



Formulation for Leibovitz's L-15 Medium ATCC® 30-2008

Inorganic Salts (g/liter)

CaCl ₂ (anhydrous)	0.14000
MgCl ₂ ·6H ₂ O	0.20000
MgSO ₄ (anhydrous)	0.09767
KCl	0.40000
KH ₂ PO ₄ (anhydrous)	0.06000
NaCl	8.00000
Na ₂ HPO ₄ (anhydrous)	0.19000

Amino Acids (g/liter)

L-Alanine	0.22500
L-Arginine (free base)	0.50000
L-Asparagine·H ₂ O	0.25000
L-Cysteine (free base)	0.12000
L-Glutamine	0.30000
Glycine	0.20000
L-Histidine (free base)	0.25000
L-Isoleucine	0.12500
L-Leucine	0.12500
L-Lysine·HCl	0.09370
L-Methionine	0.07500
L-Phenylalanine	0.12500
L-Serine	0.20000
L-Threonine	0.30000
L-Tryptophan	0.02000
L-Tyrosine·2Na·2H ₂ O	0.43000
L-Valine	0.10000

Vitamins (g/liter)

Choline Chloride	0.00100
Riboflavin-5-PO ₄ ·Na·2H ₂ O	0.00010
Folic Acid	0.00100
myo-Inositol	0.00200
Nicotinamide	0.00100
D-Pantothenic Acid (hemicalcium)	0.00100
Pyridoxine·HCl	0.00100
Thiamine·PO ₄ ·Cl·2H ₂ O	0.00100

Other (g/liter)

D-Galactose	0.90000
Phenol Red, Sodium Salt	0.01000
Sodium Pyruvate	0.55000