

## High Performance Cell Culture Media and Reagents



### Classical Media

Order No.	Description	Quantity
<b>MEM</b>		

MEM was developed by Harry Eagle, and is one of the most widely used basal media for synthetic cell culture media formulations. MEM is a modification of Basal Medium Eagle (BME), containing a higher concentration of essential nutrients than BME. MEM is frequently used to cultivate a variety of cells grown in monolayers.

01-025-1A	MEM-Eagle without L-Glutamine	500 mL
01-025-1B	MEM-Eagle without L-Glutamine	100 mL
01-025-5A	MEM-Eagle X10 without L-Glutamine	500 mL
01-040-1A	MEM-Eagle with Non-Essential Amino Acids, without L-Glutamine	500 mL
01-042-1A	MEM-Alpha, Low Glucose	500 mL
01-043-1A	MEM-Alpha, High Glucose	500 mL
01-045-1A	MEM-Eagle for Suspension cultures, 500 mL	500 mL
01-325-1B	MEM-Eagle Amino Acids Solution x 50	100 mL
01-340-1A	MEM Eagle Non-Essential Amino Acids Solution x 100	500 mL
01-340-1B	MEM Eagle Non-Essential Amino Acids Solution x 100	100 mL
01-015-5A	Basal Medium - Eagle x 10	500 mL

### DMEM

DMEM is a modification of Basal Medium Eagle (BME), with a higher concentration of amino acids and vitamins than BME and additional supplemental components. DMEM is one of the most common basal media used in cell culture, and is the basis for many specialty media and serum-free formulations.

01-050-1A	DMEM, Low Glucose	500 mL
01-052-1A	DMEM, High Glucose with L-Glutamine	500 mL
01-053-1A	DMEM, High Glucose without L-Glutamine and Phenol Red	500 mL
01-055-1A	DMEM, High Glucose without L-Glutamine	500 mL
01-055-4A	DMEM X5 High Glucose without L-Glutamine	500 mL
01-055-9A	DMEM X2, High Glucose without L-Glutamine	500 mL

Leverage a Complete Portfolio of Cell Culture Media and Reagents to Support your Research Workflow

Ensure quality cell culture with an extensive range of classical media and reagents for diverse mammalian cell cultures. From establishing and maintaining basic cell cultures to more advanced cell models, Sartorius offers a range of optimized products designed to meet your research needs.

Discover More:



01-056-1A	DMEM, High glucose with Alanine Glutamine	500 mL
01-057-1A	DMEM without Glucose and L-Glutamine	500 mL
<b>Ham's Nutrient Mixtures</b>		
Ham's F-10 and F-12 Nutrient Mixtures were originally developed for serum-free culture of CHO cells, and are now standard reagents in a cell culture lab. Ham's F-10 and F-12 are used for a variety of cell culture applications with and without the addition of a serum supplement.		
01-090-1A	F-10 (HAM) Nutrient Mixture with L-Glutamine	500 mL
01-090-1B	F-10 (HAM) Nutrient Mixture with L-Glutamine	100 mL
01-095-1A	F-12 (HAM) Nutrient Mixture	500 mL
<b>RPMI Medium</b>		
RPMI 1640 Medium was specifically developed for the long-term culture of blood cells, including normal and abnormal human leukocytes (e.g. neoplastic white blood cells). With the addition of serum, RPMI 1640 is also used as a general culture medium for hybridomas.		
01-100-1A	RPMI 1640 Medium with L-Glutamine	500 mL
01-100-1B	RPMI 1640 Medium with L-Glutamine	100 mL
01-101-1A	RPMI 1640 Medium without Glucose and L-Glutamine	500 mL
01-103-1A	RPMI 1640 Medium without L-Glutamine and Phenol Red	500 mL
01-104-1A	RPMI 1640 Medium without L-Glutamine	500 mL
01-104-1B	RPMI 1640 Medium without L-Glutamine	100 mL
01-106-1A	RPMI 1640 Medium with L-Glutamine and HEPES	500 mL
01-106-1B	RPMI 1640 Medium with L-Glutamine and HEPES	100 mL
<b>Other Classical Cell Culture Media</b>		
01-110-1A	Waymouth's MB 752/1 Medium	500 mL
01-170-1A	DMEM/F-12 (Ham) 1:1	500 mL
01-060-1A	GLASGOW MEM	500 mL
01-058-1A	Iscoe's Modified Dulbecco Medium with L-Glutamine	500 mL
01-059-1A	MCDB 153 medium (modified)	500 mL
<b>Serum Alternatives</b>		
15950-017	Ultrosor G	20 mL



