

HEK-Blue™ mDectin-2 Cells

SEAP Reporter 293 cells expressing the mouse Dectin-2 gene

Catalog code: hkb-mdect2

<https://www.invivogen.com/hek-blue-mdectin2>

For research use only

Version 20C24-MM

PRODUCT INFORMATION

Contents and Storage

• **3-7 x 10⁶ HEK-Blue™ mDectin-2 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 puromycin** (10 mg/ml). Store at 4 °C or at -20 °C.*
• **2 x 1 ml of HEK-Blue™ CLR Selection** (250x concentrate); a solution containing several selection antibiotics. Store HEK-Blue™ CLR Selection 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 pouch of HEK-Blue™ Detection**, a cell culture medium (50 ml) for real-time detection of SEAP. Store pouch at 4 °C for 6 months. Reconstituted HEK-Blue™ Detection is stable for 2 weeks at 4 °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 will 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. HEK-Blue™ mDectin-2 cells should not be passaged more than 20 times to remain fully efficient.

Quality Control

- SEAP reporter activity in response to Dectin-2 agonists and various other pathogen recognition receptor (PRR) agonists has been validated using functional assays.
- The cell surface expression of murine Dectin-2 (mDectin-2) in this cell line has been validated using fluorescence-activated cell sorting (FACS).
- The stability 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 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.

BACKGROUND

Dectin-2 is a member of the C-type lectin receptor (CLR) family and plays an important role in antifungal innate immunity^{1,2}. CLRs comprise a large family of receptors that bind to carbohydrates in a calcium-dependent manner. The lectin activity of these receptors is mediated by conserved carbohydrate-recognition domains (CRDs). Dectin-2 is a type II transmembrane CLR with a single CRD domain. It binds high mannose-type carbohydrates and has been demonstrated to be the functional receptor for α -mannans. Upon binding to its ligand, Dectin-2 signals through the kinase Syk and the adaptors CARD9/Bcl-10/MALT1 triggering the activation of NF- κ B and the subsequent production of pro-inflammatory cytokines^{3,4}. Notably, it has been demonstrated that Dectin-2 is important for the induction of Th17 in response to fungal infection⁴.

1. Hou H. et al., 2017. C-type Lectin Receptor: Old Friend and New Player. *Med Chem.* 2017;13(6):536-543. **2. Wang HL. et al., 2017.** Ligation of Dectin-2 with a novel microbial ligand promotes adjuvant activity for vaccination. *PLoS Pathog.* 213(8):e1006568. **3. Feinberg R. et al., 2017.** Mechanism of pathogen recognition by human dectin-2. *J Biol Chem.* 292(32):13402-13414. **4. Tang J. et al., 2018.** Regulation of C-Type Lectin Receptor-Mediated Antifungal Immunity. *Front Immunol.* 9:123.

PRODUCT DESCRIPTION

HEK-Blue™ mDectin-2 cells stably express the mouse Dectin-2 gene, as well as the genes of the Dectin-2 signaling pathway. They also stably express an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene. These reporter cell lines are activated by Dectin-2 ligands. They do not respond to other CLR ligands, such as trehalose- α , β -dibehenate (TDB), a Mincle ligand.

Levels of SEAP can be easily determined with HEK-Blue™ Detection, a cell culture medium that allows for real-time detection of SEAP activity. HEK-Blue™ Detection is extremely simple to use. It is applicable to high-throughput screening. HEK-Blue™ Detection contains all the nutrients necessary for cell growth and a specific SEAP color substrate. The hydrolysis of the substrate by SEAP produces a purple/blue color that can be easily detected with the naked eye or measured with a spectrophotometer.

SEAP activity can also be assessed using the alkaline phosphatase detection reagent, QUANTI-Blue™. With the QUANTI-Blue™ assay, cells are stimulated in a culture medium containing heat-inactivated fetal bovine serum. Following cell activation, QUANTI-Blue™ is used to detect SEAP in the cell supernatant. This colorimetric assay allows the same cell cultures to be repeatedly sampled for kinetic studies or further experimentation.

