# B-27™ Minus Insulin

## Description

B-27<sup>™</sup> Minus Insulin is the complete B-27<sup>™</sup> Supplement without insulin, and it is used to support growth of hippocampal and other CNS neurons in applications where the presence of insulin causes interference.

Product	Catalog no.	Amount	Storage	Shelf life*
B-27™ Minus Insulin (50X), liquid	A18956-01 A18956-02	10 mL 100 mL	–20°C to –5°C; Protect from light	12 months

\* Shelf life duration is determined from Date of Manufacture.

#### **Product use**

For Research Use Only. Not for use in diagnostic procedures.

### Safety information

Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Prepare medium

Neurobasal<sup>™</sup>-A Medium is recommended for primary neuronal cultures. Neurobasal<sup>™</sup>-A Medium requires supplementation with GlutaMAX<sup>™</sup>-I supplement (200 mM) and B-27<sup>™</sup> Minus Insulin (50X) prior to use. Thaw B-27<sup>™</sup> Minus Insulin overnight at 4°C.

- 1. As eptically add GlutaMAX  $^{\rm \tiny M}\mbox{-I}$  to 0.5 mM final concentration (2.5 mL/L) to the medium before use.
- 2. As eptically add 2% B-27  $^{\mbox{\tiny M}}$  Minus Insulin (20 mL/L) to the medium before use.

**Note:** You may aliquot the remaining B-27<sup>™</sup> Minus Insulin into working volumes and store at -20°C to -5°C. Thaw aliquots as needed. Do not freeze-thaw B-27<sup>™</sup> Minus Insulin more than twice.

- For primary rat hippocampus neuron cultures, the complete Neurobasal<sup>™</sup>-A Medium Minus Insulin (prepared from the previous steps) requires additional supplementation with 25 µM L-Glutamate up to the fourth day in culture.
- 4. Once supplemented, the complete Neurobasal<sup>™</sup>-A Medium Minus Insulin is stable for up to one week when stored in the dark at 2°C to 8°C.

#### Use

- Use B-27<sup>™</sup> Minus Insulin to supplement Neurobasal<sup>™</sup> Medium for optimal viability and long-term survival of pre-natal and embryonic neuronal cells.
- Use B-27<sup>™</sup> Minus Insulin to supplement Neurobasal<sup>™</sup>-A Medium for optimal viability and long-term survival of post-natal and adult brain neuronal cells.
- B-27<sup>™</sup> Minus Insulin when used as a supplement to RPMI 1640 has been demonstrated to support differentiation of pluripotent stem cells into cardiomyocytes.
- Areas of application include studies of neural development, neural differentiation, diabetic neuropathy, and pluripotent stem cell differentiation.

## Cell culture procedure

The following procedure has been tested on freshly isolated 18-day gestation rat hippocampal and cortical neurons, Gibco<sup>™</sup> Primary Rat Cortex Neurons, and Gibco<sup>™</sup> Primary Rat Hippocampus Neurons.

1. Coat culture surface (German glass or cell culture grade plastics) with a sterile 0.05 mg/mL solution of cold poly-D-lysine in water at 0.15 mL/cm<sup>2</sup> surface area and incubate for 1 hour at ambient temperature.

- 2. Remove poly-D-lysine solution, and rinse twice with sterile distilled water. (Rinse thoroughly, since poly-D-lysine can be toxic to the cells). Leave the plates uncovered in the cell culture hood until the wells are completely dry. Plates can be used immediately once dry, or they can be stored dry at 4°C for up to 2 weeks.
- 3. Isolate primary rat neurons or thaw cryopreserved primary rat neurons according to standard laboratory procedure or instructions supplied with the cells (See **Recovery and Culture of Cryopreserved Neurons**).
- Plate cells in pre-warmed (37°C) complete Neurobasal<sup>™</sup>-A Medium Minus Insulin (prepared as described above) at a suggested density of 160 cells/mm<sup>2</sup>, or another optimized density if required.

