

## Certificate of Origin

### Product Information

- a. Product number 21496
- b. Product name Di-2-ANEPEQ [JPW 1114]
- c. Country of raw material origin USA
- d. Country of final manufacture USA

### Manufacturing Information

- a. Is the product of synthetic origin?  Yes  No

- b. During the manufacturing process (consider all starting materials, reagents, chromatographic media, buffers, enzymes, fermentation broth, etc.) and packaging (product contact e.g. bottles caps, etc.) only synthetic origin materials were utilized?  Yes  No

If "no" is answered to either question a. or b. above, specify the origin of the materials used:  Animal  Human  Plant  
 Fermentation/cell culture origin  
 Natural\* - Inorganic  
 Natural\* - Organic  
\*Produced by nature; non-living (e.g. salts, minerals, crude oil fractions)

c. If plant or animal derived, specify the species \_\_\_\_\_

d. If animal derived, identify relevant animal health information (Provide data from test results and/or veterinarian, USDA/APHIS certificates.) \_\_\_\_\_

Animals exposed to or inoculated with any livestock or poultry disease agents exotic to the United States?  Yes  No

Animals from a facility where work with exotic disease agents affecting livestock or poultry is conducted?  Yes  No

e. If human derived, specify

Human material containing no animal material

- Preparation does contain animal derived additives  
Identify the additives \_\_\_\_\_
- Material is not of tissue culture origin
- Material is not zoonotic

Are there procedures in place to avoid cross contamination with residue of animal, human GMO origin or Allergen material that come into contact with the equipment used for manufacture of the product (e.g. other products and/or cleaning or disinfecting agents, media fills)?

- Not applicable (i.e. only synthetic origin materials and no Allergen come into contact with the equipment)
- Yes
- No

**To my best knowledge I certify that the above information is correct**

Released by:



\_\_\_\_\_  
Dr. Jennifer Liao, QC/QA Manager