

Celigo® Plate Profiles

The following table lists the Celigo plate profiles that are provided with software release v5.2.

Plate Name	Manufacturer	Well Type	Compatibles
6-Well BD Falcon™ 353046 Plate	Corning	Clear	353224, 353934, 353846, 351146, 353502
6-Well Corning™ 3516 Plate	Corning	Clear	3471, 3506, 3335
6-Well CytoOne® CC7682-7506 Plate	CytoOne	Clear	-
6-Well Greiner™ 657160 Plate	Greiner	Clear	657185, 657165
6-Well Nunc™ 140675 Plate	Thermo	Clear	-
12-Well BD Falcon™ 353043 Plate	Corning	Clear	353224, 351143, 353503
12-Well Corning™ 3513 Plate	Corning	Clear	3336, 3512
12-Well CytoOne® CC7682-7512 Plate	CytoOne	Clear	-
24-Well BD Falcon 353047 Plate	Corning	Clear	353226, 353935, 353847, 351147, 358115, 354723, 356723, 354775, 356775, 353504
24-Well Corning 3524 Plate	Corning	Clear	3337, 3526, 3527, 3573
24-Well CytoOne CC7682-7524 Plate	CytoOne	Clear	-
24-Well Greiner 662160 Plate	Greiner	Clear	662102, 622165
24-Well Revvity Visiplate 1450606 Plate	Revvity	Black	-
24-Well Seahorse XF24 Plate	Seahorse Biosciences	Clear	-
48-Well Corning 3548 Plate	Corning	Clear	-
48-Well Greiner 677180 Plate	Greiner	Clear	677102, 677165
96-Well BD Falcon 353219 Plate	Corning	Black, White	353377
96-Well BD Falcon 354640 Plate	Corning	Black, White	354650, 356650, 354651, 356651, 356701, 356693, 354649, 356649, 356640, 356700, 356692, 356717
96-Well BD Falcon 356717 Plate	Corning	Black	354717
96-Well BD Falcon 351177 U-Bottom Plate	Corning	Clear, Round Bottom ULA	-

Plate Name	Manufacturer	Well Type	Compatibles
96-Well BD Falcon 353072 Plate	Corning	Clear	351172, 353075, 354407, 354429, 353916, 353936, 354409, 354410, 354461, 354516, 354596, 354607, 354657, 354670, 354689, 356461, 356516, 356689, 356690, 356698
96-Well BD Falcon 353219 Plate	Corning	Black, White	353377
96-Well BD Falcon 354640 Plate	Corning	Black, White	354650, 354651, 356640, 356650, 356651, 354649, 356649, 356692, 356693, 356700, 356701, 356717
96-Well BD Falcon 356717 Plate	Corning	Black	354717
96-Well BD Falcon 351177 U-Bottom Plate	Corning	Clear, Round Bottom ULA	-
96-Well Corning 3596 Plate	Corning	Clear	3300, 3474, 3595, 3598, 3599, 3585, 3595, 3628, 3841
96-Well Corning 3603 Plate	Corning	Black, White	3340, 3601, 3604, 3610, 3631, 3632, 3635, 3651, 3843, 3842, 3903, 3904
96-Well Corning 3696 Plate	Corning	Half Area, Black	3686, 3688, 3690, 3693, 3694, 3695, 3696, 3697
96-Well Corning 7007 U-Bottom Plate	Corning	Clear, Round Bottom ULA	-
96-Well Greiner 655090 Plate	Greiner	Black, White	655087, 655094, 655097, 655098, 655936, 655944, 655946, 655948, 655956
96-Well Greiner 655087 Plate	Greiner	Black	655088
96-Well Greiner 655161 Plate	Greiner	Clear	655101, 655192
96-Well Greiner 655180 Plate	Greiner	Clear, chimney	655182, 655185, 655930, 655940, 655950
96-Well Greiner 675986 Plate	Greiner	Half Area, Black	67509x
96-Well Greiner 650185 U-Bottom Plate	Greiner	Clear, Round Bottom ULA	-
96-well ULA-96U Plate	-	Clear, Round Bottom ULA	-
96-Well Nunc 167008 Plate	Thermo	Clear	-
96-Well Revvity Viewplate 6005225 Plate	Revvity	Black	-
96-Well Revvity Isoplate 6005050 Plate	Revvity	Black	-
96-Well Seahorse FX96 Plate	Seahorse Bioscience	Black	-
384-well ULA-384U Plate	-	Clear, Round Bottom ULA	-
384-Well BD Falcon 353962 Plate	Corning	Clear	-
384-Well Corning 3542 Plate	Corning	Low volume, Black	3540

Plate Name	Manufacturer	Well Type	Compatibles
384-Well Corning 3680 Plate	Corning	Clear	3640, 3844, 3700, 3701, 3702, 3844
384-Well Corning 3712 Plate	Corning	Black, White	3653, 3655, 3846, 3845, 3706, 3707, 3711, 3683, 3845, 3846
384-Well Corning 3827 Plate	Corning	Low attach	-
384-Well Greiner 781182 Plate	Greiner	Clear	781185, 781186, 781061, 781940, 781930, 781950
384-Well Greiner 781091 Plate	Greiner	Black	781098, 781095, 781094, 781944, 781090, 781096, 781097, 781946, 781948, 781936, 781956
1536-Well Corning 3838 Plate	Corning	Black, White	3833, 3836, 3893
1536-Well BD Falcon 356771 Plate	Corning	Black	-
1536-Well Greiner 789866 Plate	Greiner	Black	789896
1-Well Nunc Omnitray	Thermo Fisher	Clear	242811
T25 Greiner 690175 Flask	Greiner	Clear	-
T25 Greiner 690175 Flask - Single View	Greiner	Clear	-
T25 BD Falcon 353014 Flask	Corning	Clear	-
T25 BD Falcon 353014 Flask - Single View	Corning	Clear	-
T75 BD Falcon 353136 Flask	Corning	Clear	-
T75 BD Falcon 353136 Flask - Single View	Corning	Clear	-
10cm Dish BD Falcon 353003 Dish	Corning	Clear	353803
1-Slide Holder (2/3 cover slip)	-	Clear	-
1-Slide Holder (square cover slip)	-	Clear	-
4-Slide Holder (2/3 cover slip)	-	Clear	-
4-Slide Holder (square cover slip)	-	Clear	-

For research use only. Not approved for diagnostic or therapeutic use.

