

# HEK-Blue™ IL-4/IL-13 Cells

Interleukin-4 and interleukin-13 reporter cells

Catalog code: hkb-il413

<https://www.invivogen.com/hek-blue-il4-il13>

For research use only

Version 23B14-MM

## PRODUCT INFORMATION

### Contents:

- 3-7 x 10<sup>6</sup> of HEK-Blue™ IL4/IL-13 cells in a cryovial or shipping flask.

**IMPORTANT:** If cells provided in a cryovial are not frozen upon arrival, contact InvivoGen immediately.

- 1 ml of **Blasticidin** (10 mg/ml). Store at 4 °C or at -20 °C.\*
- 1 ml of **Zeocin**® (100 mg/ml). Store at 4 °C or at -20 °C.\*
- 1 ml of **Normocin**™ (50 mg/ml), a formulation of three antibiotics active against mycoplasmas, bacteria and fungi. Store at -20 °C.\*

\*The expiry date is specified on the product label.

- 1 ml of **QB reagent** and 1 ml of **QB buffer** (sufficient to prepare 100 ml of **QUANTI-Blue™ Solution**, a SEAP detection reagent). QB reagent and QB buffer are stable for 1 year at -20 °C. QUANTI-Blue™ Solution is stable for 2 weeks at 4 °C and for 2 months at -20 °C.

### Handling Frozen Cells Upon Arrival

Cells must be thawed immediately upon receipt and grown according to handling procedures (as described on the next page) to ensure the best cell viability and proper assay performance.

**Note:** Avoid freezing cells upon receipt as it may result in irreversible damage to the cell line.

**Disclaimer:** We cannot guarantee cell viability if the cells are not thawed immediately upon receipt and grown according to handling procedures.

**IMPORTANT:** For cells that arrive in a shipping flask please refer to the enclosed 'cell recovery procedure'.

### Cell Line Stability

Cells will undergo genotypic changes over time that will result in reduced responsiveness in normal cell culture conditions. Genetic instability is a biological phenomenon that occurs in all stably transfected cells. Therefore, it is critical to prepare an adequate number of frozen stocks at early passages.

### Quality Control

- SEAP reporter activity in response to various cytokines has been validated using functional assays.
- The stability of this cell line for 20 passages following thawing has been verified.
- These cells are guaranteed mycoplasma-free.

## USE RESTRICTIONS

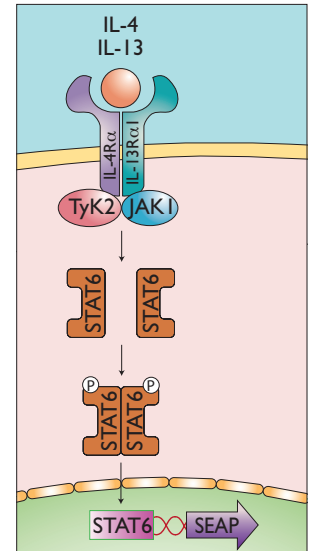
These cells are distributed for research purposes only.

This product is covered by a Limited Use License. By use of this product, the buyer agrees with the terms and conditions of all applicable Limited Use Label Licenses. For non-research use, such as screening, quality control or clinical development, contact [info@invivogen.com](mailto:info@invivogen.com).

## INTRODUCTION

The transcription factor STAT6 is activated primarily by two cytokines with overlapping biologic functions, IL-4 and IL-13. It can also be activated by IFN- $\alpha$  in a cell-specific manner.

In non-hematopoietic cells, IL-4 and IL-13 bind a receptor complex composed of the IL-4R $\alpha$  and IL-13R $\alpha$ 1. Upon ligand binding, the receptor complex activates the receptor-associated Janus kinases (JAK1 and Tyk2) leading to the recruitment of STAT6 and its phosphorylation. Activated STAT6 forms homodimers that translocate to the nucleus where they bind the promoter of responsive genes inducing gene transcription.



