

### <u>Data Sheet</u> HVEM/NF-кВ Reporter Jurkat Recombinant Cell Line Catalog # 79310

#### **Product Description**

Recombinant clonal stable Jurkat T cell line expressing firefly luciferase gene under the control of 4 copies of NF-kB response elements with constitutive expression of human HVEM (Herpes Virus Entry Mediator, Tumor Necrosis Factor Receptor Superfamily Member 14, TNFRSF14, CD270, GenBank Accession No. NM\_003820). Following activation by human HVEM ligand LIGHT, NF-kB transcription factors bind to the DNA response elements to induce transcription of the luciferase reporter gene.

#### Background

HVEM is a bidirectional switch regulating T-cell activation in a costimulatory or coinhibitory fashion whose outcome depends on the binding partner. HVEM can act as both receptor and ligand, the binding of endogenous ligand LIGHT or agonist antibodies to HVEM delivers a costimulatory signal; whereas the binding of HVEM to BTLA (IgSF) or CD160 on Effector T cells delivers a coinhibitory signal. LIGHT/HVEM axis are co-stimulatory immune checkpoint molecules extensively studied for cancer immunotherapy; LIGHT, either in the soluble monomer form or homotrimer expressed on the cell surface, can activate NF- $\kappa$ B through binding to HVEM; we have shown that LIGHT expressing cells are more potent and efficacious than the soluble LIGHT for activating HVEM expressed on T cell surface.

#### Application

- Screen for agonists or antagonists of LIGHT-HVEM signaling in a physiological relevant cellular context
- Characterize T cell-mediated immune responses of HVEM and its interactions with LIGHT
- Screen co-stimulatory immune checkpoint molecules for cancer immunotherapy

#### **Host Cell**

Jurkat T cell

#### Format

Each vial contains ~  $2 \times 10^6$  cells in 1 ml of 10% DMSO in FBS.

#### Storage

Store in liquid nitrogen immediately upon receipt.

**Thaw Medium 2 (BPS Cat. #60184):** RPMI1640 medium (Life Technologies #A10491-01) supplemented with 10% FBS (Life Technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone #SV30010.01)



**Complete Growth Medium:** Thaw Medium 2 (BPS Cat. #60184) plus 1 mg/ml G418 (Thermo Fisher, Cat. # 11811031) and 200 µg/ml of Hygromycin B (Hyclone #SV30070.01)

Cells should be grown at 37°C with 5% CO2 using complete growth medium (Thaw Medium 2 with G418 and Hygromycin B).

#### **Recommended Culture Condition**

It is recommended to quickly thaw the frozen cells from liquid nitrogen in a  $37^{\circ}$ C water-bath, then transfer the entire contents of the vial to a tube containing 10 ml of Thaw Medium 2 (no G418 and Hygromycin B). Spin down the cells, remove supernatant and resuspend cells in prewarmed Thaw Medium 2 (no G418 and Hygromycin B). Transfer the resuspended cells to a T25 flask and incubate at  $37^{\circ}$ C in a 5% CO2 incubator. This cell line tends to grow more slowly than parental WT Jurkat cells. After 24 hours of culture, add an additional 3 - 4 ml of growth medium without antibiotics. At first passage, switch to complete growth medium (contains G418 and Hygromycin B). Cells should be split before they reach  $2x10^{6}$  cells/ml.

To passage the cells, dilute cell suspension into new culture vessels at no less than 0.1x106 cells/ml. Subcultivation ratio: 1:10 to 1:20 twice a week.

#### Mycoplasma Testing

This cell line has been screened using the MycoAlert<sup>™</sup> Mycoplasma Detection Kit (Lonza, #LT07-118) to confirm the absence of Mycoplasma contamination. MycoAlert Assay Control Set (Lonza, #LT07-518) was used as a positive control.

#### Application References

- 1. Chen, *et.al.* (2013) Molecular mechanisms of T cell co-stimulation and co-inhibition. *Nat Rev Immunol.* 2013 April; 13(4): 227–242.
- 2. Steinberg, *et.al.* (2014) The Signaling Networks of the Herpesvirus Entry Mediator (TNFRSF14) in Immune Regulation. *Immunol Rev. 2011 November; 244(1): 169–187*.
- 3. Rio, *et.al.* (2014) Therapeutic blockade of LIGHT interaction with HVEM and LTβR attenuates in vivo cytotoxic allogeneic responses. *Transplantation.* 2014 December 15; 98(11): 1165–1174.
- 4. Shui *et.al.* (2014) HVEM is a TNF Receptor with Multiple Regulatory Roles in the Mucosal Immune System. *Immune Netw. 2014 Apr; 14(2): 67–72.*
- 5. Steinberg, *et.al.* (2009) Regulating the mucosal immune system: the contrasting roles of LIGHT, HVEM, and their various partners. *Semin. Immunopathol.* 31: 207-221.
- 6. Ware et.al. (2009) Targeting the LIGHT-HVEM pathway. Adv. Exp. Med. Biol. 647:146

#### Assay Principle



6042 Cornerstone Court W, Ste B San Diego, CA 92121 **Tel:** 1.858.829.3082 **Fax:** 1.858.481.8694 **Email:** info@bpsbioscience.com



**Figure 1. Expression of human HVEM validated by flow cytometry.** Flow cytometry showed PE-conjugated anti-human HVEM antibody (BioLegend, #318806) detects HVEM-positive clonal population (clone B15) (green), using wild-type Jurkat cells as a negative control (red).

