



May 26, 2023
BellBrook Labs
Madison, WI

Dear Berthold Technologies

This letter serves to certify that Tristar 5, a plate reader from Berthold Technologies has met our validation requirements for running Bellbrook Lab's Transcreener® FI assays. A Z' of 0.75 was obtained at 10% conversion when running a 10 μ M ATP/ADP standard curve. We are proud to add the Tristar 5 to our growing list of Transcreener® certified instruments.

Sincerely,
Justin Brink

Justin Brink
COO
BellBrook Labs
Madison, WI