## CELL LINE DESCRIPTION

HEK-Blue™ IL-4/IL-13 cells are specifically designed to monitor the activation of the STAT6 pathway induced by IL-4 and IL-13. These cells were generated by stably introducing the human STAT6 gene into HEK293 cells to obtain a fully active STAT6 signaling pathway. The other genes of the pathway are naturally expressed in sufficient amounts.

Furthermore, HEK-Blue™ IL-4/IL-13 cells stably express the reporter gene secreted embryonic alkaline phosphatase (SEAP) under the control of the IFN- $\beta$  minimal promoter fused to four STAT6 binding sites.

Activation of the STAT6 pathway in HEK-Blue™ IL-4/IL-13 cells induces the expression of the reporter gene. SEAP which is secreted in the supernatant is easily detectable when using QUANTI-Blue™ Solution, a medium that turns purple/blue in the presence of SEAP.

HEK-Blue™ IL-4/IL-13 cells detect human IL-4 and human/murine IL-13

- Detection range for human IL-4: 0.1 - 10 ng/ml
- Detection range for human IL-13: 1 - 100 ng/ml

HEK-Blue™ IL-4/IL-13 cells are resistant to blasticidin and Zeocin®.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

## SAFETY CONSIDERATIONS

HEK-Blue™ IL-4/IL-13 cells were derived from HEK293 cells (transformed with adenovirus 5 DNA) that require **Biosafety level 2** according to the American Center for Disease Control and Prevention (CDC) guidelines. The biosafety level may vary depending on the country. For example, in Germany HEK293 cell lines are designated Biosafety Level 1 according to the Central Committee of Biological Safety, Zentrale Kommission für die Biologische Sicherheit (ZKBS). Please check with your country's regulatory authority regarding the use of these cells.

## HANDLING PROCEDURES

### Required Cell Culture Medium

- **Growth Medium:** DMEM, 4.5 g/l glucose, 2 mM L-glutamine, 10% (v/v) heat-inactivated fetal bovine serum (FBS; 30 min at 56 °), Pen-Strep (100 U/ml-100 µg/ml), 100 µg/ml **Normocin™**
- **Freezing Medium:** DMEM with 20% FBS and 10% (v/v) DMSO
- **Test Medium:** DMEM, 4.5 g/l glucose, 2 mM L-glutamine, 10% (v/v) heat-inactivated FBS, Pen-Strep (100 U/ml-100 µg/ml) **without Blasticidin, Zeocin® and Normocin™**

### Required Selective Antibiotics

- **Blasticidin and Zeocin®**

### Initial Culture Procedure

The first propagation of cells should be for generating stocks for future use. This ensures the stability and performance of the cells for subsequent experiments.

1. Thaw the vial by gentle agitation in a 37°C water bath. To reduce the possibility of contamination, keep the O-ring and cap out of the water. Thawing should be rapid.
2. Remove the vial from the water bath as soon as the contents are thawed, and decontaminate by dipping in or spraying with 70% (v/v) ethanol.
- Note: All steps from this point should be carried out under strict aseptic conditions.*
3. Transfer cells in a larger vial containing 15 ml of pre-warmed growth medium. **Do not add selective antibiotics until the cells have been passaged twice.**
4. Centrifuge vial at 150 x g (RCF) for 10 minutes.
5. Remove supernatant containing the cryoprotective agent and resuspend cells with 1 ml of growth medium without selective antibiotics.
6. Transfer the vial contents to a 25 cm<sup>2</sup> tissue culture flask containing 5 ml of growth medium without selective antibiotics.
7. Place the culture at 37°C in 5% CO<sub>2</sub>.

