

# Frequently Asked Questions – Microbial Fermentation

---

1. 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.*

2. Are BD products GMO-Free?

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

3. Is it possible to replace bovine origin peptones used in the manufacturing process of vaccines with animal-free peptones?

*Yes, it is possible. The successful substitution is dependant on the nutritional requirements for the host bacterial cell or mammalian cell line. BD has been successful in assisting clients in this effort through our Media Optimization Service and our Custom Media Program.*

4. How safe are the bovine peptones used in the manufacturing of vaccines?

*BD animal origin peptones comply with our Animal Origin sourcing policy. Some animal origin peptones marketed in Europe also have a 'Certificate of Suitability' per EDQM. Additional custom testing or documentation can be provided on request.*

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

*No, they are not Halal or Kosher certified.*

6. Does BD perform a mycoplasma test as part of your release criteria with your 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 Biosciences**  
Advanced Bioprocessing  
800.638.8663  
[bdbiosciences.com/advbio](http://bdbiosciences.com/advbio)