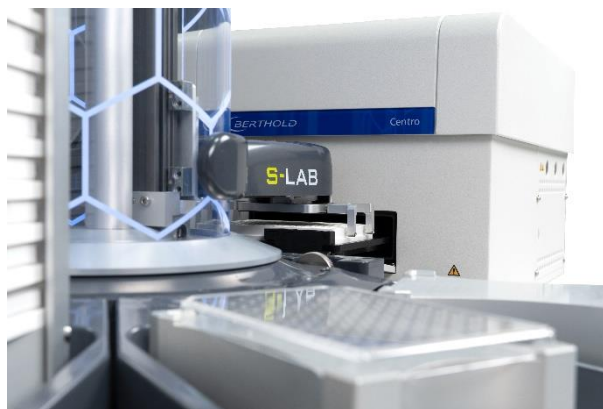


White Paper

INTEGRATION OF THE BERTHOLD TRISTAR & CENTRO WITH THE PAA S-LAB™ PRO AUTOMATED PLATE HANDLER FOR HIGH THROUGHPUT WORKFLOWS

Introduction

The need to coordinate complex tests, multiple workflows and time-consuming sample preparation pose significant challenges for laboratories. Here, laboratory automation can help by providing walk-away time, increasing throughput, and improving the effectiveness and efficiency of the testing process and reproducibility.



Berthold LB 942 Tristar 5 Multimode Reader

The Tristar 5 is a high-performance modular microplate reader. It has been specifically designed for high application flexibility and is therefore equipped with both independent, user-selectable filters and monochromators on both the excitation and emission side. The optical system thus allows great flexibility as well as high sensitivity, depending on what the application requires.

Berthold LB 963 Centro Microplate Luminometer

The Centro is a high-performance, easy-to-use, and versatile microplate luminometer for both, glow and flash luminescence assays. The optimized optical system of the instrument provides lowest background and negligible crosstalk (10^{-6}). It can be equipped with up to 3 injectors.



The S-Lab™ Pro automated plate handler attached to the Tristar 5 Multimode Reader

S-LAB™ Pro automated plate handler

The S-LAB™ Pro automated plate handler from Peak Analysis & Automation Ltd is an entry-level automation solution. It combines a compact plate handler (portrait or landscape) with the reliability of a robotic arm. The platform is capable of stand-alone operation, loading plates into automated microplate readers such as the Tristar or Centro.

The system allows unattended handling of up to 100 microplates (80 with lids). It can process lidded microplates, providing sterility for cell-based or sensitive assays and preventing evaporation. With a barcode reader and a delidding functionality, the S-LAB™ offers a reliable, flexible, and easy-to-use automation solution.

Easy setup using the Harmony™ user interface

The S-LAB™ Pro Harmony™ and Overlord™ software provide simple control via an intuitive user interface on a laptop, tablet or smartphone. Connection is via Ethernet.

PAA's Harmony™ user interface features an easy setup process for your assay. Setting up a run is straightforward, offering step-by-step guidance. Once you reviewed the entries, you just need to select run and walk away. The plates will be automatically loaded and measured into your Tristar or Centro Microplate Reader. Harmony will send you a message once the batch has been completed.

Summary

The Tristar and Centro together with the S-LAB™ Pro automated plate handler provide you with an automation platform optimised to work together

and offer a reliable solution for unattended sample analysis.

For Research Use Only. Not for use in diagnostic procedures.

© 2023 Berthold Technologies. All rights reserved. The trademarks mentioned herein are the property of Berthold Technologies or their respective owners. S-LAB™ PRO, Harmony™ and Overlord™ are a trademarks of Peak Analysis and Automation Ltd..

Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22

75323 Bad Wildbad

GERMANY

Phone: +49 7081 177 0

Email: bio@berthold.com



www.berthold.com