### Frozen Stock Preparation

1. Resuspend cells at a density of 5-7 x 10<sup>6</sup> cells/ml in freezing medium freshly prepared with cold growth medium.
- Note: A T-75 culture flask typically yields enough cells for preparing 3-4 frozen vials.*
2. Prepare 1 ml aliquots of cells in cryogenic vials.
3. Place vials in a freezing container and store at -80°C overnight.
4. Transfer vials to liquid nitrogen for long term storage.
- Note: If properly stored, cells should remain stable for years.*

## Cell Maintenance

1. HEK-Blue™ IL-4/IL-13 cells grow as adherent cells. Detach the cells using trypsin for 2-3 min at room temperature (RT).  
*Note: Prolonged action of trypsin or incubation at 37°C may alter the cell surface expression of receptors.*
2. Maintain and subculture the cells in growth medium supplemented with 10 µg/ml of **blasticidin** and 100 µg/ml of **Zeocin®**.
3. Renew growth medium twice a week.
4. Cells should be passaged when a 70-80% confluency is reached. Do not let the cells grow to 100% confluency.

## IL-4/IL-13 DETECTION

### Day 1:

1. Prepare HEK-Blue™ IL-4/IL-13 cell suspension: gently rinse cells twice with pre-warmed phosphate buffered saline (PBS), detach the cells in the presence of PBS for 2-3 min at 37°C. Tap the flask if needed. Resuspend cells in fresh, pre-warmed test medium and prepare a cell suspension at ~280,000 cells/ml.  
*Note: We recommend avoiding the use of trypsin to detach cells for the functional assays (see FAQs online).*
2. Add 20 µl of each sample per well of a flat-bottom 96-well plate.
3. Add 20 µl of **recombinant human IL-4** (1 ng/ml) or recombinant human IL-13 (10 ng/ml) as positive controls in one well.
4. Add 20 µl of TNF-α at 100 ng/ml (negative control, other cytokines can be used) in one well.
5. Add 180 µl of cell suspension (~50,000 cells) per well.
6. Incubate the plate at 37 °C in a CO<sub>2</sub> incubator for 20-24 h.

### Day 2:

1. Prepare **QUANTI-Blue™ Solution** following the instructions on the enclosed data sheet.
2. Add 180 µl of resuspended **QUANTI-Blue™ Solution** per well of a flat-bottom 96-well plate.
3. Add 20 µl of induced HEK-Blue™ IL-4/IL-13 cells supernatant.
4. Incubate the plate at 37 °C incubator for 1-3 h.
5. Determine SEAP levels using a spectrophotometer at 620-655 nm.

## RELATED PRODUCTS

Product	Description	Cat. Code
Blasticidin	Selection antibiotic	ant-bl-1
Normocin™	Antimicrobial reagent	ant-nr-1
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs
Recombinant human IL-4	Recombinant cytokine	rcyec-hil4
Zeocin®	Selection antibiotic	ant-zn-1

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

# QUANTI-Blue™ Solution

Medium for detection and quantification of alkaline phosphatase in standard and HTS assays

Catalog code: rep-qbs, rep-qbs2, rep-qbs3

<https://www.invivogen.com/quant-blue>

For research use only

Version 23A12-MM

## PRODUCT INFORMATION

**Contents:** QUANTI-Blue™ Solution is available in three pack sizes

- **rep-qbs:** 5 x 1 ml of QB reagent and 5 x 1 ml QB buffer, sufficient to prepare QUANTI-Blue™ Solution for **25 x 96-well plates** (500 ml using the standard procedure) or **20 x 1536-well plates** (85 ml using the HTS screening procedure).

- **rep-qbs2:** 10 x 1 ml of QB reagent and 10 x 1 ml QB buffer, sufficient to prepare QUANTI-Blue™ Solution for **50 x 96-well plates** (1 L using the standard procedure) or **40 x 1536-well plates** (170 ml using the HTS screening procedure).

- **rep-qbs3:** 1 x 20 ml bottle of QB reagent and 1 x 20 ml bottle of QB buffer, sufficient to prepare QUANTI-Blue™ Solution for **100 x 96-well plates** (2 L using the standard procedure) or **80 x 1536-well plates** (340 ml using the HTS screening procedure).

