

HEK-Blue™ STAT6-hSTING-R232 Cells

Human STING (variant R232)-STAT6 reporter cells

Catalog code: hkb-st6r232

<https://www.invivogen.com/sting-r232-stat6-cells>

For research use only

Version 20K02-ED

PRODUCT INFORMATION

Contents:

- 3-7 x 10⁶ HEK-Blue™ STAT6-hSTING-R232 cells in a cryovial or shipping flask. IMPORTANT: If cells provided in a cryovial are not frozen upon arrival, contact InvivoGen immediately.
- 2 x 1 ml HEK-Blue™ Selection (250x concentrate) - a solution containing several selection antibiotics. Can be stored at 4°C or -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.

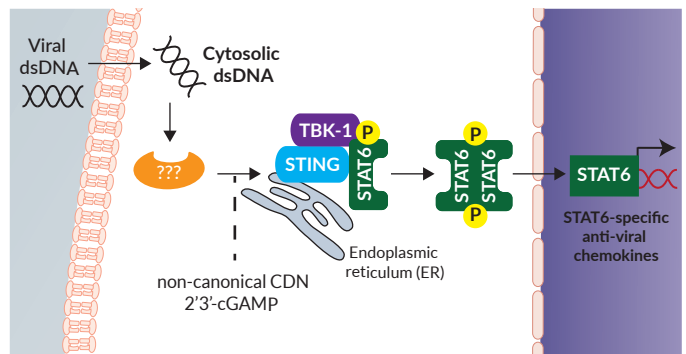
Quality Control

- STAT6-dependent SEAP reporter activity in response to various STING ligands and other cytokines has been validated.
- The stability of this cell line for 20 passages following thawing has been verified.
- These cells are guaranteed mycoplasma-free.

BACKGROUND

STING (stimulator of interferon genes) is essential for an effective response to cytoplasmic foreign or self-DNA. It directly senses cyclic dinucleotides (CDNs), which are important messengers in bacteria and innate immune agonists in mammals¹. The most prevalent variant of human STING, referred to as 'wild-type', is the R232 allele². The activation of STING causes a TANK-binding-kinase-1 (TBK1)-dependent cascade, ultimately, leading to IFN regulatory factor (IRF3)-dependent type I IFN production and NF-κB-dependent inflammatory cytokine production¹.

Additionally, signal transducer and activator of transcription 6 (STAT6) has been reported to be recruited to STING for TBK1-mediated phosphorylation during viral infection³. Ultimately, the activation of STAT6 induces a specific set of anti-viral chemokines responsible for immune cell homing, which is essential to reduce viral replication³. Notably, STING-dependent activation of STAT6 was found to be janus kinase (JAK)-independent and is thus, distinct from the 'canonical' STAT6 pathway activated by IL-4 and IL-13, which is crucial in adaptive immunity³.



1. Cheng, Z. *et al.* 2020. The interactions between cGAS-STING pathway and pathogens. *Signal Transduct Target Ther* 5, 91. 2. Yi, G. *et al.* 2013. Single nucleotide polymorphisms of human STING can affect innate immune response to cyclic dinucleotides. *PLoS One* 8, e77846. 3. Chen, H. *et al.* 2011. Activation of STAT6 by STING is critical for antiviral innate immunity. *Cell* 147, 436-446.

