

Reflects vessels supported by software version 2023A

Updated September 2023

Directions: Use CTRL-F to search by catalog number

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Agilent Seahorse	Plate	24	N/A	100777-004	24-well Seahorse XF24 V7-PS	Microplates	Standard, ATP	Yes	Yes	Yes
Agilent Seahorse	Plate	24	N/A	101037-004	24-well Seahorse XF24 V7-PET	Microplates	Standard, ATP	Yes	Yes	Yes
Aurora	Plate	384	N/A	ABB1-10100A, ABB1-11100A, ABB1-11101A, ABC1-10100A, ABC1-11100A, ABC1-11101A, ACB1-10100A, ACB1-11100A, ACB1-11101A, ACC1-10100A, ACC1-11100A, ACC1-11101A, AWB1-10100A, AWB1-11100A, AWB1-11101A, AWC1-10100A, AWC1-11100A, AWC1-11101A	384-well Aurora IQ-EB	Microplates	Standard, ATP	No	Yes	Yes
Aurora	Plate	384	N/A	ABE2-10100A, ABE2-11100A, ABE2-11101A	384-well Aurora IQ-EB Ultra Low Base	Microplates	Standard, ATP	No	Yes	Yes
BRAND	Plate	96	N/A	7816 60, 7819 00, 7819 60	96-well BRANDplate	Microplates	Spheroid	Yes	Yes	No
Cellvis	Plate	12	N/A	P12-1.5H-N	12-well Cellvis	Microplates	Standard, ATP	Yes	Yes	Yes
Cellvis	Plate	96	N/A	P96-1.5H-N	96-well Cellvis	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	6	N/A	353046, 353224, 353846, 353934, 354400, 354402, 354404, 354413, 354417, 354428, 354431, 354432, 354510, 354515, 354595, 354603, 354652, 354658, 356400, 356413, 356515, 356652	6-well Corning Falcon	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	6	N/A	3335, 3471, 3506, 3516	6-well Corning	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	12	N/A	353043, 353225, 354470, 354500, 354501, 354502, 354503, 356470	12-well Corning Falcon	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	12	N/A	3336, 3512, 3513	12-well Corning	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	24	N/A	3337, 3473, 3524, 3526, 3527	24-well Corning	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	24	N/A	353047, 353226, 353847, 353935, 354408, 354411, 354412, 354414, 354433, 354605, 354619, 354635, 354659, 356408, 356414	24-well Corning Falcon	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Plate	24	N/A	3337, 3524, 3526, 3527	24-well Corning	Microplates	OrganoidQC	Yes	No	No

