

A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells

MDA5 knockout NF-κB-SEAP & IRF-Lucia reporter lung carcinoma cells permissive to SARS-CoV-2

Catalog code: a549d-komda5-cov2r

<https://www.invivogen.com/a549dual-hace2tmprss2-cells>

For research use only

Version 21C19-NJ

PRODUCT INFORMATION

Contents

• 3-7 x 10⁶ A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells in a cryovial or shipping flask.

- 1 ml of **Blasticidin** (10 mg/ml). Store at 4°C or at -20°C.*
- 1 ml of **Hygromycin B Gold** (100 mg/ml). Store at 4°C or at -20°C.*
- 1 ml of **Puromycin** (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 & 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.

• 1 pouch of **QUANTI-Luc™**, a Lucia luciferase detection reagent. Store pouch at -20°C for 12 months. Reconstituted QUANTI-Luc™ is stable for 1 week at 4°C and for 1 month at -20°C. Protect from light. *Note: Data sheets for all components are available on our website.*

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

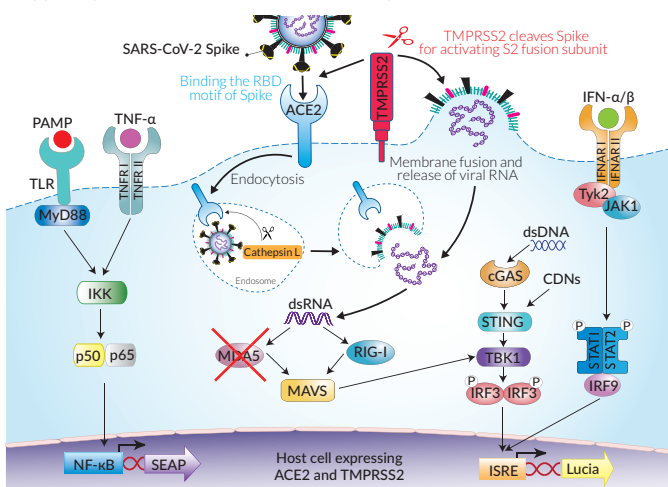
Cells undergo genotypic changes resulting in reduced responsiveness over time in normal cell culture conditions. Genetic instability is a biological phenomenon that occurs in all stably transfected cells. Therefore, it is critical to prepare an adequate number of frozen stocks at early passages. A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells should not be passaged more than 20 times to remain fully efficient.

Quality Control

- The overexpression of the human ACE2 (hACE2) gene has been verified by RT-qPCR, FACS staining, and functional assays.
- The overexpression of the human TMPRSS2 (hTMPRSS2) gene has been verified by RT-qPCR and functional assays.
- MDA5 KO has been verified by (RT-q)PCR and functional assays.
- Lucia luciferase and SEAP reporter activities have been validated using functional assays.
- The stability for 20 passages following thawing has been verified
- These cells are guaranteed mycoplasma-free.

BACKGROUND

MDA5 (melanoma-differentiation-associated gene 5) is a cytoplasmic PRR that plays an important role in antiviral response¹. It senses long double-stranded RNA (dsRNA), a replication intermediate for RNA viruses, leading to the production of type I interferons in infected cells¹. SARS-CoV-2, the causative agent of COVID-19, gains entry into target cells through ACE2 (angiotensin I-converting enzyme-2) and TMPRSS2 (transmembrane protease serine 2) surface receptors². MDA5 has been described as a major sensor of SARS-CoV-2 infection, triggering the interferon response in lung cells³.



CELL LINE DESCRIPTION

A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells were generated from A549-Dual™ KO-MDA5 lung carcinoma cells through stable integration of the human ACE2 and TMPRSS2 genes. A549-Dual™ KO-MDA5 cells feature a KO of the MDA5 gene, as well as stable integration of two inducible reporter constructs. The hACE2 and hTMPRSS2 expression allows infection by SARS-CoV-2. The secreted embryonic alkaline phosphatase (SEAP) and Lucia luciferase reporter genes allow the study of NF-κB and interferon regulatory factor (IRF) signaling pathways, respectively. Both reporter proteins are readily measurable in the culture supernatant when using QUANTI-Blue™, a SEAP detection reagent, and QUANTI-Luc™, a Lucia luciferase detection reagent. A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells are resistant to **Blasticidin**, **Hygromycin**, **Puromycin**, and **Zeocin™**.

APPLICATIONS

These cells are ideal for studying SARS-CoV-2 viral entry into host cells, the viral RNA sensing pathway, as well as screening small molecule inhibitors and neutralizing antibodies. These cells are permissive to infection by SARS-CoV-2 Spike-pseudotyped viral particles.