**Note:** For hippocampal neurons, use the complete medium supplemented with 25 µM L-Glutamate, see **Prepare Medium**.

- 5. Incubate the culture dish at  $36^{\circ}$ C to  $38^{\circ}$ C in a humidified atmosphere of 5% CO<sub>2</sub> (in air is acceptable, but 9% oxygen with 5% CO<sub>2</sub> is preferable).
- 6. After 4–24 hours of incubation, aspirate half of the medium and replace with same volume of fresh medium. Return the plate to the incubator.
- 7. **Non-hippocampal cultures:** Four days after plating, feed the cultures by aspirating half of the medium from each well and replacing with same volume of fresh medium. Repeat every three days thereafter.

**Hippocampal cultures:** Three days after plating, replace half of the medium with complete medium **without** L-Glutamate. Repeat every three days thereafter.

Note: Improved long-term survival of hippocampal neurons may be obtained by the addition of 25  $\mu M$  2-mercaptoethanol.

## Isolate primary fetal neurons

The following procedure is recommended for cultured 18-day embryonic rat hippocampal or cortical neurons.

- 1. Dissect cortex or hippocampi pairs from rat embryos at Day 18 of gestation (E18).
- 2. Collect all the tissue in a conical tube containing Hibernate-E complete medium. Leave the tissue in this tube (1 pair/2 mL) until all the dissections are completed.
- 3. Let the tissue settle to the bottom of the tubes and then carefully remove supernatant leaving only the tissue covered by a minimum amount of medium.
- Enzymatically digest the tissue in Hibernate-E, without Ca<sup>2+</sup> (BrainBits<sup>™</sup> LLC, Cat. no. HE-Ca) medium containing 2 mg/mL filter sterilized papain at 30°C for 30 minutes with gentle shaking of the tube every 5 minutes (2 pairs/mL).
- 5. Restore divalent cations with 2 volumes of Hibernate-E complete medium.

- 6. Allow non-dispersed tissue to settle for 2 minutes and then transfer the supernatant to a 15-mL tube and centrifuge for 5 minutes at  $150 \times g$ .
- Gently resuspend the pellet in 1 mL complete Neurobasal<sup>™</sup>-A Medium Minus Insulin and take an aliquot (e.g., 10 µL) for cell counting. Proceed to Recovery and Culture of Cryopreserved Neurons, steps 8–10.

Note: Gibco<sup>™</sup> Primary Rat Cortex and Rat Hippocampus Neurons (isolated from day-18 Fisher 344 rat embryos and cryopreserved in a medium containing 10% DMSO) are quality ready-to-use alternative to freshly isolated neurons (see **Recovery and Culture of Cryopreserved Neurons**).

#### Recovery and culture of cryopreserved neurons

**Important:** Primary neuronal cells will adhere to bare plastic and glassware; to maximize cell recovery and yield we recommend prerinsing all plastic and glassware with complete medium before use. **Do not** vortex or centrifuge cells at any time during this procedure as cells are extremely fragile upon recovery from cryopreservation. We recommend thawing one vial at a time. Transfer cryovial from liquid nitrogen storage to 37°C water bath minimizing handling time. A small amount of liquid nitrogen in an ice bucket can be used to transport the vials from liquid nitrogen to the water bath.