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Corning	Plate	48	N/A	3338, 3548	48-well Corning	Microplates	Standard, ATP	Yes	Yes	Yes
Corning	Plate	48	N/A	353078, 353230, 354505, 354506, 354507, 354508, 354509, 356505, 356509	48-well Corning Falcon	Microplates	Standard, ATP	Yes	Yes	Yes
Corning	Plate	48	N/A	3338, 3548	48-well Corning	Microplates	OrganoidQC	Yes	No	No
Corning	Plate	96	N/A	356259	96-well Corning Matrigel Matrix 3D Plate	Microplates	Multi Spheroid	No	Yes	No
Corning	Plate	96	N/A	3300, 3474, 3585, 3595, 3596, 3598, 3599, 3628, 3997	96-well Corning	Microplates	Organoid Assay, Embedded Multi Spheroid	Yes	No	No
Corning	Plate	96	N/A	4515, 4520	96-well Corning Black ULA Spheroid	Microplates	Spheroid	No	Yes	No
Corning	Plate	96	N/A	353219, 354640, 354649, 354650, 354651, 356640, 356649, 356650, 356651, 356692, 356693, 356700	96-well Corning Falcon Blk/Clr Imaging	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	96	N/A	356519	96-well Corning Falcon Optilux	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	96	N/A	353072, 353075, 353872, 353916, 353936, 354407, 354409, 354410, 354429, 354461, 354516, 354596, 354607, 354657, 354670, 354689, 356407, 356461, 356516, 356689, 356690, 356698	96-well Corning Falcon (All Clr)	Microplates	Standard, ATP, Whole Well, Dilution Cloning, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	96	N/A	3595	96-well Corning	Microplates	Multi Spheroid	No	Yes	No
Corning	Plate	96	N/A	7007	96-well Corning Round Bottom ULA	Microplates	Standard, Spheroid	Yes	Yes	No
Corning	Plate	96	N/A	3300, 3474, 3585, 3595, 3596, 3598, 3599, 3628, 3997	96-well Corning	Microplates	Standard, ATP, Whole Well, Dilution Cloning, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	96	N/A	4580, 4582, 4584, 4586	96-well Corning Half Area Glass Bottom	Microplates	Standard, ATP	No	Yes	Yes
Corning	Plate	96	N/A	3340, 3603, 3610, 3841, 3842, 3843, 3903, 3904	96-well Corning (Blk/Wht)	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Corning	Plate	384	N/A	353962, 353963, 354660, 354663, 354664, 354667, 356660, 356663, 356664, 356667, 356697, 356705	384-well Corning Falcon (Blk/Wht)	Microplates	Standard, ATP	No	Yes	Yes
Corning	Plate	384	N/A	353961, 354662, 354666, 356662, 356666	384-well Corning Falcon (All Clr)	Microplates	Standard, ATP	No	Yes	Yes
Corning	Plate	384	N/A	3764, 4588	384-well Corning (Blk/Wht)	Microplates	Standard, Whole Well, Dilution Cloning	Yes	Yes	Yes
Corning	Plate	384	N/A	3701	384-well Corning (All Clr)	Microplates	Standard, ATP	No	Yes	Yes
Corning	Plate	384	N/A	3985	384-well Corning (Opt Imaging)	Microplates	Standard, ATP, Whole Well, Dilution Cloning	Yes	Yes	Yes
Corning	Flask	N/A	25	353082, 353109, 354484, 354532, 354533, 354534, 354536, 356484, 356536	Corning Falcon 25 B	Tray 12	Standard, ATP	Yes	Yes	Yes

