



# Formulation for McCoy's 5A ABP-Free Medium ATCC® 30-2011

## Inorganic Salts (g/liter)

CaCl <sub>2</sub> (anhydrous)	0.10000
MgSO <sub>4</sub> (anhydrous)	0.09770
KCl	0.40000
NaHCO <sub>3</sub>	2.20000
NaCl	6.46000
NaH <sub>2</sub> PO <sub>4</sub> ·H <sub>2</sub> O	0.58000

## Amino Acids (g/liter)

L-Alanine	0.01336
L-Arginine·HCl	0.04214
L-Asparagine·H <sub>2</sub> O	0.04503
L-Aspartic Acid	0.01997
L-Cysteine·HCl·H <sub>2</sub> O	0.03514
L-Glutamic Acid	0.02210
L-Glutamine	0.21920
Glycine	0.00750
L-Histidine·HCl·H <sub>2</sub> O	0.02096
Hydroxy-L-Proline	0.01970
L-Isoleucine	0.03936
L-Leucine	0.03936
L-Lysine·HCl	0.03654
L-Methionine	0.01492
L-Phenylalanine	0.01652
L-Proline	0.01730
L-Serine	0.02630
L-Threonine	0.01790
L-Tryptophan	0.00310
L-Tyrosine·2Na·2H <sub>2</sub> O	0.02612
L-Valine	0.01760

## Vitamins (g/liter)

Ascorbic Acid	0.00050
p-Amino Benzoic Acid	0.00100
D-Biotin	0.00020
Choline Chloride	0.00500
Folic Acid	0.01000
myo-Inositol	0.03600
Nicotinamide	0.00050
Nicotinic Acid	0.00050
D-Pantothenic Acid (hemicalcium)	0.00020
Pyridoxine·HCl	0.00100
Riboflavin	0.00020
Thiamine·HCl	0.00020
Vitamin B-12	0.00200

## Other (g/liter)

D-Glucose	3.00000
Glutathione (reduced)	0.00050
Phenol Red, Sodium Salt	0.01000