**Required Material (not provided)**

- Sterile water
- Sterile screw cap tube, glass bottle or flask

**Storage and stability**

- Product is shipped at room temperature. Upon receipt, store QB reagent and QB buffer at -20°C. Product is stable for 1 year at -20°C when properly stored.

- The 20 ml bottles of QB reagent and QB buffer are designed for single use. If required, individual aliquots of QB reagent and QB buffer can be prepared upon receipt or following a single freeze-thaw cycle. Store aliquots at -20°C. **Avoid repeated freeze-thaw cycles.**

*Note:* During storage, a precipitate may form in the 20 ml bottle of QB reagent. If this occurs, vortex the product until the precipitate disappears. The formation of a precipitate does not affect the activity of the product.

- Reconstituted QUANTI-Blue™ Solution is stable for 2 weeks at 2-8°C and for 2 months at -20°C. Protect QUANTI-Blue™ from light.

**Quality Control**

Each lot is thoroughly tested to ensure the absence of lot-to-lot variation.

- Physicochemical characterization (pH, appearance).
- Functional assays using alkaline phosphatase or SEAP-expressing reporter cells.

## DESCRIPTION

QUANTI-Blue™ is a colorimetric enzyme assay developed to determine any alkaline phosphatase activity (AP) in a biological sample, such as supernatants of cell cultures. QUANTI-Blue™ Solution changes from pink to a purple-blue color in the presence of AP. Secreted embryonic alkaline phosphatase (SEAP) is a widely used reporter gene. It is a truncated form of placental alkaline phosphatase, a glycosylphosphatidylinositol (GPI)-anchored protein. SEAP is secreted into the cell culture supernatant and therefore offers many advantages over intracellular reporters.

QUANTI-Blue™ is highly sensitive for quantitative measurement. It has a higher saturation threshold than with pNPP (p-nitrophenyl phosphate) resulting in more significant differences between no, low or high AP activity. Another advantage of QUANTI-Blue™ is that it can determine secreted AP activity without disturbing cells, thus allowing the repeated sampling of cell cultures for kinetic studies.

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

## METHODS

QUANTI-Blue™ Solution has been optimized for use in 96-well plates (standard procedure) and in 1536-well plates (high throughput screening procedure).

### A. Standard procedure

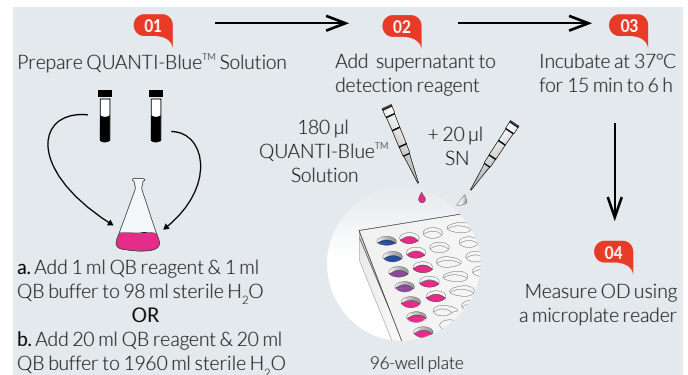


Figure 1. Standard procedure using QUANTI-Blue™ Solution.

The following protocol refers to the use of 96-well plates. Ensure QB reagent and QB buffer are completely thawed before use.

*Note:* For fast thawing, QB reagent and QB buffer can be placed at 37°C for 2 minutes. Ensure heating at 37°C does **not** exceed 5 minutes.

