

Plate Catalog Number	Plate Name	Vol. in μ L	Corning Costar Plate Number	Plate Dimensions (in mm)				
				Well Depth	Plate Length	Plate Width	A1 Row Offset *	A1 Column Offset †
X001-1EA	Black Half Area 96 Well Plate	360	3915	10.67	127.8	85.5	11.2	14.3
X001-2EA	Black Half Area 96 Well Plates	360	3915	10.67	127.8	85.5	11.2	14.3
X003-2EA	Clear Microtiter Plates	360	9017	10.67	127.8	85.5	11.2	14.3
X003-5EA	Clear Microtiter Plates	360	9017	10.67	127.8	85.5	11.2	14.3
X004-1EA	Goat anti-Mouse IgG Coated White Solid Plate	360	3922	10.67	127.8	85.5	11.2	14.3
X010-1EA	Goat anti-Mouse IgG Coated White Strip Plate	360	3923	10.67	127.64	85.6	11.3	14.22
X011-1EA	Goat anti-Mouse IgG Coated Clear Solid Plate	360	9018	10.67	127.64	85.6	11.3	14.22
X012-1EA	Goat anti-Mouse IgG Coated Clear Strip Plate	360	2592	10.67	127.64	85.6	11.3	14.22
X013-1EA	Goat anti-Rabbit IgG Coated White Solid Plate	360	3922	10.67	127.8	85.5	11.2	14.3
X014-1EA	Goat anti-Rabbit IgG Coated White Strip Plate	360	3923	10.67	127.64	85.6	11.3	14.22
X015-1EA	Goat anti-Rabbit IgG Coated Clear Solid Plate	360	9018	10.67	127.64	85.6	11.3	14.22
X016-1EA	Goat anti-Rabbit IgG Coated Clear Strip Plate	360	2592	10.67	127.64	85.6	11.3	14.22
X018-2EA	Clear Half Area 96 Well Plates	190	3695	10.54	127.8	85.5	11.2	14.3
X018-4EA	Clear Half Area 96 Well Plates	190	3695	10.54	127.8	85.5	11.2	14.3
X018-5EA	Clear Half Area 96 Well Plates	190	3695	10.54	127.8	85.5	11.2	14.3
X023-1EA	Black Half Area 96 Well Plate	190	3686	10.54	127.8	85.5	11.2	14.3
X023-5EA	Black Half Area 96 Well Plates	190	3686	10.54	127.8	85.5	11.2	14.3
X025-1EA	Black 96 Well Plate	360	3650	10.67	127.8	85.5	11.2	14.3
X025-2EA	Black 96 Well Plates	360	3650	10.67	127.8	85.5	11.2	14.3
X037-2EA	Black Half Area 96 Well Plates	190	3694	10.54	127.8	85.5	11.2	14.3
X037-4EA	Black Half Area 96 Well Plates	190	3694	10.54	127.8	85.5	11.2	14.3
X060-1EA	Donkey anti-Sheep IgG Coated Clear Solid Plate	360	9018	10.67	127.64	85.6	11.3	14.22
X061-1EA	Donkey anti-Sheep IgG Coated Clear Strip Plate	360	2592	10.67	127.64	85.6	11.3	14.22
X062-1EA	Donkey anti-Sheep IgG Coated White Solid Plate	360	3922	10.67	127.8	85.5	11.2	14.3
X063-1EA	Donkey anti-Sheep IgG Coated White Strip Plate	360	3923	10.67	127.64	85.6	11.3	14.22
X110-1EA	Black 384 Well Plate	112	3575	11.43	127.8	85.5	8.99	12.12
X115-2EA	Black Low Volume 384 Well Plates	35	3676	6.58	127.8	85.5	8.99	12.12
C033-1EA	Mouse anti-Human Cystatin C Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C075-1EA	Mouse anti-Human OPN Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C107-1EA	PKA Substrate Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C140-1EA	Mouse anti-Human Prolactin Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C161-1EA	Mouse anti-Endothelin 1 Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C187-1EA	Mouse anti-Human Insulin Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C205-1EA	Mouse anti-Human ST2 Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22
C223-1EA	Mouse anti-Myeloperoxidase Coated Plate	360	2592	10.67	127.64	85.6	11.3	14.22

* A1 Row Offset is distance from top of plate to center of well A1

† A1 Column Offset is distance from left edge of plate to center of well A1