



United States Framework for Nucleic Acid Synthesis Screening

Self-Attestation

On behalf of Revvity, Inc. and its affiliates and subsidiaries which manufacture or sell synthetic nucleic acids, we hereby self-attest to our commitment to maintaining compliance with the US Framework for Nucleic Acid Synthesis Screening in accordance with the Executive Order on the Safe, Secure, and Trustworthy Development and Use of Artificial Intelligence.

We understand that our customers and federal funding agencies rely on this attestation prior to purchasing gene synthesis products. By implementing these measures, Revvity is reinforcing its commitment to ethical standards in synthetic biology.

Should the information above change, we agree that we will promptly provide updated information to the relevant parties.

If you have any questions about our compliance with these new regulations, please don't hesitate to contact our team. We value the trust you place in us and remain dedicated to providing safe and reliable synthetic biology solutions for the advancement of research.

A handwritten signature in blue ink, appearing to read "Joel Goldberg", is written over a horizontal line.

Joel Goldberg, SVP, General Counsel and Secretary

09 October 2024