1. Reikine S, et al., 2014. Pattern Recognition and Signaling Mechanisms of RIG-I and MDA5. Front Immunol. 5:342. 2. Hoffmann M, et al., 2020. SARS-CoV-2 cell entry depends on ACE2 and TMPRSS2 and is blocked by a clinically proven protease inhibitor. Cell. 181:1-16. 3. Yin X, et al., 2021. MDA5 governs the innate immune response to SARS-CoV-2 in lung epithelial cells. Cell Reports. 34:108628.

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 1

HANDLING PROCEDURES

Required Cell Culture Medium

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

Note: Heat-inactivated FBS is also commercially available.

Required Selective Antibiotics

- **Blasticidin, Hygromycin, Puromycin, 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 T-25 culture flask containing 5 ml of growth medium.

Note: To avoid excessive alkalinity of the medium during recovery of the cells, place the tissue culture flask containing the growth medium into an incubator for at least 15 minutes before adding the vial contents.

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 1-2 frozen vials.
2. Dispense 1 ml of cell suspension into 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. A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells grow as adherent cells. To detach cells, rinse the cell layer with PBS, then incubate with 0.25% trypsin-EDTA for 2-5 minutes. Do not use a cell scraper.
2. Maintain and subculture the cells in growth medium supplemented with 10 µg/ml of **Blasticidin**, 100 µg/ml of **Hygromycin**, 0.5 µg/ml of **Puromycin**, 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.

Note: The average doubling time for the A549-Dual™ KO-MDA5-hACE2-TMPRSS2 cells is ~48 hours using the conditions described above.

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

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 to 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

REPORTER ASSAYS

Use A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells with A549-Dual™ hACE2-TMPRSS2 cells. *Note:* For best results, cells should be at 50-70% confluency on the day of the experiment.

Cell preparation

1. Rinse cell layer with PBS and detach cells with trypsin.
2. Centrifuge cells at 300 x g (RCF) for 5 min.
3. Remove supernatant and resuspend A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells at 2.8 x 10⁵ cells/ml in fresh, pre-warmed growth medium.

NF-κB induction

1. Add 20 µl of sample per well of a flat-bottom 96-well plate, including a positive control (e.g. recombinant human TNF-α) and endotoxin free water as a negative control

Note: Use new tips for each well to avoid cross-contamination.

2. Add 180 µl of cell suspension (~50,000 cells) per well.
3. Incubate the plate at 37°C in a CO₂ incubator for 18-24 h.
4. Prepare **QUANTI-Blue™ Solution** following the instructions on the enclosed data sheet.
5. Add 20 µl of A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells supernatant.
6. Add 180 µl of resuspended **QUANTI-Blue™ Solution** per well of a flat-bottom 96-well plate.
7. Incubate the plate at 37°C in a CO₂ incubator for 1-8 h.
8. Determine NF-κB-induced SEAP levels using a microplate spectrophotometer at 620-655 nm.

IRF induction

Below is a protocol for end-point readings using a luminometer, this protocol can be adapted for use with kinetic measurements.

1. Add 20 µl of sample per well including a positive control (e.g. recombinant IFN-α or IFN-β) and endotoxin free water as a negative control.

Note: Use new tips for each well to avoid cross-contamination.

2. Add 180 µl of cell suspension (~50,000 cells) per well of a flat-bottom 96-well plate.
3. Incubate the plate at 37°C in a CO₂ incubator for 18-24 h.
4. Prepare the **QUANTI-Luc™** assay solution following the instructions on the enclosed data sheet.
5. Pipet samples (20 µl per well) into a 96-well white (opaque) or black plate, or a luminometer tube.
6. Add 50 µl **QUANTI-Luc™** per well.
7. Proceed **immediately** with the measurement.

RELATED PRODUCTS

Product	Catalog Code
A549-Dual™ hACE2-TMPRSS2 Cells	a549d-cov2r
Blasticidin	ant-bl-1
Hygromycin B Gold	ant-hg-1
Puromycin	ant-pr-1
Zeocin™	ant-zn-1
3p-hpRNA/LyoVec™	tlrl-hprnalv
QUANTI-Blue™ Solution	rep-qbs
QUANTI-Luc™	rep-qlc1



Any questions about our cell lines?
Visit our FAQ page.

 **InvivoGen**
www.invivogen.com

QUANTI-Luc™

A coelenterazine-based luminescence assay reagent

Catalog code: rep-qlc1, rep-qlc2

<https://www.invivogen.com/quantiluc>

For research use only

Version 19A04-MM

PRODUCT INFORMATION

Contents

QUANTI-Luc™ is provided as packs of individually sealed pouches.

- rep-qlc1: 2 pouches of QUANTI-Luc™
- rep-qlc2: 5 pouches of QUANTI-Luc™

Each pouch contains everything needed to prepare 25 ml of reagent allowing the preparation of 500 wells of a 96-well plate.