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Corning	Flask	N/A	25	353014, 353108, 353808, 353813	Corning Falcon 25 A	Tray 20	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	25	3055, 3056	Corning 25 A	Tray 1	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	25	430168, 430372, 430639	Corning 25 B	Tray 12	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	75	3275, 3276	Corning 75 B	Tray 14	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	75	3290	Corning 75 A	Tray 15	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	75	353024, 353110, 353810, 353824	Corning Falcon 75 B	Tray 14	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	75	3290, 3814, 430641U, 430720U, 430725U, 431464U	Corning 75 U-shaped	Tray 30	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	75	353133, 353135, 353136, 354485, 354488, 354521, 354523, 354537, 356485, 356488, 356537	Corning Falcon 75 A	Tray 15	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	93	3069, 3070	RoboFlask	Microplates	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	100	3073, 3816	Corning 100 Low Profile	Tray 22	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	150	3291, 430823, 430824, 430825	Corning 150 U-shaped	Tray 32	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	150	3291, 430823, 430824, 430825	Corning 150	Tray 6	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	150	354486, 354538, 354646, 355000, 355001, 356486, 356538	Corning Falcon 150	Tray 5	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	162	3150, 3151	Corning 162	Tray 4	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	175	3292, 431079, 431080, 431085, 431306	Corning 175	Tray 5	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	175	3292, 431079, 431080, 431085	Corning 175 U-shaped	Tray 5	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	175	353028, 353112, 354487, 354526, 354528, 356487	Corning Falcon 175	Tray 5	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	225	3000, 3001	Corning 225 A	Tray 7	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	225	353138, 353139	Corning Falcon 225	Tray 10	Standard, ATP	Yes	Yes	Yes
Corning	Flask	N/A	225	431081, 431082	Corning 225 B	Tray 10	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	8	430165	Corning 35mm Dish with Holder	Tray 25	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Dish	N/A	9	430165	Corning 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	9	353001	Falcon 35mm Dish with Holder	Tray 25	Standard, ATP, Whole Well	Yes	Yes	Yes
Corning	Dish	N/A	10	353001	Corning Falcon 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	21	353002	Corning Falcon 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	22	353004	Corning Falcon Easy Grip 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	23	430166	Corning 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	63	353003, 353803, 354450, 354451, 354452, 354453, 354455, 354469, 354600, 354653, 356450, 356469, 356653	Corning Falcon 100mm Dish	Tray 9	Standard, ATP	Yes	Yes	Yes

