

Certificate of Origin

Product Information

- a. Product number 39062
- b. Product name Psoralen TMP Azide
- 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, USDAAPHIS 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 materials containing no animal material
- ☐ Preparation does contain animal derived additives
Identify the additives _____
- ☐ Material is not 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)?

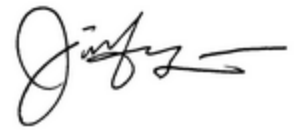
☒ Not applicable (i.e. only synthetic materials and no Allergen come into contact with the equipment)

☐ Yes

☐ No

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

Released by:



Dr. Jennifer Liao, QC/QA Manager