

RPMI 1640 Medium, HEPES

Catalog Number: 01-106-1A, 01-106-1B

Components	Concentration (mg/L)
INORGANIC SALTS	
Calcium nitrate (Ca(NO ₃) ₂ 4H ₂ O)	100.00
Potassium chloride (KCl)	400.00
Magnesium sulfate (MgSO ₄)	48.84
Sodium chloride (NaCl)	5500.00
Sodium bicarbonate (NaHCO ₃)	2000.00
Sodium Phosphate (Na ₂ HPO ₄)	800.85
OTHER COMPONENTS	
Glucose	2000.00
Glutathione Reduced	1.00
Phenol red	5.00
AMINO ACIDS	
L-Arginine	200.00
L-Asparagine-H ₂ O	50.00
L-Aspartic Acid	20.00
L-Cystine	50.00
L-Glutamic Acid	20.00
L-Glutamine	300.00
Glycine	10.00
L-Histidine	15.00
L-Hydroxyproline	20.00

Components	Concentration (mg/L)
L-Isoleucine	50.00
L-Leucine	50.00
L-Lysine hydrochloride	40.00
L-Methionine	15.00
L-Phenylalanine	15.00
L-Proline	20.00
L-Serine	30.00
L-Threonine	20.00
L-Tryptophan	5.00
L-Tyrosine	20.00
L-Valine	20.00
VITAMINS	
Biotin	0.20
D-Ca Pantothenate	0.25
Choline Chloride	3.00
Folic Acid	1.00
i-Inositol	35.00
Niacinamide	1.00
p-Aminobenzoic Acid (PABA)	1.00
Pyridoxine HCl	1.00
Riboflavin	0.20
Thiamine HCl	1.00
Vitamin B12	0.005

Components

Concentration (mg/L)

Hepes

5958.00