

ab315369 – Human Cortical Astrocytes (Male, WC-30) 96-well Astrocyte Culturing

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab315369>

Storage and Stability

Immediately transfer the vial of neurons to liquid nitrogen upon receipt. Transfer the vials of supplements to a -20°C freezer. The supplements can be stored at -20°C for up to 6 months. Alternatively, the supplements can be stored at -80°C for up to 18 months.

Materials Supplied

Item	Quantity	Storage Condition
One vial of 1 million cryopreserved human astrocytes	500 µL	-196°C
BrainFast Astro Supplement (1000x)	1 vial	-20°C or -80°C
BrainFast SK Supplement (1000x)	1 vial	-20°C or -80°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- DMEM/F12 Medium (Life Technologies #11330-032)
- Neurobasal Medium (Life Technologies #21103-049)
- N2 Supplement (Life Technologies #17502-048)
- GlutaMAX (Life Technologies #35050-061)
- Fetal Bovine Serum (Life Technologies #A3160601)
- PDL-Coated 96-Well Plates

PROCEDURE

If performing spheroid cultures, the recommended plates are either S-Bio PrimeSurface 3D Culture Spheroid (ULA) plates (CAT: MS-9096WZ) or Nunclon Sphera 96-Well, U-Shaped-Bottom Microplate (CAT: 174925). When seeding these plates cells can be seeded between 100 and 10,000 cells/well successfully.

Thawing and Seeding the Astrocytes

1. Gather the components for the Seeding Medium according to the recipe below.
2. Each vial of cells requires approximately 20 mL of seeding medium. Working in a cell culture hood (biological safety cabinet), combine all components in an appropriately sized sterile container. Allow the medium to equilibrate to room temperature for 15 minutes. Do not warm the medium in a 37°C water bath.
3. Remove the cryovial from the liquid nitrogen and place in a 37°C water bath. To minimize contamination, avoid submerging the cap. Gently move the vial within the bath to increase the rate of thawing.
4. As soon as the last of the ice melts, which will take ~75-90 seconds, remove the vial from the water bath. Disinfect the vial by spraying it with 70% ethanol and transfer it to the cell culture hood.
5. Slowly add 500 µL of seeding medium to the vial at a rate of ~1 drop/s using a 1 mL pipette tip. This process should take about 30 seconds.
6. Gently transfer all contents (1mL total) from the vial to a new sterile 50 mL conical tube.

7. To collect any residual cells, gently add another 1 mL of seeding medium to the vial and then transfer to the conical tube.
8. Slowly add an additional 3 mL of seeding medium to the 50 mL conical tube using a 10 mL serological pipette. Gently swirl the conical tube while adding the medium. This process should take about 1 minute.
9. To count the cells, gently swirl the conical tube again and remove 10 µL from the cell suspension. Count the number of viable cells per mL with a hemocytometer using the trypan blue exclusion method to identify dead/viable cells.
10. The recommended seeding density is 20,000 – 25,000 viable cells/well for a 96-well plate (~62,500 – 78,000 viable cells/cm²). Use the following equation to determine the volume of cell suspension needed for each 96-well plate:

volume of cell suspension needed (mL) = $(2.4 - 3.0 \times 10^6 \text{ cells}) / (\text{viable cells per mL})$

11. In a separate 50 mL conical tube, add the calculated volume of cell suspension needed, and then add enough medium to obtain a final volume of 12 mL. For example, if the volume of cell suspension needed is 2 mL, combine 2 mL of cell suspension with 10 mL of medium.
12. Mix completely and then plate 100 µL/well (20,000 – 25,000 cells/well) onto a PDL-coated 96-well plate using a multi-channel pipettor or liquid handler. Throughout the seeding process, be careful not to move or agitate the plate as this may lead to uneven attachment. After seeding, do not immediately transfer the plate to the incubator. Leave it in the hood for 15 minutes to allow the cells to settle to the bottom of the well. After 15 minutes, very gently transfer the plate to a humidified incubator at 37°C with 5% CO₂. Day of cell plating is designated as Day 0.

Δ Note: Entire thawing and plating process should not exceed 2 hours; post-thaw viability and overall cell health will be severely impacted and lead to an unsuccessful culture.

Day 1 Medium Change

1. On Day 1 (24 hours after seeding), prepare fresh medium according to the Day 1 Medium composition below.
2. Very gently remove the entire 100 µL from each well and gently add 100 µL of the Day 1 Medium back to each well. Take care to ensure that the astrocytes do not dry out at any point in the medium changing process. We recommend changing one row or column at a time using a multichannel pipettor. Once complete, return the plate to the incubator.

Day 4 Medium Addition

1. On Day 4 (96 hours after seeding), prepare fresh medium according to the Day 4 Medium composition below.
2. Gently add 100 µL of Day 4 Medium to each well for a final total volume of 200 µL/well.

Day 7 and Onward Medium Changes

1. Change half the medium (100 µL/well) twice weekly using the recommended Maintenance Medium.
2. Gently withdraw 100 µL/well, then add 100 µL/well fresh medium to the entire plate.
3. The astrocytes can be maintained viable and adherent in culture under the above conditions for at least 4 weeks post-seeding.

Media Compositions

Seeding Medium

	Component	Stock Conc.	Final Conc.	1 plate volumes	2 plate volumes	5 plate volumes
1	DMEM/F12 Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
2	Neurobasal Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
3	N2 Supplement	100X	1X	200 µL	400 µL	1 mL
4	GlutaMAX	200 mM	0.5 mM	50 µL	100 µL	250 µL
5	BrainFast Astro Supplement (1000x)	1000X	1X	20 µL	40 µL	100 µL
6	Fetal Bovine Serum	1X	1%	200 µL	400 µL	1 mL

Day 1 Medium

	Component	Stock Conc.	Final Conc.	1 plate volumes	2 plate volumes	5 plate volumes
1	DMEM/F12 Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
2	Neurobasal Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
3	N2 Supplement	100X	1X	200 µL	400 µL	1 mL
4	GlutaMAX	200 mM	0.5 mM	50 µL	100 µL	250 µL
5	BrainFast Astro Supplement (1000x)	1000X	1X	20 µL	40 µL	100 µL
6	BrainFast SK Supplement (1000x)	1000X	1X	20 µL	40 µL	100 µL

Day 4 Medium

	Component	Stock Conc.	Final Conc.	1 plate volumes	2 plate volumes	5 plate volumes
1	DMEM/F12 Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
2	Neurobasal Medium	1X	0.5X	9.8 mL	19.6 mL	49 mL
3	N2 Supplement	100X	1X	200 µL	400 µL	1 mL
4	GlutaMAX	200 mM	0.5 mM	50 µL	100 µL	250 µL
5	BrainFast Astro Supplement (1000x)	1000X	1X	20 µL	40 µL	100 µL

Maintenance Medium

	Component	Stock Conc.	Final Conc.	1 plate volumes	2 plate volumes	5 plate volumes
1	DMEM/F12 Medium	1X	0.5X	4.9 mL	9.8 mL	24.5 mL
2	Neurobasal Medium	1X	0.5X	4.9 mL	9.8 mL	24.5 mL
3	N2 Supplement	100X	1X	100 µL	200 µL	500 µL
4	GlutaMAX	200 mM	0.5 mM	25 µL	50 µL	125 µL

Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at the time of printing. For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)