#### **Vector and Sequence**

Human HVEM (NM\_003820.2) was cloned into pIRESHyg.



#### a.a. sequence:

MEPPGDWGPPPWRSTPKTDVLRLVLYLTFLGAPCYAPALPSCKEDEYPVGSECCPKCSPGYR VKEACGELTGTVCEPCPPGTYIAHLNGLSKCLQCQMCDPAMGLRASRNCSRTENAVCGCSPG HFCIVQDGDHCAACRAYATSSPGQRVQKGGTESQDTLCQNCPPGTFSPNGTLEECQHQTKC SWLVTKAGAGTSSSHWVWWFLSGSLVIVIVCSTVGLIICVKRRKPRGDVVKVIVSVQRKRQEAE GEATVIEALQAPPDVTTVAVEETIPSFTGRSPNH

#### Materials Required but Not Supplied

- Assay Medium: Thaw Medium 2 (BPS Cat. #60184)
- LIGHT, His-Tag (Human) (BPS Cat. #71266)
- LIGHT-CHO Recombinant Cell Line (BPS Cat. # 79262)
- 96-well tissue culture-treated white clear-bottom assay plate
- One-Step luciferase assay system (BPS bioscience # 60690) or other luciferase reagents for measuring firefly luciferase activity
- Luminometer

#### Assay Protocol

- 1. Harvest HVEM/NFκB Reporter\_Jurkat Recombinant Cell Line (effector cells) from culture in growth medium and seed cells at a density of 35,000 cells per well into white clear-bottom 96-well microplate in 90 μl of assay medium.
- Dilute the ligand (LIGHT or LIGHT/CHO cells) and agonist/antagonist (e.g. anti-LIGHT or anto-HVEM Ab) in assay medium. Add 10 μl of diluted ligand and antibody to the wells. Add 10 μl of assay medium to control wells. Add 100 μl of assay medium to cell-free control wells (for determining background luminescence). Set up each treatment in at least triplicate.
- 3. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 6 hours.
- 4. Perform luciferase assay using the ONE-Step luciferase assay system: Add 100 μl of One-Step Luciferase reagent per well and rock at room temperature for ~30 minutes. Measure luminescence using a luminometer. *If using luciferase reagents from other vendors, follow the manufacturer's assay protocol.*
- 5. Data Analysis: Obtain the background-subtracted luminescence by subtracting the average background luminescence (cell-free control wells) from the luminescence reading of all wells. The fold induction of NF-κB luciferase reporter expression = background-subtracted luminescence of stimulated well / average background-subtracted luminescence of unstimulated control wells.

#### **Figure 2.** HVEM/NF<sub>K</sub>B-Luc Reporter activities stimulated by different ligands The results are shown as fold induction of NF-κB luciferase reporter expression.



# LIGHT-HVEM Co-stimulatory cell-based reporter assay (Fold induction, n=3)



■ Series1 ■ Series2 ■ Series3

Stimulation by	Fold induction	
TNFa	5.3136	
sLIGHT	7.6818	

LIGH\_CHO 19.456

## Figure 3. Dose Response of HVEM/NF $\kappa$ B-Luc Reporter activities stimulated by different soluble LIGHT and LIGHT expresses on CHO cell surface



The results are shown as fold induction of NF-KB luciferase reporter expression.

## LIGHT-HVEM reporter assay dose response curves



License Disclosure: Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit the use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use, and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact sales@bpsbioscience.com for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

Related Products				
	Cat. #	Size		
ONE-Step <sup>™</sup> Luciferase Assay System	60690-1	10 ml		
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIA	GNOSTIC OR TH	IERAPEUTIC USE.		
To place your order, please contact us by Phone 1.858.829.	3082 Fax 1.858.	481.8694		
Or you can Email us at: info@bpsbioscience.com				
Please visit our website at: www.bpsbioscie	<u>ence.com</u>			



ONE-Step <sup>™</sup> Luciferase Assay System	60690-2	100 ml
LIGHT, His-Tag (Human)	71266	100 µg
BTLA:HVEM[Biotinylated] Inhibitor Screening Assay Kit	72008	96 rxns
Thaw Medium 2	60184	100 ml
LIGHT-CHO Recombinant Cell Line	79262	2 vials