

**Description**

B2M (Beta-2-Microglobulin) has been genetically removed by CRISPR/Cas9 genome editing from THP-1 cells .

**Background**

THP-1 cells are monocytes, derived from an acute monocytic leukemia patient. Beta-2-Microglobulin is a required component of Major Histocompatibility Complex (MHC) class 1 molecules, which present peptide fragments from within the cell to cytotoxic T cells as part of the adaptive immune system. B2M plays an essential role both in governing MHC class I molecule stability and in promoting antigen binding and presenting the antigen to CD3/TCR complex of CD8+ T cells.

**Application**

1. Study the consequences of B2M knock-down
2. Study T cell activation, antigen presentation, and immune responses

**Materials Provided**

| Components              | Format  |
|-------------------------|---|
| 2 vials of frozen cells | Each vial contains $2 \times 10^6$ cells in 1 ml of cell freezing medium (BPS Bioscience, #79796) |

**Parental Cell Line**

THP-1, human monocyte, suspension

**Mycoplasma Testing**

The cell line has been screened to confirm the absence of Mycoplasma species.

**Materials Required but Not Supplied**

These materials are not supplied with the cell line but are necessary for cell culture and cellular assays. BPS Bioscience's reagents are validated and optimized for use with this cell line and are highly recommended for best results. Media components are provided in the Media Formulations section below.

**Media Required for Cell Culture**

| Name          | Ordering Information                  |
|---------------|---------------------------------------|
| Thaw Medium 2 | <a href="#">BPS Bioscience #60184</a> |

**Storage Conditions**

Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a  $-80^{\circ}\text{C}$  freezer for long term storage. Contact technical support at [support@bpsbioscience.com](mailto:support@bpsbioscience.com) if the cells are not frozen in dry ice upon arrival.

**Media Formulations**

For best results, it is *highly recommended* to use these validated and optimized media from BPS Bioscience. Other preparations or formulations of media may result in suboptimal performance.



Note: Thaw Media do not contain selective antibiotics. Cells should be grown at  $37^{\circ}\text{C}$  with 5%  $\text{CO}_2$ . BPS Bioscience's cell lines are stable for at least 15 passages when grown under proper conditions.

### Media Required for Cell Culture

Thaw Medium 2 (BPS Bioscience, #60184):

RPMI 1640 medium supplemented with 10% FBS, 1% Penicillin/Streptomycin

### Cell Culture Protocol

#### Cell Thawing

1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. As soon as the cells are thawed (it may be slightly faster or slower than 60 seconds), quickly transfer the entire contents of the vial to a tube containing 10 ml of pre-warmed Thaw Medium 2.

**Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.**

2. Immediately spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed Thaw Medium 2.
3. Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 24 hours of culture, check for cell viability. For a T25 flask, add 3-4 ml of Thaw Medium 2, and continue growing in a 5% CO<sub>2</sub> incubator at 37°C until the cells are ready to passage.
5. Cells should be passaged before they reach a density of 2 x 10<sup>6</sup> cells/ml. At first passage and subsequent passages, use Thaw Medium 2.

#### Cell Passage

Dilute the cell suspension into new culture vessels before they reach a density of 2 x 10<sup>6</sup> cells/ml, at no less than 0.2 x 10<sup>6</sup> cells/ml of Growth Medium 2. The sub-cultivation ratio should maintain the cells between 0.2 x 10<sup>6</sup> cells/ml and 2 x 10<sup>6</sup> cells/ml.

#### Cell Freezing

1. Spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cell pellet in 4°C Freezing Medium (BPS Bioscience #79796, or 10% DMSO + 90% FBS) at a density of ~2 x 10<sup>6</sup> cells/ml.
2. Dispense 1 ml of cell aliquots into cryogenic vials. Place the vials in an insulated container for slow cooling and store at -80°C overnight.
3. Transfer the vials to liquid nitrogen the next day for storage.



Note: It is recommended to expand the cells and freeze at least 10 vials at an early passage for future use.

**A. Validation Data**



Figure 1. Genomic Sequencing of B2M in the B2M Knockout THP-1 Cell Line. Genomic DNA from the B2M Knockout THP-1 cells were isolated and sequenced. The PAM (Protospacer Adjacent Motif) is shown in blue, the sgRNA (synthetic guide RNA) is shown in green, and the Indels (Insertions / Deletions) in the two B2M alleles are highlighted in red. The B2M genomic DNA is labeled as Ref.

