm x 12 mm), Sarstedt, Leicester, UK (Product No. 555.476).

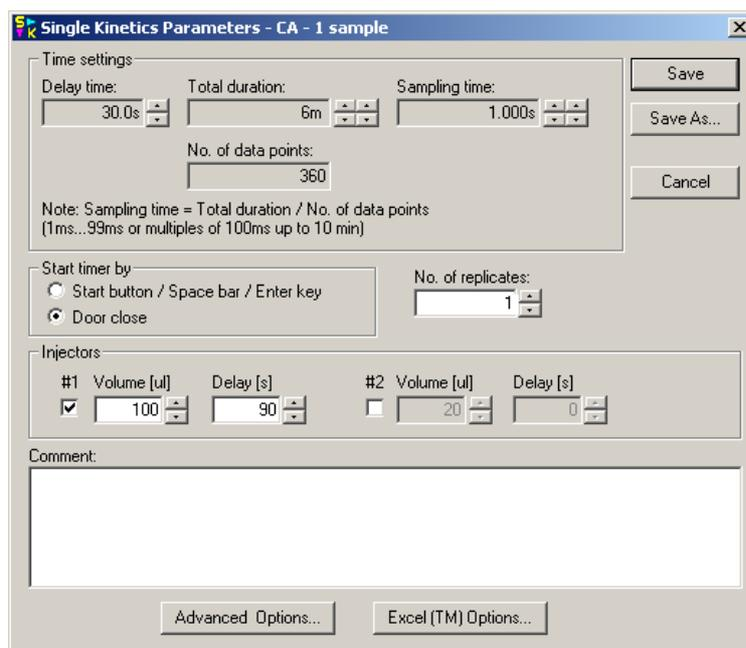
### Method:

For more detailed protocols please refer to the kit instructions supplied by the manufacturer. A product support document for the ABEL® Cell Activation assay can be downloaded from <http://www.knightscientific.com/downloads>.

1. Reconstitute reagents according to kit instructions.
2. Prime automatic reagent injectors (each injector has to be primed with at least 3 x 2.5mL of the respective solution):
  - a. Prime injector 1 with fMLP
  - b. Prime injector 2 with PMA
3. Dilute whole blood to 1:100 in 2 mL by adding 20µL of anti-coagulated blood (preferably EDTA) to 2mL of blood dilution buffer. For controls, plasma is used in place of blood.
4. To each cuvette add: 100 µL diluted blood (or diluted plasma control), 450 µL assay buffer, 250 µL Pholasin® (2.5 µg), 100 µL Adjuvant-K™ (1 enhancement unit).
5. Incubate the cuvette in a water bath at 37°C for 6 minutes.
6. Create a protocol in the FB12/Sirius software (Fig.1.).

Select Single Kinetics and set the following parameters:

- Delay = 30 seconds (allows background light to equilibrate)
- Sampling time = 1.000 seconds
- Inject 100 µL fMLP or PMA after 1 minute of reading
- Continue reading for 5 minutes (if using fMLP) or 7 minutes (if using PMA)



#1	Volume [ul]	Delay [s]	#2	Volume [ul]	Delay [s]
<input checked="" type="checkbox"/>	100	90	<input type="checkbox"/>	20	0

Figure 1. Parameter settings for assay using fMLP stimulant in Sirius Protocol Manager

## Sirius L Tube Luminometer FB12/Sirius Software V2.0

### Example:

Leucocyte activation was measured in whole blood from three healthy adult donors. Capillary blood samples were taken by finger prick using a safety lancet (Sarstedt) and stored in Microvette® tubes containing EDTA anticoagulant (Sarstedt). Blood plasma was used as a negative (no cell) control. The assay was performed as described above with samples assayed in duplicate to show the reproducibility of the assay.

### Results:

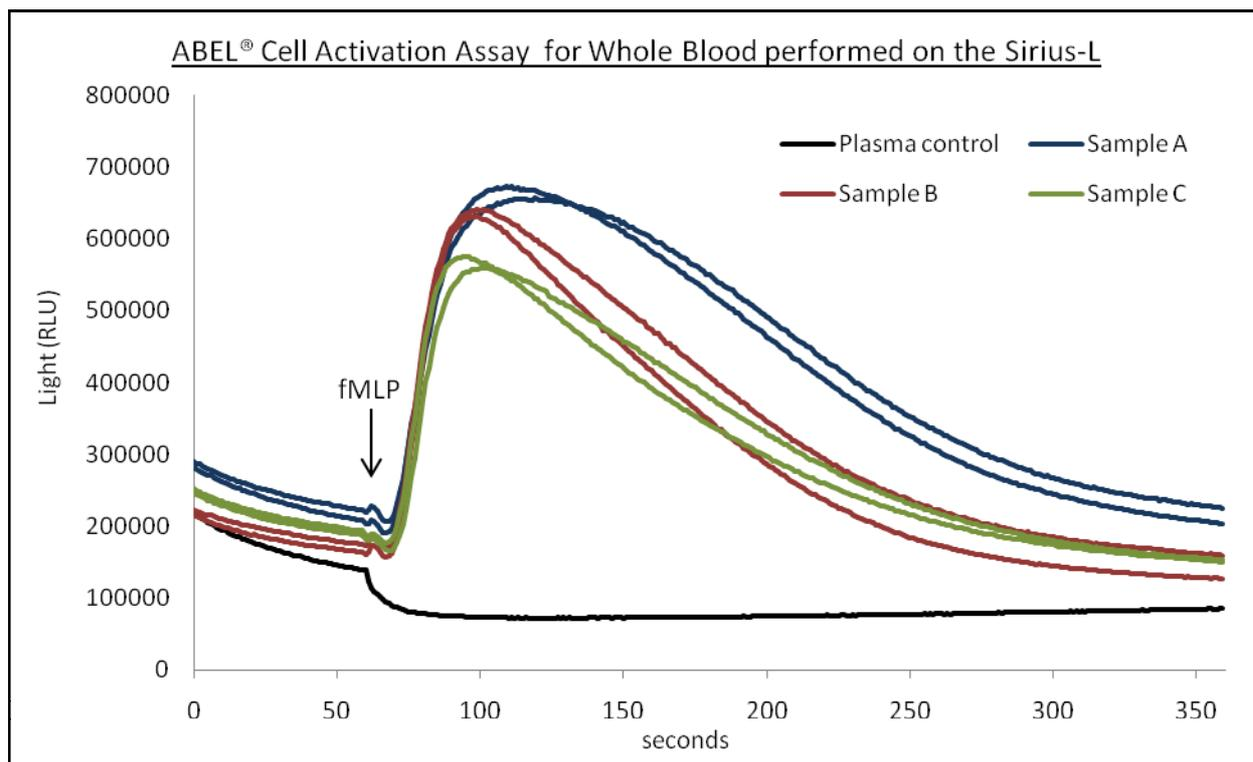


Figure 2: Free radical production by activated leukocytes in whole blood samples from 3 individuals monitored by the luminescence of Pholasin®. The cells are activated by the injection of the stimulant fMLP after 1 minute. The black line represents the plasma (no cell) control.

### Summary:

The activation of the NADPH oxidase in leucocytes in the blood can be monitored by the luminescence of Pholasin® enhanced with Adjuvant-K™.

The Sirius L tube luminometer showed an excellent ability to detect the production of reactive oxygen species produced by living cells from a small sample (20 µL) of capillary blood.

### Acknowledgement:

Data provided by Knight Scientific Ltd., UK  
Figures and text by courtesy of Knight Scientific Ltd., UK