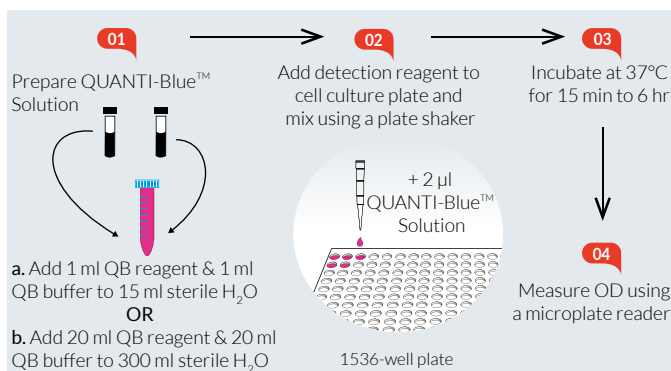
1. In a sterile bottle or flask, prepare QUANTI-Blue™ Solution by adding:
  - a. 1 ml of QB reagent and 1 ml of QB buffer to 98 ml of sterile water.
  - b. 20 ml of QB reagent and 20 ml of QB buffer to 1960 ml of sterile water.
2. Mix by vortexing and incubate at room temperature for 10 min before use.
3. Use QUANTI-Blue™ Solution immediately or store at 2-8°C or -20°C.
4. Dispense 180 µl of QUANTI-Blue™ Solution per well into a flat-bottom 96-well plate.
5. Add 20 µl of the sample (supernatant of SEAP-expressing cells) or negative control (cell culture medium).
6. Incubate at 37°C for 15 min to 6 h.
7. Measure optical density (OD) at 620-655 nm using a microplate reader.

*Note:* If the negative control turns purple/blue, it means the fetal bovine serum (FBS) contains alkaline phosphatase. We recommend heating FBS at 56°C for 30 min to inactivate the alkaline phosphatase activity.

For different cell culture plate formats, please refer to the table below:

	96-well plate	24-well plate	12-well plate
QUANTI-Blue™	180 µl	450 µl	900 µl
Supernatant	20 µl	50 µl	100 µl

## B. High Throughput Screening (HTS) procedure



**Figure 2. High throughput screening procedure using QUANTI-Blue™ Solution.**

This procedure has been optimized for use in HTS screening procedures in 1536-well plates. QUANTI-Blue™ Solution is added directly to the cell suspension to reduce liquid handling.

Ensure QB reagent and QB buffer are completely thawed before use.  
*Note:* For fast thawing, QB reagent and QB buffer can be placed at 37°C for 2 minutes. Ensure heating at 37°C does **not** exceed 5 minutes.

1. Dispense cell suspension and test compounds into a 1536-well plate in a volume that does not exceed **5 µl** per well. Incubate cells with test compounds for the desired period of time.
2. Prepare QUANTI-Blue™ Solution by adding:
  - a. **1 ml** of QB reagent and **1 ml** of QB buffer to **15 ml** of sterile water in a sterile 50 ml screw cap tube.
  - b. **20 ml** of QB reagent and **20 ml** of QB buffer to **300 ml** of sterile water in a sterile glass bottle or flask.
3. Mix well by vortexing and incubate at room temperature for 10 minutes before use.
4. Use QUANTI-Blue™ Solution immediately or store at 2-8°C or -20°C.
5. Dispense **2 µl** of QUANTI-Blue™ Solution to the wells containing  $\leq 5 \mu\text{l}$  of cell culture in a 1536-well plate.
6. Mix using a plate shaker.
7. Incubate at 37°C for 15 min to 6 h.
8. Measure OD at 620-655 nm.

*Note:* If the negative control turns purple/blue, it means the fetal bovine serum (FBS) contains alkaline phosphatase. We recommend heating FBS at 56°C for 30 min to inactivate the alkaline phosphatase activity.

## RELATED PRODUCTS

Product	Catalog Code
pNifTy2-SEAP (Zeo®)	pnifty2-seap
pSELECT-zeo-SEAP	psetz-seap
HEK-Blue™ Detection	hb-det2
Recombinant SEAP Protein	rec-hseap
<b>Reporter cells</b>	
HEK-Blue™ hTLR2	hkb-htlr2
HEK-Blue™ hTLR4	hkb-htlr4
RAW-Blue™ Cells	raw-sp
THP1-Blue™ NF-κB Cells	thp-nfkb
THP1-Blue™ ISG Cells	thp-isg

For a complete list of InvivoGen's Reporter Cell Lines visit <https://www.invivogen.com/reporter-cells>

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)