



# Lyphocheck<sup>®</sup> Hemoglobin A<sub>2</sub> Control

**BIO-RAD**