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Corning	Dish	N/A	63	3262, 3296, 430167, 430293	Corning 100mm Dish	Tray 9	Standard, ATP	Yes	Yes	Yes
Corning	Dish	N/A	153	430599	Corning 150mm Dish	Tray 11	Standard, ATP	Yes	Yes	Yes
CytoOne	Plate	6	N/A	CC7682-7506	6-well CytoOne	Microplates	Standard, ATP	Yes	Yes	Yes
CytoOne	Plate	12	N/A	CC7682-7512	12-well CytoOne	Microplates	Standard, ATP	Yes	Yes	Yes
CytoOne	Plate	24	N/A	CC7682-7524	24-well CytoOne	Microplates	Standard, ATP	Yes	Yes	Yes
CytoOne	Plate	48	N/A	CC7682-7548	48-well CytoOne	Microplates	Standard, ATP	Yes	Yes	Yes
CytoOne	Plate	96	N/A	CC7672-7596, CC7682-7596	96-well CytoOne	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Eppendorf	Plate	6	N/A	0030 720.016, 0030 720.113, 0030 720.121	6-well Eppendorf	Microplates	Standard, ATP	Yes	Yes	Yes
Eppendorf	Plate	12	N/A	0030 721.012, 0030 721.110	12-well Eppendorf	Microplates	Standard, ATP	Yes	Yes	Yes
Eppendorf	Plate	24	N/A	0030 722.019, 0030 722.116	24-well Eppendorf	Microplates	Standard, ATP	Yes	Yes	Yes
Eppendorf	Plate	48	N/A	0030 723.015, 0030 723.112	48-well Eppendorf	Microplates	Standard, ATP	Yes	Yes	Yes
Eppendorf	Plate	96	N/A	0030 730.011, 0030 730.119, 0030 730.127	96-well Eppendorf	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Eppendorf	Flask	N/A	25	0030 710.010, 0030 710.029, 0030 710.118, 0030 710.126	Eppendorf 25	Tray 26	Standard, ATP	Yes	Yes	Yes
Eppendorf	Flask	N/A	75	0030 711.017, 0030 711.025, 0030 711.114, 0030 711.122	Eppendorf 75	Tray 28	Standard, ATP	Yes	Yes	Yes
Eppendorf	Flask	N/A	175	0030 712.110, 0030 712.129	Eppendorf 175	Tray 29	Standard	Yes	Yes	Yes
Essen BioScience	Slide	1	N/A	95004380	FLR Calibration Slide	MicroSlides	Standard, ATP	Yes	Yes	Yes
Essen BioScience	Plate	96	N/A	4582, 4599, 4648	Incucyte® 96-Well Clear-view Plate	Microplates	Chemotaxis (Top/Bot)	No	Yes	No
Essen BioScience	Plate	96	N/A	4379	Incucyte® 96-Well Image-lock Plate	Microplates	Scratch Wound, Standard (with lock)	Yes	Yes	Yes
Essen BioScience	Plate	96	N/A	4600, 4601	Incucyte® 96-Well Clear-view Reservoir	Microplates	Standard, ATP	Yes	Yes	Yes
Greiner	Plate	6	N/A	657160	6-well Greiner	Microplates	Standard, ATP	Yes	Yes	Yes
Greiner	Plate	24	N/A	662102, 662160, 662165	24-well Greiner	Microplates	Standard, ATP	Yes	Yes	Yes
Greiner	Plate	96	N/A	655087, 655088, 655090, 655098, 655976	96-well Greiner micro-clear	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Greiner	Plate	96	N/A	655160, 655162, 655180, 655182, 655980, 655982	96-well Greiner	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Greiner	Plate	384	N/A	781090, 781091, 781092, 781093, 781094, 781095, 781096, 781097, 781098, 781801, 781936, 781944, 781946, 781948, 781956	384-well Greiner micro-clear	Microplates	Standard, ATP	No	Yes	Yes
Greiner	Plate	384	N/A	781165, 781182	384-well Greiner (All Clr)	Microplates	Standard, ATP	No	Yes	Yes
Greiner	Flask	N/A	25	690160	Greiner 25	Tray 20	Standard, ATP	Yes	Yes	Yes
Greiner	Flask	N/A	75	658170, 658175, 658195, 658940, 658950	Greiner 75	Tray 14	Standard, ATP	Yes	Yes	Yes
Greiner	Flask	N/A	182	660160, 660175, 661160, 661175, 661195, 661940, 661950	Greiner 182	Tray 6	Standard, ATP	Yes	Yes	Yes