Storage and Stability

- Store QUANTI-Luc™ pouches at -20°C for 12 months.
- Reconstituted QUANTI-Luc™ is stable for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

Note: This product is photosensitive and should be protected from light.

DESCRIPTION

QUANTI-Luc™ is an assay reagent containing all the components required to quantitatively measure the activity of Lucia luciferase and other coelenterazine-utilizing luciferases. QUANTI-Luc™ contains the coelenterazine substrate and stabilizing agents for the luciferase reaction. The light signal produced is quantified using a luminometer and expressed as relative light units (RLU). The signal produced correlates to the amount of luciferase protein expressed, indicating promoter activity in the reporter assay.

QUANTI-Luc™ is optimized for use with Lucia luciferase reporter cell lines. Lucia luciferase is a secreted coelenterazine luciferase encoded by a synthetic gene. As Lucia luciferase is secreted, it can be directly measured in the cell culture medium using bioluminescent assays.

InvivoGen provides a recombinant Lucia luciferase protein (see Related Products) which is a positive control for QUANTI-Luc™. A dilution series of the recombinant Lucia luciferase protein can also be used to determine the linear range of the assay.

METHODS

Preparation of QUANTI-Luc™

1. Pour the pouch contents into a 50 ml screw cap tube.
2. Add 25 ml of sterile water.
3. Swirl product gently until powder is completely dissolved.
4. Use QUANTI-Luc™ assay solution immediately or store until required for use. Reconstituted QUANTI-Luc™ can be stored for 1 week at 4°C and for 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles.

Note: This product is photosensitive and should be protected from light.

Detection of luciferase activity from cell culture medium

To obtain **end-point readings** using a luminometer **with an injector**.

1. Set the luminometer with the following parameters: 50 µl of injection, end-point measurement with a 4 second start time and 0.1 second reading time.
2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Prime the injector with the QUANTI-Luc™ assay solution and proceed **immediately** with the measurement.

To obtain **end-point readings** using a luminometer **without injectors**.

1. Set the luminometer with a 0.1 second reading time.
2. Pipet 10-20 µl of sample per well into a 96-well white (opaque) or black plate, or a luminometer tube.
3. Add 50 µl of QUANTI-Luc™ assay solution to each well or tube.
4. Gently tap the plate several times to mix (do **not** vortex).
5. Proceed **immediately** with the measurement.

RELATED PRODUCTS

Product	Catalog Code
QUANTI-Luc™ Gold (For standard and HTS assays)	rep-qlcg1
pSelect-zeo-Lucia™ (expression plasmid)	psetz-lucia
Recombinant Lucia luciferase protein	rec-lucia
Reporter Cells	
THP1-Dual™ (IRF-Lucia/NF-κB-SEAP) Cells	thpd-nfis
THP1-Lucia™ NF-κB Cells	thp1-nfkb

For a complete list of InvivoGen's Lucia luciferase Reporter Cell Lines visit <https://www.invivogen.com/lucia-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



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 20C16-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 (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 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 Hong Kong: +852 3622-3480

E-mail: 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

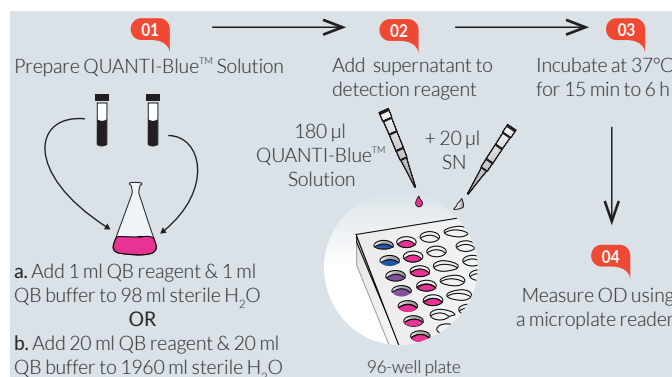


Figure 1. Standard procedure using QUANTI-Blue™ Solution.

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

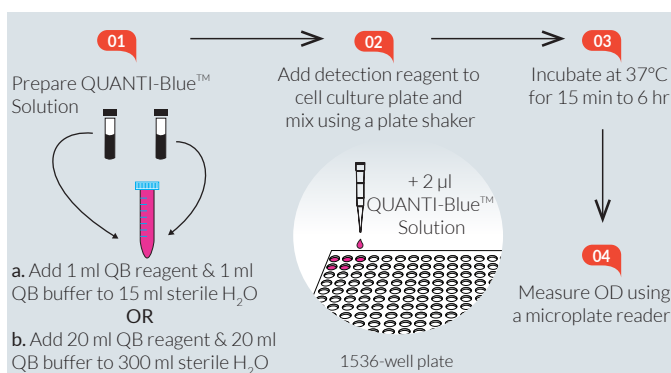


Figure 2. High throughput screening procedure using QUANTI-Blue™ Solution.

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
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