

Agilent BioTek Synergy H1 Shake Settings Optimized for Volume and Plate Type

The Agilent BioTek Synergy H1 multimode reader offers a broad range of shaking speeds and amplitudes. If the wells of a microplate are almost full, shaking can result in spillage inside the instrument. The following table is designed to help prevent this. Select your microplate type and assay volume, and you will get the acceptable shake amplitude on your instrument.

Sample Volume (mL)	Linear (mm)	Orbital Slow (mm)	Orbital Fast (mm)	Double Orbital Slow (mm)	Double Orbital Fast (mm)
6-Well Microplate					
0 to 3	1 to 6	1 to 6	1 to 6	1 to 6	1 to 6
3 to 4	1 to 6	1 to 6	1 to 6	1 to 6	1
4 to 5	1	1 to 6	1 to 6	1 to 6	1
5 to 7	1	1 to 6	1 to 3	1	No
7 to 8	1	No	1 to 3	No	No
9	1	No	No	No	No
12-Well Microplate					
0 to 1.5	1 to 6	1 to 6	1 to 6	1 to 6	1 to 6
1.5 to 2.5	1 to 2	1 to 6	1 to 6	1 to 6	1
2.5 to 3	1 to 2	1 to 6	No	No	1
3 to 4	No	1 to 6	No	No	No
24-Well Microplate					
0 to 0.75	1 to 6	1 to 6	1 to 6	1 to 6	1 to 6
0.75 to 1	1	1 to 6	1 to 6	1 to 6	1 to 6
1 to 1.5	1	1 to 6	1 to 6	1 to 6	No
1.5 to 2	1	1 to 6	No	1 to 6	No
48-Well Microplate					
0 to 0.5	1 to 6	1 to 6	1 to 6	1 to 6	1 to 6
0.5 to 1	1	1 to 6	1 to 6	1 to 6	1 to 6
1 to 1.3	No	1 to 6	1 to 6	1 to 6	No
96-Well Microplate					
0 to 0.25	1 to 6	1 to 6	1 to 6	1 to 6	1 to 6

www.agilent.com/lifesciences/biotek

DE44335.0669675926

This information is subject to change without notice.