For more information, visit <https://www.invivogen.com/quanti-blue>.

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 3622-34-80

E-mail: info@invivogen.com



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



SAFETY CONSIDERATIONS

Biosafety Level: 2

HEK-Blue™ mDectin-2 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 Media

- **Growth Medium:** DMEM (4.5 g/l glucose), 10% (v/v) fetal bovine serum (FBS), Pen-Strep (100 U/ml-100 µg/ml), 100 µg/ml Normocin™, 2 mM L-glutamine

- **Freezing Medium:** DMEM, 20% (v/v) FBS and 10% (v/v) DMSO

Required Selective Antibiotics

- **Blasticidin, puromycin and HEK-Blue™ CLR 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 must 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% ethanol. *Note: All of the operations 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 1000-1200 RPM (RCF = 200-300 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 T-25 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 3-7 x 10⁶ cells/ml in freezing medium prepared extemporaneously with cold DMEM.

Note: A T-75 culture flask typically yields enough cells for preparing 3-4 frozen vials.

2. Aliquot 1 ml cells 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. Maintain and subculture the cells in growth medium supplemented with 30 µg/ml **blasticidin**, 1 µg/ml **puromycin** and 1X HEK-Blue™ CLR Selection.
2. Renew growth medium 2 times a week.
3. Cells should be passaged when a 70-80% confluency is reached. Detach the cells in presence of PBS by tapping the flask or by using a cell scraper. Do not let the cell grow to 100% confluency.

Note: The response of HEK-Blue™ mDectin-2 cells can be altered by the use of trypsin. Avoid trypsin to detach HEK-Blue™ mDectin-2 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

REPORTER ASSAY

HEK-Blue™ Detection is a cell culture medium that allows the detection of SEAP as the reporter protein is secreted by the cells. Prepare HEK-Blue™ Detection following the instructions on the enclosed data sheet.

Note: Before the test, the cells should be 50-80% confluent.

1. Coat several wells of a 96-well plate with a Dectin-2 ligand, such as Furfurman (10-1 µg/ml), a cell wall preparation of *Malassezia furfur*. These wells serve as positive controls for the reporter assay.

- 1.a - To coat wells, resuspend the Dectin-2 ligand in water, dispense a small volume (20 µl) of the Dectin-2 ligand per well of a 96-well plate and allow to dry for 1 hour at 70 °C.

- 1.b - Alternatively, resuspend the Dectin-2 ligand in isopropanol, dispense a small volume (20 µl) per well of a 96-well plate and allow to dry for 1 hour at room temperature.

2. Add 20 µl of each test compound per well of a 96-well plate. *Note: To coat the wells with the Dectin-2 ligand present in your samples, one of the methods described above may be used.*

3. Add 20 µl of a negative control (e.g. sterile endotoxin-free water) in one well.

4. Remove HEK-Blue™-mDectin-2 cells from the incubator and discard growth medium.

5. Gently rinse cells with pre-warmed 5-10 ml PBS (for a T-75 flask).

6. Add 2-5 ml pre-warmed PBS (for a T-75 flask) and place the cells at 37 °C for 1-2 min, detach the cells by tapping the flask or by using a cell scraper. Dissociate cell clumps by gently pipetting up and down.

Note: Do not use trypsin to detach HEK-Blue™-mDectin-2 cells.

7. Count cells which have been resuspended in pre-warmed PBS.

Note: Avoid centrifugation of HEK-Blue™-mDectin-2 cells.

8. Prepare a cell suspension ~280,000 cells per µl in HEK-Blue™ Detection medium and immediately add 180 µl of the cell suspension (~50,000 cells) per well.

Note: Avoid prolonged incubation of cells at room temperature in HEK-Blue™ Detection medium as it may lead to high background or false positive readings.

9. Incubate the plate at 37°C in 5% CO₂ for 16-24h. Hydrolysis of SEAP substrate can be observed with naked eye and determined using a spectrophotometer at 620-655 nm.

