

# 293XL-hTLR8A

293XL cell line expressing the human TLR8a gene

Catalog # 293xl-htr8

For research use only

Version # 10D14-MM

## PRODUCT INFORMATION

### Contents and Storage

- 1 vial of 293XL-hTLR8A Cells (5-7 x 10<sup>6</sup> cells) in Freezing Medium
- IMPORTANT:** Cells are shipped frozen. If cells are not frozen upon arrival, contact InvivoGen immediately.
- 100 µl Blastidicin selective antibiotic (10 mg/ml). Store at -20°C. Product is stable for 1 year when stored at -20°C.
  - 1 ml Normocin™ (50 mg/ml). Normocin™ is a formulation of three antibiotics active against mycoplasmas, bacteria and fungi. Store at -20°C. Product is stable for 18 months when stored at -20°C.

## PRODUCT DESCRIPTION

293XL-hTLR8A cells are designed for studying the stimulation of human TLR8a (hTLR8a). 293XL-hTLR8A cells were obtained by co-transfection of the hTLR8 and the human antiapoptotic Bcl-XL genes. HEK293 cells express endogenous levels of TLR3, TLR5, and NOD1. *Note: The control cell line for 293XL-hTLR8A cells is 293XL/null cells (which do not express hTLR8).*

TLR8 was identified together with TLR7 and TLR9 and is expressed more abundantly in PBL and lung<sup>1</sup>. Human TLR8 and TLR7 were reported to independently confer responsiveness to R848, an imidazoquinolone with antiviral activity. R848 induces the activation of NF-κB in HEK293 cells transfected with human TLR8 but not in HEK293 cells expressing murine TLR8, suggesting a species specificity<sup>2</sup>. Sequence-specific single-stranded RNA (ssRNA) was identified as the natural ligand of human TLR8<sup>3</sup>. GU-rich ssRNA oligonucleotides derived from HIV-1 were shown to activate DC and macrophages. TLR8 recognizes ssRNA in the endosomal compartment and leads to the secretion of cytokines<sup>3</sup>.

1. **Chuang TH, and RJ. Ulevitch, 2000.** Cloning and characterization of a sub-family of human toll-like receptors: hTLR7, hTLR8 and hTLR9. *Eur Cytokine Netw*, 11(3):372-8. 2. **Jurk M, et al., 2002.** Human TLR7 or TLR8 independently confer responsiveness to the antiviral compound R-848. *Nat Immunol*, 3(6):499. 3. **Heil F, et al., 2004.** Species-specific recognition of single-stranded RNA via toll-like receptor 7 and 8. *Science*. 303(5663):1526-9.

### Handling Cells Upon Arrival

We strongly recommend that you propagate the cells, using the provided procedure, as soon as possible. This will ensure the best cell viability and assay performance. Frozen cells may be placed in liquid nitrogen until you are ready to thaw and propagate them, however, this may reduce cell viability.

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873  
Outside US: (+1) 858-457-5873  
Europe: +33 562-71-69-39  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)  
Website: [www.invivogen.com](http://www.invivogen.com)

### Product Warranty

InvivoGen warrants that cells shall be viable upon shipment from InvivoGen for a period of thirty days, provided they have been properly stored and handled during this period.

### 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.

293XL-hTLR8A cells should not be passaged more than 20 times to remain fully efficient. 293XL-hTLR8A cells should be maintained in Growth Medium as described below in the presence of Normocin™ (100 µg/ml) and the selective antibiotic, Blastidicin (10 µg/ml). Antibiotic pressure with Blastidicin is required to maintain the plasmid coding for hTLR8a.

### Quality control

Expression of hTLR8a gene was confirmed by RT-PCR. These cells were stimulated with various TLR8 Ligands. 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](mailto:info@invivogen.com)

### SAFETY CONSIDERATIONS

**Biosafety Level:2**

## HANDLING PROCEDURES

### Required Cell Culture Medium

- Growth Medium: DMEM, 4.5 g/l glucose, 10% (v/v) fetal bovine serum, 50 U/ml penicillin, 50 µg/ml streptomycin, 100 µg/ml Normocin™, 2 mM L-glutamine
- Freezing Medium: DMEM, 4.5 g/l glucose, 20% (v/v) fetal bovine serum, 50 U/ml penicillin, 50 µg/ml streptomycin, 100 µg/ml Normocin™, 2 mM L-glutamine, 10% (v/v) DMSO



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA

### 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 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 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- Aliquot 1 ml cells into cryogenic vials.

