

SmartNotes

What is the appropriate refrigeration unit to support storage of freshly expressed breastmilk?

Q
A

The TSG205 is designed to support the requirements of freshly expressed breastmilk.

There is nothing more important than ensuring the safe and effective storage of expressed breast milk. The storage of breastmilk needs to be managed in nourishment rooms, mother/baby rooms and neonatal intensive care (NICU) situations. The CDC recommends very specific storage conditions:

https://www.cdc.gov/breastfeeding/recommendations/handling_breastmilk.htm

The recommended storage temperature for breastmilk is +1 to +4°C, meaning it has an even more stringent temperature recommendations than vaccines or even blood. Breastmilk is often stored in close proximity to patients in the NICU and mother/baby suites, so noise from a storage unit should also be considered.





TSG205
countertop refrigerator

TSG205 with Phononic™ solid state technology allows for the most precise and consistent temperature control

The temperature range at which breastmilk should be stored (+1 to +4C) is very difficult for most types of refrigeration systems to manage given the cyclical nature of vapor compression-based refrigeration systems. However, the solid state technology used in the TSG205—which is powered by Phononic ensures that units can not only be set at the appropriate point but they can also maintain the proper range during normal operation—something other refrigerators cannot claim.

TSG205 with Phononic solid state technology allows for whisper quiet operation

The AAP recommends that the noise level in a NICU environment not exceed 45 dba (insert citation here from the AAP). The solid state technology that drives the TSG205 ensures operating noise levels less than 35 dba during operation, meaning you are always providing the safest environment for the most sensitive ears.

Summary: The TSG205 is the designed specifically for the safe storage and management of expressed breastmilk, supporting recommended temperature and noise conditions ensuring an optimum environment.



Capitol Vial Snappies™
Breast Milk
Container System

Visit thermofisher.com/whisperquiet for more information.

Australia: +61 39757 4300
Austria: +43 1 801 40 0
Belgium: +32 53 73 42 41
China: +800 810 5118 or
+400 650 5118
France: +33 2 2803 2180
Germany national toll-free: 0800 1 536 376
Germany international: +49 6184 90 6000

India toll free: 1800 22 8374
India: +91 22 6716 2200
Italy: +39 02 95059 552
Japan: +81 3 5826 1616
Netherlands: +31 76 579 55 55
New Zealand: +64 9 980 6700
Nordic/Baltic/CIS Countries:
+358 9 329 10200

Russia: +7 812 703 42 15
Spain/Portugal: +34 93 223 09 18
Switzerland: +41 44 454 12 12
UK/Ireland: +44 870 609 9203
USA/Canada: +1 866 984 3766

Other Asian Countries: +852 2885 4613
Countries not listed: +49 6184 90 6000