CELL LINE DESCRIPTION

HEK-Blue™ STAT6-hSTING-R232 cells are specifically designed to monitor the activation of the STAT6 signaling pathway induced by STING. These cells were generated by stable transfection and overexpression of the wild-type human (h)STING variant R232 in a human embryonic kidney (HEK)293-derived cell line that expresses human STAT6, and a STAT6-inducible secreted embryonic alkaline phosphatase (SEAP) reporter. The other genes of this signaling pathway are naturally expressed at sufficient levels. Numerous studies have indicated that type I interferons (IFNs) can activate STAT6, thus, HEK-Blue™ STAT6-hSTING-R232 cells have IFNα-R2 knocked out and subsequently, do not respond to type I IFNs, such as IFNα and IFNβ. HEK-Blue™ STAT6-hSTING-R232 cells stably express SEAP under the control of the IFN-β minimal promoter fused to four STAT6 binding sites. Thus, activation of the STAT6 pathway in these cells by hSTING agonists (i.e. 2'3'-cGAMP) induces the expression of SEAP, which is easily detectable in the cell culture supernatant with QUANTI-Blue™ Solution, a SEAP detection reagent. HEK-Blue™ STAT6-hSTING-R232 cells are resistant to Blasticidin, Hygromycin B, 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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com



Any questions about our cell lines?

Visit our FAQ page

 **InvivoGen**
www.invivogen.com

SAFETY CONSIDERATIONS

Biosafety Level 2

HEK-Blue™ STAT6-hSTING-R232 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 HEK-Blue™ Selection and Normocin™**

Required Selective Antibiotics

- **HEK-Blue™ Selection**

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 200-300 x g for 5 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 cm² tissue culture flask containing 5 ml of growth medium without selective antibiotics.
7. Place the culture at 37°C in 5% CO₂.

Frozen Stock Preparation

1. Resuspend cells at a density of 5-7 x 10⁶ 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

- Maintain and subculture the cells in growth medium supplemented with **HEK-Blue™ Selection**.
- Renew growth medium twice a week.
- Cells should be passaged when a 70-80% confluency is reached. Do not let the cells grow to 100% confluency.

REPORTER ASSAY

Day 1:

1. Add 20 µl of each sample per well of a flat-bottom 96-well plate.
2. 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.
3. Add 20 µl of TNF-α at 100 ng/ml (negative control, other cytokines can be used) in one well.
4. Prepare a cell suspension of HEK-Blue™ STAT6-hSTING-R232 cells (~3 x 10⁵ cells per ml) in test medium (containing 10% v/v heat-inactivated FBS).

Note: Some FBS may contain alkaline phosphatases that can interfere with SEAP quantification. We recommend to use heat-inactivated FBS to inactivate these enzymes which are thermosensitive.

5. Add 180 µl of cell suspension (~50,000 cells) per well.
6. Incubate the plate at 37 °C in a CO₂ 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™ STAT6-hSTING-R232 cell supernatant.
4. Incubate the plate at 37 °C for 1-3 h.
5. Determine SEAP levels using a spectrophotometer at 620-655 nm.

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.

RELATED PRODUCTS

| Product | Catalog Code |
|---------------------------|----------------|
| 2'3'-cGAMP | tlrl-nacga23 |
| 2'3'-cGAM(PS)2 (Rp/Sp) | tlrl-nacga2srs |
| 3'3'-cGAMP | tlrl-nacga |
| cAIM(PS)2 Difluor (Rp/Sp) | tlrl-nacairs |
| HEK-Blue™ Selection | hb-sel |
| Normocin™ | ant-nr-1 |
| QUANTI-Blue™ Solution | rep-qbs |
| Recombinant human IL-4 | rcyec-hil4 |
| THP1-Dual™ KI-hSTING-R232 | thpd-r232 |
| THP1-Dual™ KO-STING | thpd-kostg |

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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com



Any questions about our cell lines?

Visit our FAQ page

 **InvivoGen**
www.invivogen.com

QUANTI-Blue™ Solution

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

Catalog code: rep-qbs, rep-qbs2

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

For research use only

Version 19F11-MM

PRODUCT INFORMATION

Contents

QUANTI-Blue™ Solution is available in two pack sizes:

- **rep-qbs** containing 5 x 1 ml of QB reagent and 5 x 1 ml QB buffer to prepare 500 ml of QUANTI-Blue™ Solution sufficient for 25 x 96-well plates (standard procedure) or 20 x 1536-well plates (HTS screening)
- **rep-qbs2** containing 10 x 1 ml of QB reagent and 10 x 1 ml QB buffer to prepare 1 liter of QUANTI-Blue™ Solution sufficient for 50 x 96-well plates (standard procedure) or 40 x 1536-well plates (HTS screening)

Required Material (not provided)

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

Storage and Stability

- Store QB reagent and QB buffer at -20°C. Product is stable for 1 year at -20°C when properly stored.
- 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 (including pH, solubility).
- 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 GPI-anchored protein. SEAP is secreted into cell culture supernatant and therefore offers many advantages over intracellular reporters.

