



## Formulation for McCoy's 5A Peptone-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<br>(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 |