

Certificate of Origin

Product Information

- a. Product number 22002
- b. Product name Calcein, AM *CAS 148504-34-1*
- 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