

Anti Human Trafficking Modern Slavery Statement July 24th, 2024

Arbor Assays is committed to preventing modern slavery and human trafficking in our business operations and supply chains. As a small employee owned company, we recognize our responsibility to act ethically and with integrity in all our business dealings. This statement outlines the steps we have taken to ensure that slavery and human trafficking do not occur within our business or supply chains.

OUR COMMITMENT TO PREVENTING MODERN SLAVERY

While we may not have extensive resources for enforcement, we are committed to taking practical steps to minimize the risk of modern slavery in our operations and supply chains. Our approach includes:

Ethical Conduct: Promoting a culture of ethical conduct and integrity within our company.

Supplier Engagement: Communicating our commitment to preventing modern slavery to our suppliers and seeking their cooperation in adhering to our standards.

DUE DILIGENCE

We aim to address the risk of modern slavery through the following measures:

Supplier Vetting: We conduct basic vetting of our suppliers to ensure they align with our values and commitment to ethical practices.

Contractual Agreements: Where possible, we include clauses in our supplier contracts that require compliance with anti-slavery and human trafficking laws.

REPORTING CONCERNS

We encourage our employees, suppliers, and stakeholders to report any concerns or suspicions of modern slavery or human trafficking. Reports can be made directly to Braden@ArborAssays.com



COMMITMENT TO IMPROVEMENT

We are committed to continuously improving our practices to combat modern slavery. We will:

Review Practices: Regularly review our practices and seek feedback from employees and suppliers to identify areas for improvement.

Stay Informed: Stay informed about best practices and legislative changes related to modern slavery.

This statement has been approved by the Board of Directors of Arbor Assays.

Braden Robison

President & CEO