FEATURES AND ADVANTAGES

- **Requires small samples of cell supernatants** - 20 µl is sufficient.
- **No need to process samples** - Preparation of cell lysates or heating of samples is not required.
- **Determine secreted AP activity without disturbing cells** - The same cell cultures can be repeatedly sampled for kinetic studies.
- **Assay can be completed in 30 min** - Hands-on time no longer than 10 min. The enzymatic activity can be detected as early as 15 min after incubation of the samples in QUANTI-Blue™ Solution.
- **Wide dynamic range allows to detect low and high levels of AP** - No need to perform multiple sample dilutions.
- **Highly sensitive for quantitative measurement** - Higher saturation threshold than with pNPP (p-nitrophenyl phosphate) resulting in more significant differences between no, low or high AP activity.
- **Extremely simple to use** - 1) Prepare solution with water, 2) add sample to detection reagent, 3) incubate at 37°C, and 4) assess AP activity.

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

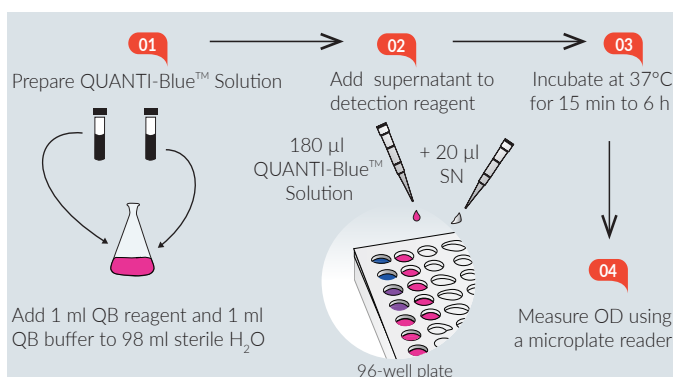


Figure 1. Standard procedure using QUANTI-Blue™ Solution.

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. Prepare 100 ml of QUANTI-Blue™ Solution by adding 1 ml of QB reagent and 1 ml of QB buffer to 98 ml of sterile water in a sterile glass bottle or flask.
 2. Mix well 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 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 to heat 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 |

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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

B. High Throughput Screening (HTS) procedure

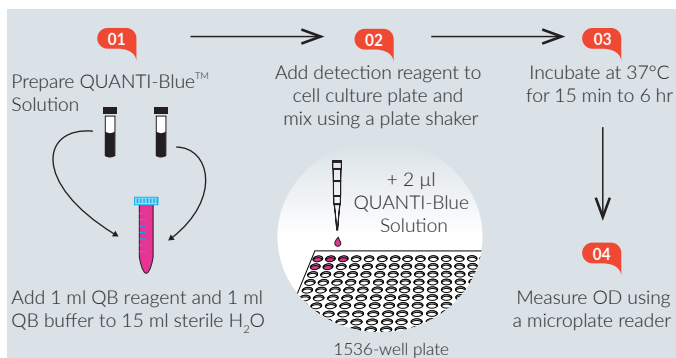


Figure 2. High throughput screening procedure using QUANTI-Blue™ Solution.

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 17 ml of QUANTI-Blue™ Solution by adding 1 ml of QB reagent and 1 ml of QB buffer to 15 ml of sterile water in a 50 ml screw cap tube.
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 ≤ 5 µl 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 to heat 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 |
| Reporter cells | |
| 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 Hong Kong: +852 3622-3480
E-mail: info@invivogen.com