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Greiner	Dish	N/A	8	627160	Greiner 35mm Dish with Holder	Tray 25	Standard, ATP, Whole Well	Yes	Yes	Yes
Greiner	Dish	N/A	9	627160	Greiner 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Greiner	Dish	N/A	23	628160	Greiner 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Greiner	Dish	N/A	64	664160, 664940, 664950	Greiner 100mm Dish	Tray 9	Standard, ATP	Yes	Yes	Yes
Greiner	Dish	N/A	153	639160	Greiner 150mm Dish	Tray 11	Standard, ATP	Yes	Yes	Yes
Greiner/GNF	Flask	N/A	84	779160	CELLSTAR AutoFlask	Microplates	Standard, ATP	Yes	Yes	Yes
IBIDI	Slide	8	N/A	80821, 80822, 80823, 80824, 80825, 80826	8-well IBIDI MicroSlide	Tray 18	Standard, ATP	Yes	Yes	Yes
IBIDI	Slide	36	N/A	80601, 80602, 80604, 80606	IBIDI MicroSlide VI	Tray 18	Standard, ATP	Yes	Yes	Yes
IBIDI	Flask	N/A	2	80171, 80172, 80176	IBIDI MicroSlide I	Tray 18	Standard, ATP	Yes	Yes	Yes
IBIDI	Flask	N/A	3	80121, 80126	IBIDI MicroSlide Y	Tray 18	Standard, ATP	No	Yes	Yes
IBIDI	Dish	N/A	3	81151, 81156	IBIDI 35mm Dish	Tray 21	Standard, ATP	No	Yes	Yes
Iwaki	Plate	6	N/A	3810-006, 4810-010, 4810-020	6-well Iwaki	Microplates	Standard, ATP	Yes	Yes	Yes
Iwaki	Plate	12	N/A	3815-012	12-well Iwaki	Microplates	Standard	Yes	Yes	Yes
Iwaki	Plate	24	N/A	5826-024	24-well Iwaki Glass Bot	Microplates	Standard	Yes	Yes	Yes
Iwaki	Plate	24	N/A	1820-024, 3820-024	24-well Iwaki	Microplates	Standard	Yes	Yes	Yes
Iwaki	Plate	48	N/A	1830-048, 3830-048	48-well Iwaki	Microplates	Standard	Yes	Yes	Yes
Iwaki	Plate	96	N/A	3860-096, 3861-096	96-well Iwaki	Microplates	Standard, Neuronal Activity	Yes	Yes	Yes
Iwaki	Plate	384	N/A	3721-384	384-well Iwaki	Microplates	Standard	No	Yes	Yes
Iwaki	Flask	N/A	25	3100-025	Iwaki 25	Tray 12	Standard, ATP	Yes	Yes	Yes
Iwaki	Flask	N/A	75	3110-075	Iwaki 75	Tray 15	Standard, ATP	Yes	Yes	Yes
Iwaki	Dish	N/A	11	1000-035, 3000-035	Iwaki 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Iwaki	Dish	N/A	23	1010-060, 3010-060	Iwaki 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Matrix	Plate	384	N/A	4313, 4336	384-well Matrix Screen-Mates (AllClr)	Microplates	Standard, ATP	No	Yes	Yes
Matrix/Greiner	Plate	384	N/A	4329, 4330, 4331, 4332	384-well Matrix Screen-Mates (Blk/Wht)	Microplates	Standard, ATP	No	Yes	Yes
MatTek	Plate	6	N/A	P06G-1.5-20-F	6-well MatTek	Microplates	Standard, ATP	Yes	Yes	Yes
MatTek	Plate	12	N/A	P12G-1.5-14-F	12-well MatTek	Microplates	Standard, ATP	Yes	Yes	Yes
MatTek	Plate	24	N/A	P24G-1.5-13-F	24-well MatTek	Microplates	Standard, ATP	Yes	Yes	Yes
MatTek	Dish	N/A	9	P35G-1.0-14-C	MatTek 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Nanofiber Solutions	Plate	96	N/A	9601, 9602	96-well Nanofiber Solutions	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Nunc	Slide	8	N/A	154534	8-well Nunc Lab-Tek Slide	MicroSlides	Standard, ATP	Yes	Yes	Yes
Nunc	Plate	6	N/A	140675, 140685	6-well Nunc	Microplates	Standard, ATP	Yes	Yes	Yes
Nunc	Plate	8	N/A	167064	8-well Nunc	Microplates	Standard, ATP	Yes	Yes	Yes
Nunc	Plate	12	N/A	150628	12-well Nunc	Microplates	Standard, ATP	Yes	Yes	Yes
Nunc	Plate	24	N/A	142475, 142485	24-well Nunc	Microplates	Standard, ATP	Yes	Yes	Yes