3- Place vials in a freezing container (Nalgene) 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 10 µg/ml Blasticidin.

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 cells grow to 100% confluency.

*Note: The response of 293XL-hTLR8A cells can be altered by the action of trypsin. Do not use trypsin to detach 293XL-hTLR8A cells.*

### TLR8 Stimulation

TLR8 stimulation can be assessed by determining the levels of IL-8 using an ELISA kit or by measuring the activation of NF-κB. InvivoGen has developed a simple and convenient method to evaluate TLR stimulation through NF-κB activation based on the use of an NF-κB-inducible SEAP reporter system (pNiFty-SEAP) and QUANTI-Blue™, a SEAP detection medium. Alternatively, InvivoGen provides HEK-Blue™ hTLR8 cells (cat. code hkb-htr8), a SEAP reporter cells line expressing the hTLR8 gene. Determine SEAP levels using a spectrophotometer at 620-655 nm.

### Day 1: Transfection of 293XL-hTLR8A cells with pNiFty-SEAP

1- Prepare pNiFty-SEAP/LyoVec™ complexes following the instructions provided in the technical data sheet of LyoVec™.

*Note: If using another transfection reagent, perform transfection according to the manufacturer's recommendations.*

2- Seed 50,000 cells per well of a flat-bottom 96-well plate in 200 µl Growth Medium.

3- Add 10 µl of pNiFty(2)-SEAP/LyoVec™ complexes per well.

4- Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 20-24 h.

### Day 2: TLR8 Stimulation

- Remove medium and replace with 180 µl of fresh Growth Medium which contains 10% (v/v) heat-inactivated FBS.

*Note: Some fetal bovine serum (FBS) may contain alkaline phosphatases that can interfere with SEAP quantification. To ensure that these thermosensitive enzymes are inactive, use heat-inactivated FBS (30 min at 56°C). Heat-inactivated FBS is also commercially available.*

- Add 20 µl of each sample per well of a 96-well plate.

- Add 20 µl of a positive control (such as ssRNA40/ LyoVec™, 50 µg/ml) in one well.

- Add 20 µl of a negative control (such as sterile, endotoxin-free water) in one well.

- Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 20-24 h.

### Day 3: Detection and Quantification of SEAP

- Prepare QUANTI-Blue™ following the instructions on the pouch.

- Add 180 µl of resuspended QUANTI-Blue™ per well of a 96-well plate.

- Add 20 µl of induced 293XL-hTLR8A Cells supernatant.

- Incubate the plate at 37°C incubator for 1-3 h.

- Determine SEAP levels using a spectrophotometer at 620-655 nm.

*Note: For faster reading or high-throughput applications we recommend the use of the one step HEK-Blue™ Detection growth medium. This medium allows for the combined growth of your cells and reading of SEAP activity.*

### Specificity of 293XL-hTLR8A Cells

As HEK293 cells express endogenous levels of TLR3, TLR5, and NOD1, 293XL-hTLR8A cells will respond to TLR3, TLR5 and NOD1 ligands. To ensure the specificity of the hTLR8 activation, we recommend that you perform experiments with the control cell line 293XL-null cells. This will avoid misleading results, due to direct activation of NF-κB via a non-hTLR8 pathway (e.g. TNFα activation of NF-κB).

## RELATED PRODUCTS

Product	Description	Catalog Code
Blasticidin (100 mg)		ant-bl-1
Normocin™		ant-nr-1
293XL-null (Control cell line)		293xl-null
pNiFty2-SEAP (NF-κB inducible reporter plasmid)		pnifty2-seap
LyoVec™ (Transfection reagent)		lyec-1
QUANTI-Blue™ (5 pouches)		rep-qb1
HEK-Blue™ Detection (2 pouches)		hb-det1
ORN02 / LyoVec™ (TLR8 ligand)		trl-orn2
ORN06 / LyoVec™ (TLR8 ligand)		trl-orn6
ssRNA40 / LyoVec™ (TLR8 ligand)		trl-rna40
ssPoly(U) / LyoVec™ (TLR8 ligand)		trl-lpu
CL075 (TLR7/8 ligand)		trl-c75
R848 (TLR7/8 ligand)		trl-r848

### TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

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

Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA