

We currently have laboratory equipment at reduced rates. This equipment will be sold on a first come, first served basis. This equipment may have been returned from a trade show, returned from a customer and refurbished, or acquired a scratch during production. Hurry, these items are popular and go fast.

List effective April 1, 2019

Model	Job/Serial No.	Product Number	Sales Price	Description
Shakers - MaxQ	300000571	SHKE6000-R	4635	- LIKE NEW - LIKE NEW -
LN2 Cryo -	300000863	7403TF-R	4800	- - -
LN2 Cryo -	300000864	7406-R	4800	- - -
LN2 Cryo -	300006202	7404-R	4800	- SCRATCHES - SCRATCHES FRONT, BACK & SIDES -
Shakers - 0	300011536	SHKE7000-R	3838	NONE - SCUFFS/BLACK MARKS - LIKE NEW -
Shakers - MaxQ	300052501	SHKE6000-R	4635	NONE - BACK WALL SCUFFED - LEFT SIDE DENT -
Shakers - MaxQ	300052535	SHKE6000-R	4635	NONE - RIGHT SIDE WALL SCUFFED - AND PAINT SCUFFED OFF TOP - DOOR DENTED LEFT FRONT CORNER
LN2 Cryo - 0	300144053	7404-R	4800	NONE - LIKE NEW CONDITION - MARKS ON OUTSIDE OF UNIT - SEVERAL DENTS, SCRATCHES AND
Shakers - 0	300160920	SHKE6000-R	4635	NONE - BACK WALL SCUFFED - DENT LEFT SIDE -
CO2 Incubators - 0	300100419	3310-R	7847	NONE - LIKE NEW CONDITION - LIKE NEW CONDITION -
CO2 Incubators - 0	300185719	3310-R	7847	PENDING - PENDING - PENDING -
CO2 Incubators - 0	300185720	3310-R	7847	NONE - SCUFFED - LIKE NEW CONDITION -
CO2 Incubators - 0	300192028	3307-R	7062	NONE - LIKE NEW CONDITION - CONTROL PANEL DENTED - RIGHT CORNER DENTED
CO2 Incubators - 0	300210545	3310-R	7847	NONE - SCUFFED - RIGHT SIDE SCUFFED AND - SCRATCHED ALONG BOTTOM
CO2 Incubators - 0	300170793	3310-R	7847	NONE - LIKE NEW CONDITION - AND SMALL DENT IN FRONT NEAR - TOP OF CONTROL PANEL SCUFFED
CO2 Incubators - 0	300210544	3310-R	7847	NONE - LIGHT SCUFFS AND SCRATCHES - DOOR LIGHTLY SCRATCHED. - TOUCHUP
CO2 Incubators - 0	300170794	3310-R	7847	NONE - LIKE NEW CONDITION - CONTROL PANEL LEFT SIDE LARGE - DECAL ON BOTTOM OF DOOR DAMAGE
CO2 Incubators - 0	300210543	3310-R	7847	NONE - LIKE NEW CONDITION - OF DOOR. PAINT CHIPPED & TOUCH - SMALL DENT LOWER LEFT CORNER
CO2 Incubators - 0	300217486	370-R	3715	PENDING - PENDING - PENDING -
CO2 Incubators - 0	300190795	370-R	3715	PENDING - DENTS & SCUFFS INNER WALLS & - PENDING -
Shakers - 0	300239784	SHKE6000-R	4635	NONE - SHELVES SCUFFED - LIKE NEW CONDITION - SCRATCHES ON RIGHT AND
ULT Freezers - 0	300235364	TSE400A-R	7276	NONE - PENDING - DOOR DENTED BY CONTROL PANEL - LEFT SIDE SCRATCHED - TOUCH UP
CO2 Incubators - 0	300183317	3310-R	7847	NONE - LIKE NEW CONDITION - TOP LIGHTLY SCRATCHED -