

SARTORIUS

DMEM, High Glucose

Catalog Number: 01-055-1, 01-055-1A

Contains 4500 mg/L D-glucose without Sodium Pyruvate. This preparation is without L-Glutamine. Add 584.0 mg/L L-Glutamine.

Components	Concentration (mg/L)
INORGANIC SALTS	
CaCl ₂ ·2H ₂ O	264.92
Ferric Nitrate (Fe(NO ₃) ₃ ·9H ₂ O)	0.10
Potassium Chloride (KCl)	400.00
MgSO ₄ ·7H ₂ O	200.00
Sodium Chloride (NaCl)	6400.00
Sodium Bicarbonate (NaHCO ₃)	3700.00
Sodium Phosphate (NaH ₂ PO ₄ ·H ₂ O)	125.00
OTHER COMPONENTS	
D-Glucose	4500.00
Phenol Red	15.00
AMINO ACIDS	
L-Arginine-HCl	84.00
L-Cystine	48.00
Glycine	30.00
L-Histidine HCl-H ₂ O	42.00
L-Isoleucine	105.00
L-Leucine	105.00
L-Lysine-HCl	146.00
L-Methionine	30.00

SARTORIUS

Components	Concentration (mg/L)
L-Phenylalanine	66.00
L-Serine	42.00
L-Threonine	95.00
L-Tryptophan	16.00
L-Tyrosine	72.00
L-Valine	94.00
VITAMINS	
D-Calcium pantothenate	4.00
Choline Chloride	4.00
Folic Acid	4.00
i-Inositol	7.20
Niacinamide	4.00
Pyridoxine HCl	4.00
Riboflavin	0.40
Thiamine HCl	4.00