

CX & CXR Series



CX & CXR Series

Cryo Express (CX) "dry shippers" are designed to safely transport a variety of materials at cryogenic temperatures. The unique absorbent material prevents a liquid spill if the unit is tipped over. Storage temperature inside the shipping cavity remains at approximately -190°C until the liquid nitrogen evaporates from the absorbent material.



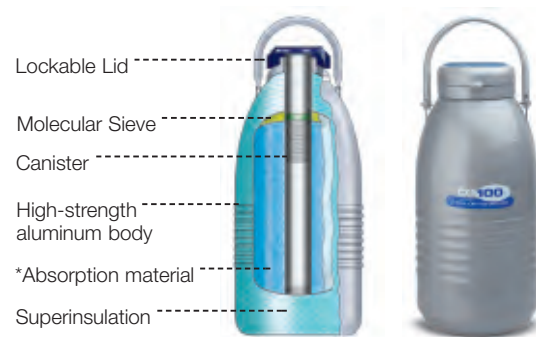
CX & CXR Series

Models	CX100	CXR100	CXR500
Static Holding Time days (1)	22	16	11
Working Time days (2)	17	11	7
Evaporation Rate (1) liters/days	0.19	0.23	0.64
Liquid Nitrogen Adsorbed - liters	4.1	3.7	7
Weight Empty	- lbs. 11.7 - kg. 5.3	11.7 5.3	30 13.6
Weight Full (3)	- lbs. 19 - kg. 8.6	18.3 8.3	42.5 19.3
Neck Diameter	- in. 2.78 - mm. 71	3.58 91	8.5 216
Overall Height	- in. 18.4 - mm. 467	19.4 493	26.9 683
Overall Diameter	- in. 9.2 - mm. 234	9.2 234	15.5 391
Number of Canisters	1	1	N/A
Canister Dimensions	- in. 2.64x11.0 - mm. 67x2.79	2.64x11.0 67x2.79	N/A N/A
Number of 1.2 ml & 2.0 ml vials (5/cane) (4)	85	85	500
Number of 1.2 ml & 2.0 ml vials (6/cane) (4)	102	102	500
Number of 1/2 straws (10/cane)	280	280	N/A
Number of 1/2 cc straws - Bulk (2 levels)	980	980	N/A
Padded carton	3701-9277	N/A	N/A
5 shelf rack	N/A	N/A	RS30-9C44
Poly Carb 100 box	N/A	N/A	R24K-9C44
Shipping Case	CX10-8C00	CX10-8C00	CP19-8C00

- (1) Evaporation rate and static holding time are nominal. Actual rate may be affected by the nature of the contents, atmospheric conditions, container history, and manufacturing tolerances.
 (2) Work time is an arbitrary, reference-only value to estimate container performance under the actual operating conditions.
 (3) Without canisters or racks
 (4) CXR500 vials are stored in 100 cell boxes

Features

- (1) Advanced Concept Absorbent enables faster charging
- (2) Complies with IATA regulations
- (3) New rugged construction includes – dependable, stronger necktube design, ribbed high strength aluminum body, and durable paint system
- (4) Superior vacuum performance with super insulation provides maximum holding times
- (5) Optional storage accessories include a five shelf rack box system and Blood Bag rack for the CXR500
- (6) Temperature loggers are available for all models
- (7) Lockable lids



*CXR Exclusive Feature: Removable absorption material



Taylor-Wharton

A legacy of excellence. A vision of promise.



Jena Bioscience GmbH
 Loebstedter Strasse 71
 07749 Jena, Germany

Ph.: +49-3641-6285000
 Fax: +49-3641-6285100
www.jenabioscience.com