

Microbe Acrostic

Instructions: Print the puzzle or fill out in Adobe Acrobat. Enter the answers to the clues, one letter per square. Then enter each letter in its corresponding spot in the grid. When you are done, the grid will spell out a quote. The first letters of the clue answers will spell the author and the title of the work.

1 E	2 V	3 U	4 C	5 J	6 I		7 U	8 S	9 K		10 S	11 L	12 J	13 A	14 V	15 Q		16 B	17 T		18 J	19 W	20 J	21 D		22 S	23 V	24 T		25 V	26 E	27 A
28 Q	29 O	30 V		31 I	32 J		33 P		34 H	35 R	36 E	37 S		38 I	39 P		40 J	41 D	42 U		43 I	44 C	45 F		46 X	47 Q		48 M	49 A	50 D	51 U	
52 N	53 L	54 U	55 D	56 S		57 X	58 J	59 V	60 F	61 A		62 Q	63 H	64 V		65 X	66 U	67 K	68 H	69 O	70 N	71 C		72 L	73 T	74 X		75 P	76 F	77 R	78 X	79 U
	80 I	81 Q	82 O	83 D	84 G	85 H		86 U	87 M		88 F	89 U	90 V	91 U	92 G	93 E		94 N	95 K	96 G	97 R		98 I	99 F	100 N		101 D	102 U	103 C	104 U		105 H
106 B	107 G		108 O	109 R	110 C	111 M		112 P	113 V	114 B	115 X		116 V	117 D	118 F	119 N	120 D	121 S	122 K	123 P	124 W		125 T	126 H	127 N		128 J	129 H	130 Q	131 T	132 P	

Tips for filling out with Adobe Reader/Acrobat: The cursor will advance to the next clue space after typing in the clue spaces or the grid spaces; hit shift+tab to backspace.

A. Name of a bacterial cell wall stain, such as CF[®] Dye WGA

61 13 49 27

B. Number of colonies to pick to isolate a clone

16 106 114

C. Common wavelength range of spectrophotometers (hyphen)

44 103 110 4 71

D. Workspace for many scientists (2 wds)

101 55 83 120 21 117 50 41

E. Old-fashioned contest to settle a point of honor

1 26 93 36

F. Particle of light

88 118 76 45 99 60

G. Toxic heavy metal

92 84 96 107

H. One of the three domains of life

63 85 34 126 129 68 105

I. What you're using to solve this puzzle

98 31 6 43 38 80

J. Wastewater; common sample for vPCR testing

128 20 32 18 12 58 5 40

K. Skilled lab worker

122 9 67 95

L. Not in

11 53 72

M. Toll or charge

87 48 111

N. Defrosted

94 52 119 100 127 70

O. Iron-containing porphyrin compound

108 29 82 69

P. Sweet-smelling hydrocarbon

33 75 132 112 39 123

Q. Solution that resists pH change

28 81 47 62 15 130

R. Metric prefix for a tiny number

109 77 97 35

S. Fungal polysaccharide; Calcofluor staining target

10 8 37 22 121 56

T. Intellectual property concern

125 73 24 17 131

U. Tough cookie of the microbial world

42 66 7 3 89 54 86 91 51 102 79 104

V. Possible cold cause

64 23 116 25 90 113 59 30 2 14

W. One of three elements of the psyche named by Freud

19 124

X. Organism that grows in the presence of oxygen

78 65 115 46 57 74