## Buffers & Lab Waters

Order No.	Description	Quantity
<b>Earle's Balanced Salt Solutions (EBSS)</b>		
02-010-1A	Earle's Balanced Salt Solution	500 mL
02-011-1A	Earle's Balanced Salt Solution without phenol red	500 mL
<b>Hanks' Balanced Salt Solutions (HBSS)</b>		
02-015-1A	Hanks' Balanced Salt Solution	500 mL
02-015-1B	Hanks' Balanced Salt Solution	100 mL
02-015-5A	Hanks' Balanced Salt Solution X10	500 mL
02-016-1A	Hanks' Balanced Salt Solution without phenol red	500 mL
02-017-1A	Hanks' Balanced Salt Solution without Calcium and Magnesium	500 mL
02-017-1B	Hanks' Balanced Salt Solution without Calcium and Magnesium	100 mL
02-017-5A	Hanks' Balanced Salt Solution X10 without Calcium and Magnesium	500 mL
02-018-1A	Hanks' Balanced Salt Solution without Calcium, Magnesium and Phenol Red	500 mL

Dulbecco's Phosphate-Buffered Saline (DPBS)		
02-020-1A	Dulbecco's Phosphate Buffered Saline with Calcium and Magnesium	500 mL
02-020-1B	Dulbecco's Phosphate Buffered Saline with Calcium and Magnesium	100 mL
02-023-1A	Dulbecco's Phosphate Buffered Saline without Calcium and Magnesium	500 mL
02-023-1B	Dulbecco's Phosphate Buffered Saline without Calcium and Magnesium	100 mL
02-023-5A	Dulbecco's Phosphate Buffered Saline X10 without Calcium and Magnesium	500 mL
03-015-1B	EDTA Solution 0.05%	100 mL
HEPES Buffer Solution		
03-025-1B	HEPES Buffer 1M	100 mL
03-025-1C	HEPES Buffer 1M	20 mL
Other Buffers & Salt Solutions		
02-045-1A	Alsever's Solution	500 mL
01-862-1B	EDTA 0.5M	100 mL
01-870-1A	TAE Electrophoresis Buffer X50	500 mL
01-871-1A	TBE Electrophoresis Buffer X5, 500 mL	500 mL
01-888-1B	Red Blood Cell (RBC) Lysis Solution	100 mL
Lab Waters		
01-869-1A	Molecular Biology Grade Water (DNase and RNase-free)	500 mL
01-869-1B	Molecular Biology Grade Water (DNase and RNase-free)	100 mL
03-055-1A	Cell Culture Grade Water	500 mL



## Reagents, Supplements & Kits


Order No.	Description	Quantity
20-700-20	EZ-PCR Mycoplasma Detection Kit <i>Only available in North America</i>	1 kit
01-357-1B	Yeastolate Solution X50	100 mL
03-020-1A	L-Glutamine Solution 200 mM	500 mL
03-020-1B	L-Glutamine Solution 200 mM	100 mL
03-020-1C	L-Glutamine Solution 200 mM	20 mL
03-022-1B	L-Alanyl-L-Glutamine (Stable Glutamine) 200 mM	100 mL
03-042-1B	Sodium Pyruvate Solution 100 mM (11.0 mg/mL)	100 mL
03-040-1B	Sodium Bicarbonate Solution (7.5%)	100 mL
12-005-1A	Potassium Chloride Solution 0.075M	500 mL
12-005-1B	Potassium Chloride Solution 0.075M	100 mL

## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)