

EX-CELL[®] ANTIFOAM

Frequently Asked Questions

CATALOG NO. 59920C

What is the density of EX-CELL ANTIFOAM?

1.0273

What is the metabolic fate of EX-CELL ANTIFOAM?

Polydimethylsiloxanes (PDMS) are diverse polymetric compounds that are commonly known as silicones. PDMS make up a significant percentage of the raw material (Dow Corning[®] Q7-2587) used in EX-CELL ANTIFOAM. This compound does not bioaccumulate in living organisms because it is too large to be absorbed by biological membranes.

Why is EX-CELL ANTIFOAM sterilized by gamma irradiation versus autoclaving or sterile filtering?

The packaging used for this product is polyethylene (PE) media bags and it cannot tolerate the extreme temperature or pressure used in autoclaving. Antifoam is too thick to filter through a sterile filter. The PE media bags used during manufacture are not sterile prior to fill and the product is terminally sterilized so as to prevent the film from cross-linking due to two irradiations.

Are the polyethylene (PE) media bags used for EX-CELL ANTIFOAM considered as animal component free (ACF)?

Yes, the product contact layer in polyethylene bag is considered as animal component free.

How is the final product concentration tested and what test is used?

EX-CELL ANTIFOAM is tested with the use of a quantification assay and allows for the reporting of the active ingredient simethicone in the final product. The test name is Simethicone Emulsion and is based on the USP method. EX-CELL ANTIFOAM is the industry's only antifoam tested for accurate final product concentration.

Are there any special preparation instructions for use of EX-CELL ANTIFOAM?

EX-CELL ANTIFOAM is pre-diluted with water-for-injection (WFI) grade water, packaged into polyethylene (PE) media bags and sterilized by gamma irradiation. This product comes prepared and ready-to-use. The only recommendation is to vigorously mix the product prior to addition in cell culture. One to five minutes using a manual or mechanical rocking motion should achieve a homogenous mixture.

How is it proven that EX-CELL ANTIFOAM is sterile by gamma irradiation?

Spiking studies were performed to demonstrate a >6 log reduction of microorganisms. Dose mapping was also performed on the product to determine that the product is completely irradiated.

Are the PE media bags used for EX-CELL ANTIFOAM already irradiated prior to fill?

No, the bags are not irradiated prior to fill. This prevents the bags from being irradiated twice.

What is the storage temperature and shelf life of this product?

EX-CELL ANTIFOAM can be stored at ambient temperatures (15 to 30°C) and has a current shelf of 24 months.