Nunc	Plate	96	N/A	156545, 161093, 167008, 168055	96-well Nunc	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Nunc	Plate	96	N/A	167311, 167314, 267312, 267313	96-well Nunc Edge	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Nunc	Plate	96	N/A	152028, 152036, 152037, 152040, 165305, 165306	96-well Nunc opt bottom (Blk/Wht)	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
Nunc	Plate	384	N/A	142761, 142762, 152029, 152041	384-well Nunc opt bottom (Blk/Wht)	Microplates	Standard, ATP	No	Yes	Yes
Nunc	Plate	384	N/A	164555, 164688	384-well Nunc	Microplates	Standard, ATP	No	Yes	Yes
Nunc	Flask	N/A	25	136196, 163371	Nunc 25	Tray 20	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	25	132703, 132706, 156340, 156367	Nunc EasYFlask 25	Tray 13	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	75	132704, 132707, 156472, 156499	Nunc EasYFlask 75	Tray 17	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	75	153732, 178905	Nunc 75	Tray 14	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	175	132705, 132708, 159910, 159920	Nunc EasYFlask 175	Tray 5	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	175	156502, 178883, 178983	Nunc 175	Tray 5	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	175	132867, 132913, 132920	Nunc TripleFlask	Tray 5	Standard, ATP	Yes	Yes	Yes
Nunc	Flask	N/A	225	159933, 159934	Nunc 225	Tray 10	Standard, ATP	Yes	Yes	Yes
Nunc	Dish	N/A	9	150318, 153066	Nunc 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Nunc	Dish	N/A	18	174888	Nunc 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Nunc	Dish	N/A	63	150350, 150679, 172958	Nunc 100mm Dish	Tray 9	Standard, ATP	Yes	Yes	Yes
Nunc	Dish	N/A	153	168381	Nunc 150mm Dish	Tray 11	Standard, ATP	Yes	Yes	Yes
Nunclon	Plate	48	N/A	150687, 152640	48-well Nunc	Microplates	Standard, ATP	Yes	Yes	Yes
ORGANOGENIX	Plate	96	N/A	NCP-HH-96-10, NCP-HH-96-2, NCP-HS-96-10, NCP-HS-96-2, NCP-LH-96-10, NCP-LH-96-2, NCP-LS-96-10, NCP-LS-96-2, NCP-LSH-96-2	96-well ORGANOGENIX	Microplates	Standard, Neuronal Activity	Yes	No	No
PerkinElmer	Plate	96	N/A	6055300, 6055302, 6055308, 6055500, 6055508	96-well PerkinElmer CellCarrier-96 Ultra	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
PerkinElmer	Plate	96	N/A	6055300, 6055302, 6055308	96-well PerkinElmer CellCarrier-96 Ultra	Microplates	Multi Spheroid	No	Yes	No
PerkinElmer	Plate	96	N/A	6005181, 6005182, 6005225	96-well PerkinElmer ViewPlate	Microplates	Standard, ATP, Neuronal Activity	No	Yes	No
PerkinElmer	Plate	96	N/A	6005450, 6005458, 6005550, 6005558, 6005920, 6005928	96-well PerkinElmer CellCarrier	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes
PerkinElmer	Plate	384	N/A	6007480, 6007490	384-well PerkinElmer ViewPlate	Microplates	Standard, ATP	No	Yes	Yes
PerkinElmer	Plate	384	N/A	6057300, 6057302, 6057308	384-well PerkinElmer CellCarrier-384 Ultra	Microplates	Standard, ATP	No	Yes	Yes
Porvair	Plate	384	N/A	311003, 311503, 312003, 312030, 312503	384-well Porvair Krystal	Microplates	Standard, ATP	No	Yes	Yes
Sarstedt	Plate	6	N/A	83.3920, 83.3920.300, 83.3920.500	6-well Sarstedt	Microplates	Standard, ATP	Yes	Yes	Yes
Sarstedt	Plate	12	N/A	83.3921, 83.3921.300, 83.3921.500	12-well Sarstedt	Microplates	Standard, ATP	Yes	Yes	Yes
Sarstedt	Plate	24	N/A	83.3922, 83.3922.300, 83.3922.500	24-well Sarstedt	Microplates	Standard, ATP	Yes	Yes	Yes
Sarstedt	Plate	96	N/A	83.3924, 83.3924.300, 83.3924.500	96-well Sarstedt	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes