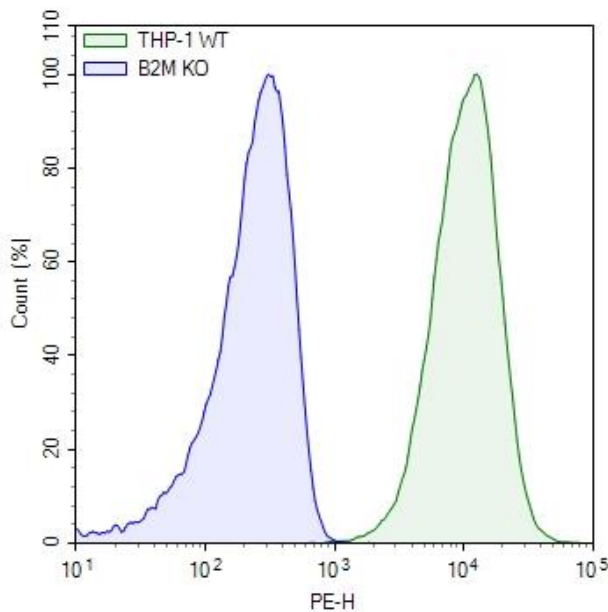


Figure 2. B2M Expression in B2M knockout THP-1 cells. Flow cytometry was performed using a PE-labeled anti-human B2M antibody (BioLegend, #395703). Parental THP-1 cells (green) were compared to B2M Knockout THP-1 cells (blue). The Y-axis is the % cell number. The X-axis is the intensity of PE.

**Sequence**

Human beta-2-microglobulin (B2M), mRNA, NCBI Reference Sequence: NM\_004048.4, with the sgRNA targeting sequence underlined:

ATTCCTGAAGCTGACAGCATTTCGGGCCGAGATGTCTCGCTCCGTGGCCTTAGCTGTGCTCGCGCTACTCTCTCTTTCTGGCCT  
 GGAGGCTATCCAGCGTACTCCAAAGATTCAGGTTTACTCACGTCATCCAGCAGAGAATGGAAAGTCAAATTCCTGAATTGC  
 TATGTGTCTGGGTTTCATCCATCCGACATTGAAGTTGACTTACTGAAGAAATGGAGAGAGAATTGAAAAAGTGGAGCATTCA  
 GACTTGTCTTTAGCAAGGACTGGTCTTTCTATCTCTTGTACTACACTGAATTCACCCCCACTGAAAAAGATGAGTATGCCTG  
 CCGTGTGAACCATGTGACTTTGTACAGCCCAAGATAGTTAAGTGGGATCGAGACATGTAAGCAGCATCATGGAGGTTTGA  
 AGATGCCGCATTTGGATTGGATGAATTCAAATCTGCTTGCTTGCTTTTAAATATTGATATGCTTATACACTTACACTTTATG  
 CAAAAATGTAGGGTTATAATAATGTTAACATGGACATGATCTTCTTTATAATTCTACTTTGAGTGCTGTCTCCATGTTTATG  
 GTATCTGAGCAGGTTGCTCCACAGGTAGCTCTAGGAGGGCTGGCAACTTAGAGGTGGGGAGCAGAGAATTCTCTTATCCAA  
 CATCAACATCTTGGTCAGATTTGAACTCTTCAATCTCTTGCCTCAAAGCTTGTTAAGATAGTTAAGCGTGCATAAGTTAACT  
 TCCAATTTACATACTGCTTAGAATTTGGGGGAAAATTTAGAAATATAATTGACAGGATTATTGGAAATTTGTTATAATGAA  
 TGAAACATTTTGTATATAAGATTCATATTTACTTCTTATACATTTGATAAAGTAAGGCATGGTTGTGGTTAATCTGGTTTATT  
 TTTGTTCCACAAGTTAAATAAATCATAAACTTGA

**License Disclosure**

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

**Troubleshooting Guide**

Visit [bpsbioscience.com/cell-line-faq](https://bpsbioscience.com/cell-line-faq) for detailed troubleshooting instructions. For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

**Related Products**

| <i>Products</i>  | <i>Catalog #</i> | <i>Size</i> |
|--|------------------|-------------|
| TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line | 78364            | 2 vials     |
| B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line     | 78363            | 2 vials     |
| B2M Knockout Jurkat Cell Line                              | 78342            | 2 vials     |
| B2M (Human) CRISPR/Cas9 Lentivirus (Integrating)           | 78340            | 500 µl x 2  |
| B2M (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)       | 78341            | 500 µl x 2  |