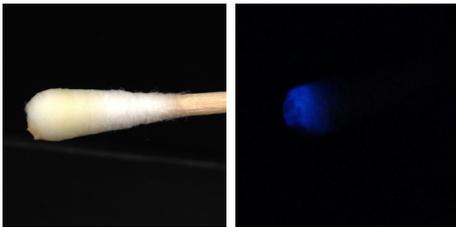


PRODUCT SPOTLIGHT

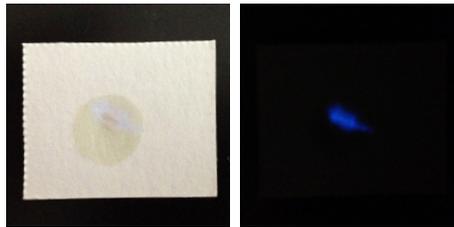
LABELED CONTROL STRAIN DETECTION

COTTON SWAB



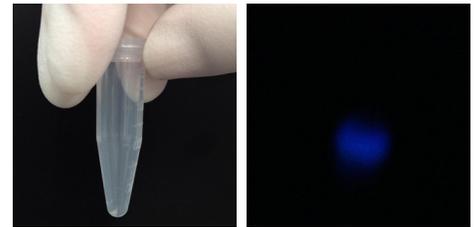
- 1 Prepare the reaction solution by mixing 1 μ L of Nano-Glo[®] Luciferase Assay Substrate (white cap) with 49 μ L of Nano-Glo[®] Luciferase Assay Buffer (yellow cap)
- 2 Swipe a cotton swab across the surface of a plate to collect colonies
- 3 Saturate the cotton swab with the reaction solution
- 4 Visualize in a dark area

PAPER



- 1 Prepare the reaction solution by mixing 1 μ L of Nano-Glo[®] Luciferase Assay Substrate (white cap) with 49 μ L of Nano-Glo[®] Luciferase Assay Buffer (yellow cap)
- 2 Saturate laboratory paper (e.g. Whatman[™]) with the reaction solution
- 3 Scoop 3-5 colonies onto a sterile loop
- 4 Scratch cells onto the saturated paper and visualize in a dark area

BROTH



- 1 Prepare the reaction solution by mixing 1 μ L of Nano-Glo[®] Luciferase Assay Substrate (white cap) with 49 μ L of Nano-Glo[®] Luciferase Assay Buffer (yellow cap)
- 2 Centrifuge 500 μ L of broth culture and remove the supernatant
- 3 Resuspend the pellet in reaction solution
- 4 Visualize in a dark area