- 1. Rinse a sterile 15-mL conical culture tube with complete Neurobasal<sup>™</sup>-A Medium Minus Insulin and leave in the hood prior to thawing cells.
- 2. If removing vial from liquid nitrogen storage, twist cap slightly to release pressure and then retighten cap.
- 3. Rapidly thaw (<2 minutes) frozen vial by gently swirling in a 37°C water bath. Remove from water bath when only one tiny ice crystal is left (vial should still be cold to the touch).
- 4. Transfer the vial into the hood and disinfect with 70% isopropyl alcohol. Collect the liquid to the bottom of the vial by gently tapping the vial on the hood's surface.
- 5. Use a pre-rinsed P-1000 pipette tip to very gently transfer the cells to the pre-rinsed 15-mL conical tube.
- 6. Rinse the cryovial with 1 mL of pre-warmed complete Neurobasal<sup>™</sup>-A Medium Minus Insulin and extremely slowly add to the cells in the 15-mL tube at the rate of one drop per second. Mix by gentle swirling after each drop. Do not add the full amount of media to the tube at once. This may lead to decreased cell viability due to osmotic shock.
- 7. Slowly (dropwise) add an additional 2 mL of pre-warmed complete Neurobasal<sup>™</sup>-A Medium Minus Insulin to the tube (for a total suspension volume of 4 mL). Mix the suspension very gently with P-1000 pipette without creating any air bubbles.
- 8. Add 10  $\mu$ L of cell suspension to a microcentrifuge tube containing 10  $\mu$ L of 0.4% Trypan blue, using a pre-rinsed tip. Mix only by gently tapping the tube. Determine the viable cell density using a manual (i.e., hemocytometer) counting method. The viability of thawed cells should be >50%.
- Plate ~1 × 10<sup>5</sup> cells (see Cell Culture Procedure, steps 4–5) (or desired cell density) per well in a poly-D-lysine coated 48-well plate (see Cell Culture Procedure, steps 1–2). Dilute cell suspension to 500 µL per well by adding pre-warmed complete Neurobasal<sup>™</sup>-A Medium Minus Insulin.
- 10. Follow **Cell Culture Procedure**, steps 6–7, to maintain neuronal cell cultures. Incubate at  $36^{\circ}$ C to  $38^{\circ}$ C in a humidified atmosphere of 5% CO<sub>2</sub> (in air is acceptable but 9% oxygen with 5% CO<sub>2</sub> is preferable).

## **Related products**

Product	Cat. no.
B-27 <sup>™</sup> Supplement (50X), liquid	17504
Neurobasal <sup>™</sup> Medium (1X), liquid	21103
Neurobasal <sup>™</sup> -A Medium (1X), liquid	10888
RPMI 1640 Medium (1X), liquid	11875
GlutaMAX <sup>™</sup> -I (100X), liquid	35050
Hibernate E	A12476
Primary Rat Cortex Neurons, $1 \times 10^6$ viable cells/vial $4 \times 10^6$ viable cells/vial	A10840-01 A10840-02
Primary Rat Hippocampus Neurons, $1 \times 10^6$ viable cells/vial	A10841
2-mercaptoethanol (1000X), liquid	21985
Countess <sup>™</sup> II Automated Cell Counter	AMQAX1000
Trypan Blue Stain	15250

#### Explanation of symbols and warnings

The symbols present on the product label are explained below:

5	1	1	1	
$\wedge$		×	STERILE A	Í
Caution, consult accompanying documents	Temperature Limitation	Keep away from light	Sterilized using aseptic processing techniques	Consult instructions for use
LOT	REF	444		Read SDS
Batch Code	Catalog number	Manufacturer	Use By:	Read Safety Data Sheet

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **www.lifetechnologies.com/termsandconditions**. If you have any questions, please contact Life Technologies at **www.lifetechnologies.com/support**.

#### Important licensing information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

#### References

- 1. Singh *et al.* (2012) Resistance to trophic neurite outgrowth of sensory neurons to insulin. J Neurochem *121*, 263–276.
- Sekine *et al.* (2012) Highly efficient generation of definitive endoderm lineage from human induced pluripotent stem cells. Transpl Proceed 44, 1127–1129.
- 3. Lian *et al.* (2012) Robust cardiomyocyte differentiation from human pluripotent stem cells via temporal modulation of canonical Wnt signaling. PNAS *109*, E1848–1857.
- Lian *et al.* (2013) Directed cardiomyocyte differentiation from human pluripotent stem cells by modulating Wnt/β-catenin signaling under fully defined conditions. Nature Protocols 8, 162–175.

For additional technical information such as Safety Data Sheets (SDS), Certificates of Analysis, visit **www.thermofisher.com/support**. For further assistance, email **techsupport@lifetech.com** 

© 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. B-27 is a registered trademark of Southern Illinois University. BrainBits is a registered trademark of BrainBits LLC.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

