

For service purposes only

INTENDED USE

LANCE™ Controls are intended for *in vitro* use for testing instrument performance in LANCE measurements. The functioning of Wallac multilabel plate readers and microplate imagers can be verified with these controls.

PACKAGE CONTENTS

1 bottle (20 mL) of LANCE Positive Control (+)
1 bottle (20 mL) of LANCE Negative Control (-)

The Positive Control is a ready-for-use solution with Eu-W1024-biotin and SA-APC in polarization buffer.

The Negative Control is a ready-for-use solution with Eu-W1024-biotin and SA in polarization buffer.

The manufacturing date of the control is stated on the bottle label. Store in dark at +2 - +8°C.

PRODUCT USE

Allow the bottle to reach room temperature (+20 - +25°C) before use.

Take a clean black 96-well plate and dispense 200 µL of Positive Control to each well of one strip (e.g. 12 replicates) with a clean dispenser (not a glass pipette). Similarly dispense 200 µL of Negative Control to each well of another strip (e.g. 12 replicates). With a 384-well plate use a 50 µL dispensing volume, and with a 1536-well plate use a 5 µL dispensing volume. Measure the fluorescence within 30 minutes. When using the 1420 VICTOR™ Multilabel Counter, use a 665 nm filter and a typical window of 50-100 µs. When using the 1430 ViewLux™ Ultra HTS Microplate Imager, use a 671 nm filter. Count the signal-to-noise ratio (positive +/-negative -).

The avoidance of contamination and resulting high fluorescent background demands high standard pipetting technique. The same pipette tip must not be used for pipetting any other reagent. In case of test solution contamination, firstly use Enhancement Solution (prod. no. 1244-105) for cleaning, then wipe with a soft cloth or tissue paper soaked in clean water.

PRODUCT SPECIFICATION

The signal-to-noise ratio is typically over 100, when measured using a black standard microplate (prod. no. 1132 0129 or 1508-0010) and the VICTOR Multilabel Counter. Typically, the LANCE Positive Control gives a signal higher than 50000 counts, and the LANCE Negative Control gives less than 1000 counts.

November 2001