Specificity of HEK-Blue™ mDectin-2 cells

As their parental cell line, HEK293 cells, express endogenous levels of TLR3, TLR5 and NOD1, HEK-Blue™ mDectin-2 cells will respond to TLR3, TLR5 and NOD1 agonists, such as **poly(I:C)**, **flagellin** and **C12-iE-DAP**, respectively. In order to identify Dectin-2-specific responses, we recommend to use **HEK-Blue™ Null1-v cells** as a control cell line.

RELATED PRODUCTS

Product	Description	Cat. Code
Blasticidin	Selection antibiotic	anti-bl-1
Furfurman	Dectin-2 agonist	t1rl-ffm
HEK-Blue™ CLR Selection	Selection antibiotics	hb-csm
HEK-Blue™ Detection	SEAP detection medium	hb-det2
HEK-Blue™ Null1-v cells	Parental cell line	hkb-null1v
Puromycin	Selection antibiotic	ant-pr-1
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs



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



HEK-Blue™ Detection

Cell culture medium for the real-time detection of secreted alkaline phosphatase

Catalog code: hb-det2, hb-det3

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

For research use only

Version 23L22-MM

PRODUCT INFORMATION

Contents

HEK-Blue™ Detection is provided in sealed pouches and is available in two quantities:

- hb-det2: 5 pouches
- hb-det3: 10 pouches

Each pouch contains everything needed to prepare 50 ml of medium for the colorimetric detection of secreted embryonic alkaline phosphatase (SEAP).

Storage and stability

- Store sealed pouches at 2-8°C. Unopened pouches are stable for at least 6 months when stored properly.

Important: For the exact expiry date please see the corresponding CoA.

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

DESCRIPTION

HEK-Blue™ Detection is a cell culture medium developed to provide a fast and convenient method to monitor SEAP expression. Detection of SEAP occurs as the reporter protein is secreted by the cells grown in HEK-Blue™ Detection, which will change to a purple/blue color in the presence of alkaline phosphatase activity.

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. It allows the determination of reporter activity without disturbing the cells, does not require the preparation of cell lysates, and can be used for kinetic studies. Using HEK-Blue™ Detection, SEAP expression can be observed visually, and unlike fluorescent or luminescent reporters can be easily quantified using a microplate reader or spectrophotometer.

HEK-Blue™ Detection is applicable for high-throughput screening.

METHODS

Preparation of HEK-Blue™ Detection

1. Pour the contents of one pouch of HEK-Blue™ Detection into a sterile vial/bottle.
2. Solubilize the powder with 50 ml of cell culture grade water.
3. Vortex vigorously until powder is completely dissolved.
4. Warm reconstituted HEK-Blue™ Detection to 37°C for at least 3 hours.
5. Filter the medium through a 0.2 µm bottle-top vacuum filter into a sterile vial/bottle.
Note: We recommend using filter units providing a large filter area to facilitate filtration.
6. Keep the HEK-Blue™ Detection medium at 37°C before use or store at 2-8°C for up to 2 weeks.

Detection of SEAP activity

The following protocol is for the use of HEK-Blue™ Detection in 96-well plates. This will vary slightly depending on the volume of reagents needed, based on different plate sizes.

1. Prepare the cell suspension by detaching the cells and resuspending in a small volume of PBS.
2. Count the cells.
3. Add an appropriate amount of PBS-resuspended cells in HEK-Blue™ Detection to obtain a cell suspension at the expected concentration.
4. Add 20 µl of SEAP-inducer compound or negative control (such as PBS) per well.
5. Add 180 µl of cell suspension per well.
Note: To obtain more consistent results, we recommend to mix the SEAP-inducer and cell suspension by pipetting up and down.
6. Incubate overnight at 37°C, in 5% CO₂.
7. Determine SEAP activity with the naked eye or by reading the optical density (OD) at 620-655 nm.

RELATED PRODUCTS

Product	Description	Cat. Code
pSELECT-zeo-SEAP	SEAP reporter gene	psetz-seap
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs
Recombinant SEAP Protein	Control for SEAP assays	rec-hseap

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

