

Toxic Substance Control Act (TSCA) Certification

Date:

Waybill or reference number:

Check only one

Positive Certification

I certify that all chemical substances in this shipment comply with all applicable rules or orders under TSCA and that I am not offering a chemical substance for entry in violation of TSCA or any applicable rule or order thereunder.

or

Negative Certification

I certify that all chemicals in this shipment are not subject to TSCA.

Company name: Reaction Biology

Company address: One Great Valley Parkway, Suite 2

Certifier name: Haiching Ma

Certifier title: Chief Science Officer

Certifier phone number: 610 722 0247

Certifier email address: info@reactionbiology.com

Certifier signature:



Product description

1.
2.
3.
4.
5.
6.
7.
8.
9.
10.

If the certifier is unsure if their chemical substance is subject to TSCA compliance, contact the Environmental Protection Agency, TSCA Assistance Office at 1.202.554.1404 between 8:30 a.m. and 5:00 p.m.