Manufacturer	Vessel Type	Wells	Area (cm ²)	Catalog #s	Vessel Name	Tray Name	Scan Types	4x	10x	20x
Sarstedt	Flask	N/A	25	83.3910, 83.3910.002, 83.3910.300, 83.3910.302, 83.3910.500, 83.3910.502	Sarstedt 25	Tray 13	Standard	Yes	Yes	Yes
Sarstedt	Flask	N/A	75	83.3911, 83.3911.002, 83.3911.300, 83.3911.302, 83.3911.500, 83.3911.502	Sarstedt 75	Tray 17	Standard	Yes	Yes	Yes
Sarstedt	Dish	N/A	9	83.3900, 83.3900.300, 83.3900.500	Sarstedt 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
Sarstedt	Dish	N/A	23	83.3901, 83.3901.300, 83.3901.500	Sarstedt 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
Sarstedt	Dish	N/A	58	83.3902, 83.3902.300, 83.3902.500	Sarstedt 100mm Dish	Tray 9	Standard, ATP	Yes	Yes	Yes
Sartorius	Plate	96	N/A	BA-04855, BA-04856, BA-04857	Incucyte® 96-Well Imago-lock Plate	Microplates	Scratch Wound, Standard (with lock)	Yes	Yes	Yes
S-BIO	Plate	6	N/A	MS-80060	6-well S-BIO	Microplates	Standard, ATP	Yes	Yes	Yes
S-BIO	Plate	12	N/A	MS-80120	12-well S-BIO	Microplates	Standard, ATP	Yes	Yes	Yes
S-BIO	Plate	24	N/A	MS-80240	24-well S-BIO	Microplates	Standard, ATP	Yes	Yes	Yes
S-BIO	Plate	48	N/A	MS-80480	48-well S-BIO	Microplates	Standard, ATP	Yes	Yes	Yes
S-BIO	Plate	96	N/A	MS-9096UZ	96-well S-BIO PrimeSurface	Microplates	Standard, Spheroid	Yes	Yes	No
S-BIO	Plate	96	N/A	MS-3096F, MS-8096F, MS-8196F5	96-well S-BIO	Microplates	Standard, ATP	Yes	Yes	Yes
S-BIO	Plate	384	N/A	MS-9384UZ	384-well S-BIO PrimeSurface	Microplates	Spheroid	No	Yes	No
SPL	Plate	6	N/A	30006, 31006, 32006	6-well SPL	Microplates	Standard, ATP	Yes	Yes	Yes
SPL	Plate	12	N/A	30012, 31012, 32012	12-well SPL	Microplates	Standard, ATP	Yes	Yes	Yes
Thermo Scientific	Plate	12	N/A	130185	12-well Thermo BioLite	Microplates	Standard, ATP	Yes	Yes	Yes
Thermo Scientific	Flask	N/A	75	130190, 130193	Thermo BioLite 75	Tray 17	Standard, ATP	Yes	Yes	Yes
TPP	Plate	6	N/A	92006, 92406	6-well TPP	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
TPP	Plate	12	N/A	92012, 92412	12-well TPP	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
TPP	Plate	24	N/A	92024, 92424	24-well TPP	Microplates	Standard, ATP, Whole Well	Yes	Yes	Yes
TPP	Plate	96	N/A	92096, 92696	96-well TPP	Microplates	Standard, ATP, Whole Well, Neuronal Activity	Yes	Yes	Yes
TPP	Flask	N/A	25	90025, 90026	TPP 25	Tray 13	Standard, ATP	Yes	Yes	Yes
TPP	Flask	N/A	75	90075, 90076	TPP 75	Tray 17	Standard, ATP	Yes	Yes	Yes
TPP	Flask	N/A	150	90150, 90151	TPP 150	Tray 5	Standard, ATP	Yes	Yes	Yes
TPP	Dish	N/A	9	93040	TPP 35mm Dish	Tray 21	Standard, ATP	Yes	Yes	Yes
TPP	Dish	N/A	23	93060	TPP 60mm Dish	Tray 19	Standard, ATP	Yes	Yes	Yes
TPP	Dish	N/A	153	93150	TPP 150mm Dish	Tray 11	Standard, ATP	Yes	Yes	Yes
WPI	Dish	N/A	4	FD35-100, FD35PDL-100	WPI 35mm Fluorodish	Tray 21	Standard, ATP	Yes	Yes	Yes
zell-kontakt	Plate	96	N/A	5241-20	96-well zell-kontakt Imaging Plate CG	Microplates	Standard, ATP, Neuronal Activity	Yes	Yes	Yes

North America

Phone +1 734 769 1600

Europe

Phone +44 1763 227400

Email:

Japan

Phone +81 3 6478 5202

China

Phone +86 21 6878 2300

Rest of Asia Pacific and other countries around the world:

Phone +65 6872 3966