

Frequently Asked Questions – Mammalian Cell Culture

Peptones

1. Are BD peptones processed in dedicated facilities?

BD has a dedicated Animal-Free (AF) processing area for select animal-free products (TC Yeastolate UF, Soy Peptone UF, and Yeast Extract UF) at our Sparks, MD facility. AF supplements will be manufactured at BD's dedicated animal-free/antibiotic-free facility in Miami, FL beginning in 2011.

2. What are potential sources of endotoxin?

Potential sources include

i) endotoxin present in the raw materials or ingredients used to manufacture the products,

ii) bacterial growth during the digestion process. To minimize endotoxin levels, BD offers ultra-filtered peptones as part of our product portfolio.

3. Do carbohydrates vary in composition with different peptones? Do they have an effect on glycosylation?

Each substrate has a unique carbohydrate structure profile. Glycosylation should be evaluated during and/or after the media development process.

4. Does BD provide trending data on your current peptones? What is the variation in the typical analysis data?

BD has size exclusion data and typical analysis information in the Bionutrients Technical Manual.

5. Are BD peptones GMO-Free?

BD does not certify GMO-free on our standard products.

6. Do increased concentrations of peptones in the basal medium contribute to increase in Osmolality?

Peptones are protein hydrolysates and consist of a mixture of peptides, vitamins, carbohydrates, salts, minerals and trace elements. Each peptone is unique in composition and may contribute to an increase in osmolality if supplemented incrementally in concentration because of rise in the total solute level in the medium.

7. Are BD peptones Halal and (or) Kosher certified?

No, they are not Halal or Kosher certified.

8. Does BD perform a mycoplasma test as part of the release criteria for animal free peptones?

Mycoplasma testing is not currently included as part of standard release testing, but may be requested as either a fee for service test or through our custom program.

BD Cell™ Medium

1. Does BD Cell MAb Media contain glucose? How much?

Yes, BD Cell media contains glucose, at 6.5 g/L.

2. Do BD Cell Serum Free and BD Cell Animal-Free media contain insulin?

Yes, it contains human recombinant insulin at 10 µg/mL.

3. Do BD Cell media contain L-glutamine?

Yes, BD Cell media do contain L-glutamine.



BD Select™ CHO Medium

1. Does Select CHO medium contain Insulin?
What expression system is used to produce Insulin?
Yes, it contains hr Insulin. The expression system is Yeast.
2. How much glucose is contained in Select CHO?
5-7 g/L of free glucose.
3. How much total protein is contained in Select CHO?
The only protein present is hr-Insulin, in the range of 9-10 mg/L.
4. Why does Select CHO contain HT?
Is there any specific reason?
The majority of the cell culture media contain Hypoxanthine and Thymidine, and nucleosides which act as intermediates for RNA/DNA synthesis.
5. Does Select CHO work well without Pluronic F68?
Yes, Select CHO will work in the absence of pluronic but single cell suspension culture performs well in an aerated system and pluronic helps cells stay in suspension.
6. Does Select CHO contain any chelating agent such as EDTA or EGTA.
There are no chelating agents present in Select CHO.

BD Select™ CD1000 Medium

1. Is there Pluronic F68 in the Select CD1000 medium?
No, it should be supplemented prior to use. A recommended concentration is 0.05-0.1%.
2. What is the protocol for preparing Select CD1000 from powder?
Please refer to the label or the technical sheet for preparation instructions.
3. What is the buffering system used in Select CD1000 medium?
The buffering system used in Select CD1000 medium is a bicarbonate and HEPES system.



BD Biosciences
Advanced Bioprocessing
800.638.8663
bdbiosciences.com/advbio